

USER MANUAL

CELLOMETER MATRIX™ SOFTWARE



8004707 Rev A

For Research Use Only. Not for use in
diagnostic procedures.



Cellometer Matrix™ Software User Manual

8004707 Rev A

Updated for Matrix v6.0.1 Software Release

May 2024

© 2024 Revvity, Inc. All rights reserved.

No part of this work may be reproduced in any form or by any means – graphic, electronic or mechanical, including photocopying, taping or information storage and retrieval systems – without the written permission of Revvity, Inc.

While every precaution has been taken in the preparation of this document, Revvity assumes no responsibility for errors or omissions, or for damages resulting in the use of information contained in this document or from the use of programs and source code that may accompany it. In no event shall the publisher and author be liable for the loss of profit, or any other commercial damage caused or alleged to have caused directly or indirectly by this document.

Cellometer™ and Cellaca™ are trademarks of Revvity, Inc.

Other products referred to in this document may be either trademarks and/or registered trademarks of their respective owners. Revvity makes no claim to these trademarks.

Table of Contents

CHAPTER 1. OVERVIEW	1
Matrix Software Overview.....	1
Available Software Run Modes	2
Live Mode	2
Data Analysis Mode	2
Simulated Mode	2
Minimum Computer Requirements	3
Matrix 21 CFR Part 11 Module Overview	3
Understanding Screen Variations between Instruments	4
Variations in Display of Consumable Formats	4
Variations in Load Process for Consumables	5
Variations in Preview Focus Controls	5
Variations in Default Assays/Report Templates	5
About this User Manual.....	6
How Screen Variations are Identified in this Manual	6
CHAPTER 2. MATRIX SCREEN ELEMENTS	7
Launching the Matrix Software	7
Matrix Screen Elements.....	8
Navigation Bar	8
Workflow Tabs.....	9
Simplified Workflow	11
User Favorites.....	11
Streamlined User Experience	11
Focus Methods	12
Manual focus.....	12
Image-Based Auto Focus	12
Slide-based Auto Focus	12
Viewing Software Version	13
CHAPTER 3. ACQUIRING SAMPLE DATA.....	15
Entering Setup Details	15
Entering Parameter Settings.....	16
Entering Well Details	17
Setting Auto Export Location.....	18

Previewing the Sample	19
Previewing Live Images.....	19
Adjusting Focus.....	19
Performing a Count	20
CHAPTER 4. ANALYZING SCAN RESULTS	21
Selecting Scan Results	21
Searching for a Scan Result.....	22
Viewing a Scan Result.....	23
Managing the Results List.....	23
Viewing Count Results	24
Understanding Default Report Tabs	24
Changing Well View Image Display	25
Varying Well View Channels/Counted Overlay Display	25
Understanding Custom Reporting	26
Verifying Auto Exports.....	26
Exporting and Printing Scan Results (Manual Settings).....	27
Performing a Recount.....	28
Refining Assay Details/Selecting New Assay	28
Managing Channel Mappings	29
Modifying Auto Export Options	30
Clicking the Recount Button	30
CHAPTER 5. CUSTOMIZING SCAN RESULT REPORTS.....	31
Viewing Custom Reports	31
Setting Reports and Exports for an Assay	32
Changing Template used in Well View Report Display.....	33
Editing Assay Exports.....	33
Managing the Assay Reports List.....	33
Adding Report Templates	34
Creating a Time Course Series	35
Best Practices and Workflow Tips	37
Creating a Time Course Report Template	38
CHAPTER 6. MANAGING FAVORITES	39
Viewing the Favorites List.....	39
Searching for Favorites	40
Maintaining the Favorites List	41
Importing Favorites	41
Exporting Favorites	42

Renaming Favorites	42
Deleting Favorites	42
Showing/Hiding Favorites in Selection Panel	42
Editing a Favorite	43
Modifying Favorite Details.....	44
Choosing a Consumable Type.....	45
Managing Favorite Reports and Exports	45
Creating a New Favorite	46
CHAPTER 7. MANAGING ASSAYS	47
Viewing the Assays List.....	47
Searching for Assays	48
Maintaining the Assays List	49
Importing Assays	49
Exporting Assays	50
Renaming Assays	50
Deleting Assays	50
Showing/Hiding Assays in Dropdown	51
Editing an Assay	51
Modifying Basic Assay Details.....	52
Defining Assay Imaging and Analysis Parameters	53
Imaging Mode Options	53
Analysis Mode Options.....	54
Dilution Factor	55
Channel Imaging and Cell Type Parameters	55
Managing Assay Reports and Exports.....	57
Creating a New Assay	60
CHAPTER 8. MANAGING CELL TYPES	61
Viewing the Cell Types List	61
Searching for Cell Types	62
Maintaining the Cell Types List.....	63
Importing Cell Types	63
Exporting Cell Types	64
Renaming Cell Types.....	64
Deleting Cell Types	64
Showing/Hiding Cell Types in Dropdown	64
Editing a Cell Type	65
Basic Cell Type Details	66

Brightfield Parameters.....	67
Fluorescence Parameters	69
Creating a New Cell Type.....	70
CHAPTER 9. MANAGING ACS TEMPLATES.....	71
Viewing the ACS Templates List	71
Searching for ACS Templates.....	72
Maintaining the ACS Templates List.....	73
Importing ACS Templates	73
Exporting ACS Templates.....	74
Renaming ACS Templates.....	74
Deleting ACS Templates.....	74
Showing/Hiding ACS Templates in Dropdown.....	75
CHAPTER 10. MANAGING REPORT TEMPLATES	77
Viewing the Report Templates List.....	77
Searching for Report Templates.....	78
Maintaining the Report Templates List	80
Importing Report Templates	80
Exporting Report Templates	80
Renaming Report Templates	81
Deleting Report Templates.....	81
Showing/Hiding Report Templates in Dropdown.....	81
Editing a Report Template.....	82
Editing Basic Report Template Details.....	83
Editing Report Elements	85
Creating A New Report Template.....	86
CHAPTER 11. USING THE 21 CFR PART 11 MODULE	87
Functionality Overview.....	87
Establishing Your User Hierarchy	88
Monitoring User Actions.....	88
Controlling User Access via Roles	89
Understanding Role Privileges.....	89
Data Privileges	90
Settings Privileges.....	90
Users Privileges	91
Roles Privileges.....	92
Audits Privileges	92

First-Time Login as Default Admin User	93
Creating an Administrator User	94
Managing Roles	95
Creating a Role.....	95
Copying a Role	96
Renaming a Role	96
Deleting a Role.....	96
Guidelines for Managing Roles.....	96
Managing Users	97
Creating a User	97
Disabling/Enabling a User	98
Resetting a User Password	99
Deleting a User	99
Guidelines for Managing Users.....	99
e-Signing Count Results	100
Viewing Audit Trail	101
Searching the Log	101
Exporting Log Details	101
CHAPTER 12. CONTACTING SUPPORT	103
Scope of Support Services	103
Contact Methods	103
Reporting an Issue to Support.....	103
Generating Diagnostic Reports.....	104
APPENDIX A. REPORT DESIGNER FOR WPF REFERENCE	105
APPENDIX B. SOFTWARE LICENSE DETAILS	107
Terms and Conditions.....	107

This page intentionally left blank.

Chapter 1. Overview

This chapter presents introductory concepts about the Matrix software including an understanding of available software run modes, minimum computer requirements, screen variations between instruments, and the intended scope of information contained in this manual.

MATRIX SOFTWARE OVERVIEW

The Matrix software used to run Revvity automated cell counting instruments is shared by the Cellometer (e.g., Ascend/K2) and Cellaca (e.g., MX/PLX) product families. In addition, default assays and report templates (both identified with instrument prefixes such as *ASD_*, *K2_*, *MX_*, or *PLX_*) are included for *all* supported instrument types. *Available assays and report templates for all instruments can be viewed in the **Manage** workflow tab.*

Note: When entering setup details, only assays for your specific instrument type will be displayed for selection. After a count is complete, report templates used in the presentation of scan results will be dynamically adjusted based on instrument consumable format variations (i.e., whether running multiple samples or a single sample).

Key elements in Matrix software functionality are described below:

- *Acquiring Sample Data* – The **Acquire** tab launches the two sequential screens – *Setup* and *Preview* – in the data acquisition workflow. Users can either select a *Favorite* (containing previously saved counting parameter settings), or enter setup details, select at least one well, and click the **Preview** or **Count** (displayed if *Skip Preview* feature is enabled) button. *Preview mode allows users to preview live channel images and adjust instrument focus/fluorescent exposure for each channel, while Count mode engages the camera to acquire sample images as specified by the assay.* The Matrix imaging and pattern-recognition software declusters, identifies, and counts individual cells according to defined cell type parameters.
- *Analyzing Scan Results* – The **Data** tab launches the three sequential screens – *Select*, *Results*, and *Recount* – in the data analysis workflow. Once a scan result is selected, users can vary the image displayed in the viewing pane by clicking available *Channel* buttons (e.g., **BR**, **FL1**, **FL2**, etc.) appearing across the top of the **Well View** tab or by choosing another well in the map. In addition, users can increase the magnification of the image and/or enable a graphic overlay to highlight counted cells. Data associated with the well appears below the image or can be viewed in a consolidated report including data from *all* wells by clicking the **Summary** tab.

To further refine the analysis of a scan, click the **Recount** button and either select a new assay, or click the **View** button to view parameters for the current assay and edit settings as necessary. *Parameters for the Last Used Assay are provided for reference.*
- *Setting Up Auto Exports* – The Matrix auto export feature allows you to select the *Images* (e.g., *Raw Images* and *Colorized Images*), *Data* (e.g., *Well Level CSV*, *Object Level CSV*, and *Object Level ACS*), and *Archive* (e.g., *Data Set*) output file types to be exported for an assay after completion of scan analysis. In addition, ACS templates can be imported into the Matrix software and assigned to an assay. On export, data will auto populate into the specified ACS template and can launch De Novo Software *FCS Express* if the *Auto Open* option is also selected and the application is installed.

- *Customizing Reports* – The Matrix custom reporting feature allows you to assign and/or modify report templates to be used by assays when generating scan results. You can change the report template used for displaying **Well View** tab data and manage the *Reports List* associated with an assay. You can also enable up to five additional tabs or specify output file types (e.g., *CSV*, *Excel*, *PDF*, and *Word*) to be automatically exported, opened or printed after completion of scan analysis.
- *Managing Favorites, Assays, Cell Types, ACS Templates, and Report Templates* – The **Manage** tab > *Favorites, Assays, Cell Types, ACS Templates, and Report Templates* options display lists of these entities currently loaded in your instrument system. From these screens you can import/export, rename, delete, or show/hide an entity in its list. For most entities you can also view, create, or use locked entities as templates to create new ones.

AVAILABLE SOFTWARE RUN MODES

In scenarios where an Operating Computer cannot always remain connected to its instrument, you can run the Matrix software in multiple modes to continue your analysis of captured sample images. In addition, you can export scan results to an alternative computer (e.g., workstation or personal laptop on which the Matrix software has also been installed) and perform sample analysis in a different location.

Note: Computers used to import and analyze scan result data when *not* connected to an instrument must meet minimum requirements (i.e., *Processor, Memory, and Operating System*) as noted below to run Matrix software in *Data Analysis* or *Simulated* modes. See the user manual for your instrument regarding technical specifications.

The following software *Run Modes* are available.

Live Mode

If the instrument and camera are physically connected to the Operating Computer via the Revvity-provided USB Connector Cable, the Matrix software runs in *Live Mode*.

Running the software in *Live mode* is necessary when acquiring sample data. Some **Acquire** tab functionality (e.g., adjusting focus and fluorescent exposure) is *not* available unless the software is running in *Live mode*.

Data Analysis Mode

If the instrument and camera are *not* physically connected to the Operating Computer *or* an alternative computer is used to continue your analysis of captured sample images, the Matrix software runs in *Data Analysis Mode*.

Simulated Mode

If the instrument and camera are *not* physically connected to the Operating Computer, a *Simulated Mode* feature uses previously stored simulated images to mimic camera functionality. *Simulated Mode* is only available to Revvity Sales/Field Application Specialists to demonstrate and test instrument functionality.

MINIMUM COMPUTER REQUIREMENTS

The following *minimum* requirements apply to Operating Computers used for running the Matrix software while connected to an instrument. *Some instruments include a Revvity-provided Laptop which may be more advanced.*

- Windows 10 PC
- Intel® i7, (2.10 – 4.80 GHz) Processor
- 16 Core, 24 Threads
- 32 GB RAM
- NVIDIA RTX A2000 8GB GDDR6 Graphics Card
- 1080p Display Resolution
- 1 TB+ Hard Drive
- USB 2.0 Port

In addition, these requirements also apply to any alternative computer (e.g., workstation or personal laptop) used to run the Matrix software in *Data Analysis* mode (i.e., while *not* connected to an instrument). *Scan results can be exported from the Operating Computer and imported onto an alternative computer for further analysis.*

MATRIX 21 CFR PART 11 MODULE OVERVIEW

The Matrix software offers a *21 CFR Part 11* module that is compliant with the *Code of Federal Regulations (CFR) Title 21 Part 11 – Electronic Records: Electronic Signatures* published by the U. S. Food and Drug Administration (FDA). In summary, this module ensures that an organization’s use of electronic records and digital signatures in place of traditional paper-based documentation is in compliance with current FDA regulations.

Key elements in Matrix *21 CFR Part 11* module functionality are described below:

- *Electronic Signatures* are captured during the counting and analysis workflow using **e-Sign/e-Unsign** buttons, and include the name of an authorized user, date/time when the signature was executed, reason for the action performed, and meaning of the signature (i.e., for the *Review, Approval, or Rejection* of an action). If a user’s assigned role does not have permissions to approve count results, a supervisor can log in to e-Sign the record on behalf of the user.
- *Electronic Records* are created as the result of linking electronic signatures to user actions ensuring that records have not been copied or falsified in any manner.
- An *Audit Trail* provides assurance regarding the integrity of an electronic record and continually monitors all users performing actions, the type of actions performed, and the date/time associated with user actions.

This module may be enabled for your instrument at time of purchase or implemented as an upgrade in the field by our Support team. See *Chapter 11. Using the 21 CFR Part 11 Module* starting on page 87 for a full description of module functionality.

UNDERSTANDING SCREEN VARIATIONS BETWEEN INSTRUMENTS

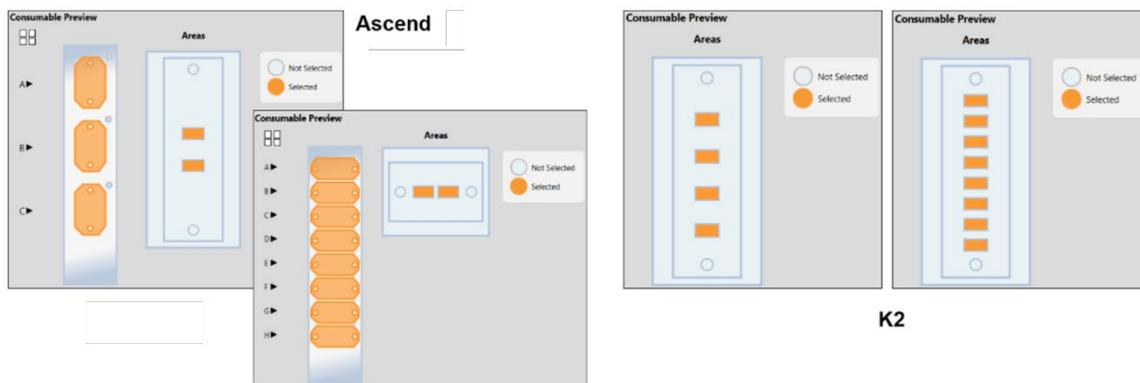
As the Matrix software is shared by multiple Cellometer products, it must account for physical differences between instruments, such as:

- Display of consumable formats (i.e., multi-well plates vs. chambered slides)
- Load process for consumables (i.e., clickable software buttons vs. physical insertion/removal process)
- Preview focus controls (i.e., automatic internal focus vs. physical external knob) and auto focus methods
- Default nature of assays and report templates (i.e., running and reporting on multiple samples vs. a single sample, and specifying the number of images to be captured per well)

To accommodate these differences, variations appear in the *Navigation Bar* and **Acquire** workflow tab screens, in assay file names, and in the display of generated reports based on the needs of each instrument. Although the visual presentation of these differences varies, the core functionality driving these tasks remains the same.

Variations in Display of Consumable Formats

The software display of Cellometer consumable formats is presented in the *Consumable Details* area of the **Acquire** tab Setup screen. The Matrix software supports multi-chambered slides for Ascend/K2 as shown below.



The visualization of “wells” on the chambered slides is depicted by *Images Per Well* (i.e., 1, 4, or 8) to be captured during the scan, with each image representing a cross-section of the counting chamber. *While Ascend slides are inserted completely into the instrument, only one chamber can be inserted at a time into the K2.*

Variations in Load Process for Consumables

For instruments that have an automatic feed mechanism (e.g., Cellometer Ascend), the *Navigation Bar* visible across the top of the screen displays an **Eject** or **Load** button that reflects the current position of the consumable. *The Eject button will become disabled while the instrument performs a count.*

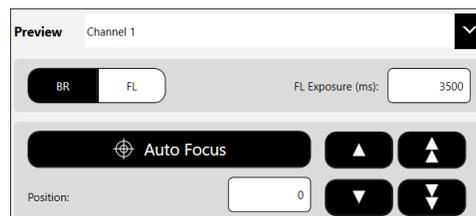


For Ascend, click the **Eject** button  to eject consumable or the **Load** button  to retract consumable. *Clicking the Eject or Load button toggles its display to the opposite action.*

No buttons are necessary for instruments that use a manual *Sample Slot* (e.g., Cellometer K2) as users can easily insert or remove consumable chambered slides from the slot located on the front of the instrument. *Only one chamber of a K2 slide can be positioned in the Sample Slot at a time.*

Variations in Preview Focus Controls

For instruments that use automatic internal focus (e.g., Cellometer Ascend), *Focus* controls are displayed while you are in the **Acquire** tab Preview screen. You can use the **Auto Focus** button to perform automatic focus adjustment or the *Coarse Focus/Fine Focus* controls to move through the focal planes of sample cells. In addition, you can specify a numerical value for the vertical position of the objective lens.



For instruments that have a physical *Focus Knob* located on the right side of the instrument (e.g., Cellometer K2), users can slowly turn the knob to manually adjust sample focus. Once good focus has been achieved during initial setup, the instrument should perform most counting operations with only minor adjustments.

Variations in Default Assays/Report Templates

Default assays and report templates that were common between instruments in previous releases have been optimized to accommodate for the display needs for specific instruments (i.e., whether you are running and reporting on multiple samples vs. a single sample). Although the core nature of these assays and report templates remains the same, minor variations exist in their presentation as noted below:

- Default assay and report template names include a prefix (e.g., “ASD_”, “MX_”, “PLX_”, or “K2_”) to identify instrument type.
- Default report templates are able to automatically adjust their presentation to display results from either multiple samples or a single sample based on your instrument type.

ABOUT THIS USER MANUAL

The intended scope of this *Cellometer Matrix Software User Manual* is to provide full software-related details regarding shared functionality as it has been applied across Cellometer products (e.g., Ascend/K2). This manual provides information on the following topics:

- [Matrix Overview and Screen Variations](#)
- [Screen Elements and Home Tab](#)
- [Acquire Tab Functionality](#)
- [Data Tab Functionality](#)
- [Custom Reporting Functionality](#)
- [Managing Assays and Assays List](#)
- [Managing Cell Types and Cell Types List](#)
- [Managing ACS Templates](#)
- [Managing Report Templates and RTs List](#)
- [21 CFR Part 11 Module Functionality](#)
- [Contacting Support and Reporting Issues](#)
- [Report Designer for WPF Reference](#)
- [Software License Details](#)

For hardware-related details regarding the setup, safety, care, and use of your instrument, refer to the specific *User Manual* and *Quick Start Guide* available for your instrument.

The following *Precaution Signifiers* are used in conjunction with the  symbol in this user manual:



IMPORTANT: Note indicating that to skip or move past *<content_of_note>* may result in improper functionality of the instrument.



CAUTION: Note indicating that *<content_of_note>* may damage instrument to the point where it will no longer function as expected.



WARNING: Note indicating that *<content_of_note>* may permanently damage instrument and cause personal injury or harm.

How Screen Variations are Identified in this Manual

As the Matrix software is shared by multiple product families, it must account for physical differences between instruments. To accommodate for these differences, screen variations are noted throughout this guide using instrument-specific sections introduced with a special statement that is *highlighted in italic font and orange text*. Be sure to choose the section containing information applicable to your instrument.

Chapter 2. Matrix Screen Elements

This chapter describes launching the Matrix software and presents basic graphical user interface (GUI) elements common to all instruments.

The following notes regarding instrument setup must be addressed *prior* to launching the software:

- Be sure to verify your system is set up as recommended in the unboxing and site preparation sections of the instrument's User Manual, and that the Operating Computer and instrument are both connected/powering on. In addition, follow all equipment safety protocols while using the instrument and keep the area around it clean before, during, and after operation.
- When connecting the computer to the instrument, users must wait until the instrument makes an audible click (i.e., indicating the instrument motors are communicating with the computer) before launching the software. Not waiting for this click can result in errors during the startup sequence. Keep in mind this note will apply each time the computer and instrument are disconnected/connected, or powered off/on again.

LAUNCHING THE MATRIX SOFTWARE

The Matrix software is shared by the Cellometer (e.g., Ascend/K2) and Cellaca (e.g., MX/PLX) product families.

If the Operating Computer running the Matrix software is a touchscreen device, you can interact directly with the graphical user interface (GUI) by tapping gently on screen elements (e.g., tabs, dropdowns, or buttons) using a finger or stylus. As an alternative, a USB mouse may be connected. *If you choose to use touchscreen functionality, the term "click" as it appears in this guide may be replaced with "tap" interchangeably.*

From desktop of Operating Computer, launch the Matrix software by double-clicking the **Matrix** icon. *The instrument will run through a startup sequence that includes connecting to the database.* The default screen for your instrument is displayed (e.g., the **Acquire** tab Setup screen).



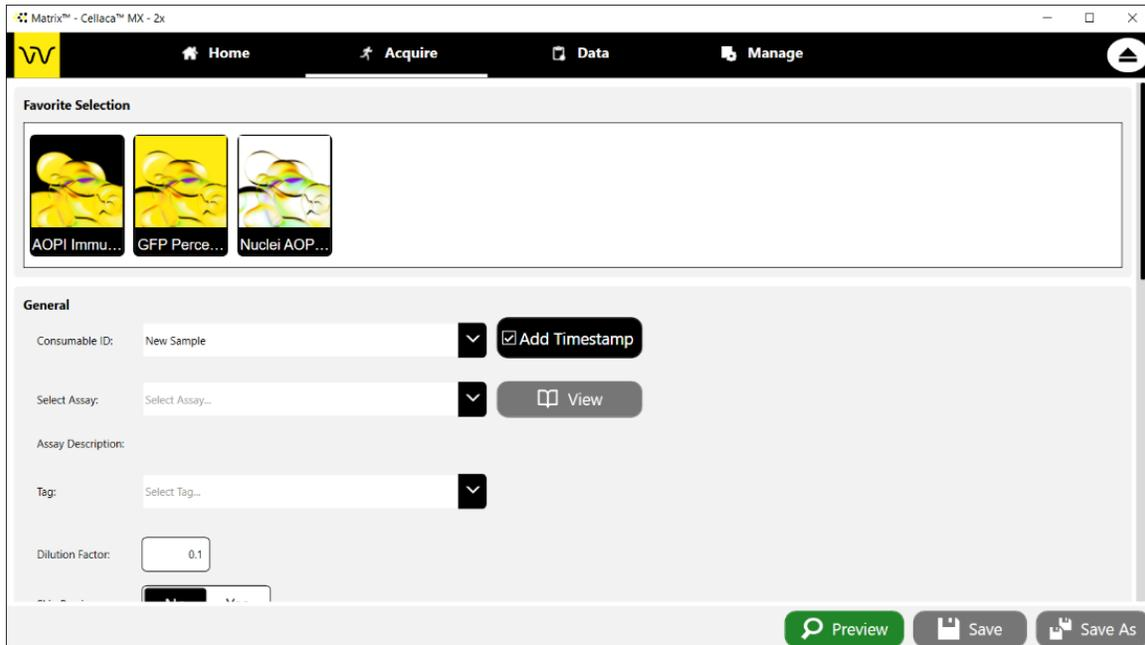
Note: If the Matrix 21 CFR Part 11 module has been enabled for your system, users must log in *before* they can use the software. See *Chapter 11. Using the 21 CFR Part 11 Module* starting on page 87 for module functionality.

After initial setup of an instrument, it must be calibrated using the Matrix software prior to first-time use. *You will be prompted with a message if you attempt to acquire data before calibration has been performed.*

The calibration process takes a background image that will be used to normalize the cell counter for each installed filter pair *without* a consumable counting loaded in the instrument. For assistance, contact Support by visiting <https://www.revivity.com/contact-us> or send email to: CellC-support@revivity.com

MATRIX SCREEN ELEMENTS

Upon launching the Matrix software you are presented with the default screen for your instrument (e.g., the **Acquire** tab Setup screen). Basic screen elements common to all instruments and their display variations based on instrument type are described below.



Navigation Bar

The *Navigation Bar* visible across the top of the screen is always displayed. *Elements displayed in this bar may vary based on instrument type.*



Functionality of Navigation Bar elements is described below.

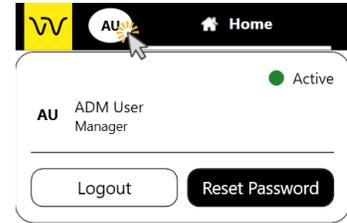
Workflow Tabs: The **Home**, **Acquire**, **Data**, and **Manage** tabs represent key areas of functionality as described in *Workflow Tabs* on page 9. Clicking a tab may display options for that tab (if available) in a sub navigation bar.

If the *Matrix 21 CFR Part 11* module is enabled, additional **Roles**, **Users**, and **Audit Trail** tabs are displayed.

Eject/Load Buttons: Available on *Ascend* only. Control movement of the automatic feed mechanism (i.e., ejects slide from instrument). *Eject* button becomes temporarily disabled while the instrument performs the count process.



User ID/User ID Card: If the Matrix 21 *CFR Part 11* module is enabled, the identity of the currently logged in user  is displayed. Clicking the User ID displays the User ID Card which contains the **Logout** and **Reset Password** buttons. See the *Matrix Software User Manual* for details on using module functionality.



Workflow Tabs

Functionality associated with each workflow tab is described below. Note that clicking the **Acquire** and **Data** tabs launches a series of screens that guide you through a sequential process, while options in the **Manage** tab may be individually selected to perform distinct tasks.



Home Tab: Displays the logo for the instrument to which you are connected and the Matrix software logo. Powered By  Matrix™

Also contains the **About Matrix** button which displays software version details and Revvity contact information, as well as the **Generate Diagnostic Report/Clear All Logs** buttons used to assist Support with troubleshooting technical issues. See *Viewing Software Version* on page 13 and *Generating Diagnostic Reports* on page 104 for more information.



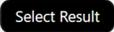
Acquire Tab: Selects the *Data Acquisition* workflow. In the **Acquire** tab you can either select a favorite (with all assay settings pre-defined), or enter *General* and *Consumable Details* information to select an assay, before advancing to preview the sample and confirm focus/ fluorescent exposure for each channel. See *Performing a Count* starting on page 20 for more information.

Note: The **Skip Preview** field selection is used to toggle the display of the **Preview** (i.e., **Skip Preview** is set to *No*) and **Count** (i.e., **Skip Preview** is set to *Yes*) buttons.

Use the **Back** button  located in bottom left corner of the screen to return to the previous screen in the process. *Clicking the **Acquire** tab while already in the workflow no longer returns to the previous screen as it did in earlier releases.*



Data Tab: Selects the *Data Analysis* workflow. You must first select a scan result to view its count results. You can then fine-tune assay parameter settings to perform a recount and create a new scan result. See *Chapter 4. Analyzing Scan Results* starting on page 21 for more information.

Use the **Select Result** button  located in bottom left corner of the screen to return to the previous screen in the process. *Clicking the **Data** tab while already in the workflow no longer returns to the previous screen as it did in earlier releases.*



Manage Tab: Contains the various system lists of favorites, assays, cell types, ACS templates, and report templates used in the *Data Acquisition* and *Data Analysis* workflows.

- In the *Favorites* screen you can view the library of favorites available in your system, import or export favorites, manage the list of favorites, as well as create new favorites or modify existing favorite details.

Favorites enable users to select previously saved counting parameter settings and instantly run samples without having to choose an application, thus bypassing the setup process and possibly the Preview screen (if *Skip Preview* feature is enabled). Scans that are performed frequently can be run with minimal software interaction.

- In the *Assays* screen you can view the library of assays available in your system, import or export assays, manage the list of assays, as well as create new assays or modify existing assay details.
- In the *Cell Types* screen you can view the library of cell types available in your system, import or export cell types, manage the list of cell types, as well as create new cell types or modify existing cell type details.
- In the *ACS Templates* screen you can view the library of ACS templates available in your system, import or export ACS templates, and manage the list of ACS templates. *ACS templates cannot be created or viewed in the Matrix software.*

ACS Templates can be used for populating data into specified formats when files are generated on export. If the *Auto Open* export option is also selected, the De Novo Software FCS Express application is launched for viewing the output file.

- In the *Report Templates* screen you can view the library of report templates available in your system, import or export report templates, manage the list of report templates, as well as create new report templates or modify existing report template details.

If the *Matrix 21 CFR Part 11* module has been enabled for your system, a user hierarchy must be established requiring the creation of roles that control varying levels of access to functionality available in each of these tabs. See *Chapter 11. Using the 21 CFR Part 11 Module* starting on page 87 for details.

Simplified Workflow

Simplified workflow features have been integrated into the Matrix software graphical user interface (GUI) to streamline user input in the most common data acquisition and analysis workflows.

USER FAVORITES

A key feature introduced in the v6.0 release is the creation of user favorites. A *Favorite* is a collection of count parameters such as an assay, consumable name/type, and reports/exports settings that can be saved for quick re-use. Users can assign images to be associated with favorites for quick reference, specify a prefix to be used for a series of consumables within an experiment, and enable the option of skipping the Preview screen (in favor of proceeding directly to the Count screen).

Favorites enable users to select previously saved counting parameter settings and instantly run samples without having to choose an application, thus bypassing the setup process and possibly the Preview screen (if *Skip Preview* feature is enabled). Scans that are performed frequently can be run with minimal software interaction.

Favorite Selection Panel: The *Favorite selection* panel is displayed across the top of the Setup screen and highlights either the last used favorite or the system default. To run the selected favorite, simply click the **Preview** (if enabled) or **Count** button to proceed.



To run a different assay, users can either scroll across the panel to select a different favorite or create a new setup by selecting an assay and/or entering new parameter settings.

STREAMLINED USER EXPERIENCE

Additional simplified workflow features built into the Matrix software may be used to streamline the overall user experience.

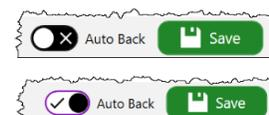
Add Timestamp Button: When entering setup details, clicking the **Add Timestamp** button automatically appends a timestamp to the Consumable ID when the scan result is created. *This button is enabled by default.*



Skip Preview Feature: When entering setup details, the **Skip Preview** field selection is used to toggle the display of the **Preview** (i.e., **Skip Preview** is set to **No**) and **Count** (i.e., **Skip Preview** is set to **Yes**) buttons.



Auto Back Button: If enabled, the **Auto Back** feature toggles functionality of the **Save** button to be **Save and Back**, automatically returning users to the previous screen when clicked. Click the **Auto Back** button  (located to left of **Save** button) to enable or disable this feature.



Single Folder Button: When entering setup details in the *Reports and Exports* area, the **Single Folder** button toggles between exporting files using the hierarchical folder structure traditionally available in earlier Matrix releases (button is *not* selected) or to a single folder (i.e., button is selected). *This button is also available when exporting scan results and is disabled by default.*



Back Button: While using the **Acquire** workflow tab, the Matrix software launches a series of screens that guide you through a sequential process. Use the **Back** button located in bottom left corner of the screen to return to the previous screen in the process. *Clicking the **Acquire** tab while already in the workflow no longer returns to the previous screen as it did in earlier releases.*



Select Result Button: While using the **Data** workflow tab, the Matrix software launches a series of screens that guide you through a sequential process. Use the **Select Result** button located in bottom left corner of the screen to return to the previous screen in the process. *Clicking the **Data** tab while already in the workflow no longer returns to the previous screen as it did in earlier releases.*



Focus Methods

Matrix software offers the following focusing methods.

MANUAL FOCUS

The traditional *Manual* focusing operation allows users to make adjustments to the initial Z-position using focusing controls (i.e., **Coarse Focus Manual Offset** and **Fine Focus Manual Offset** buttons) in the Preview Screen. This method uses an integrated *Focus Map* option to make small adjustments in the Z plane based on the consumable.

IMAGE-BASED AUTO FOCUS

The image-based auto focusing operation relies on the contrast within an image. Z-stacking is performed on the initial image, then the software analyzes objects in the image, selects images with the best contrast, and applies an algorithm to sharpen the focus further. Users have the option to perform image-based auto focusing for either the brightfield or fluorescent channels.

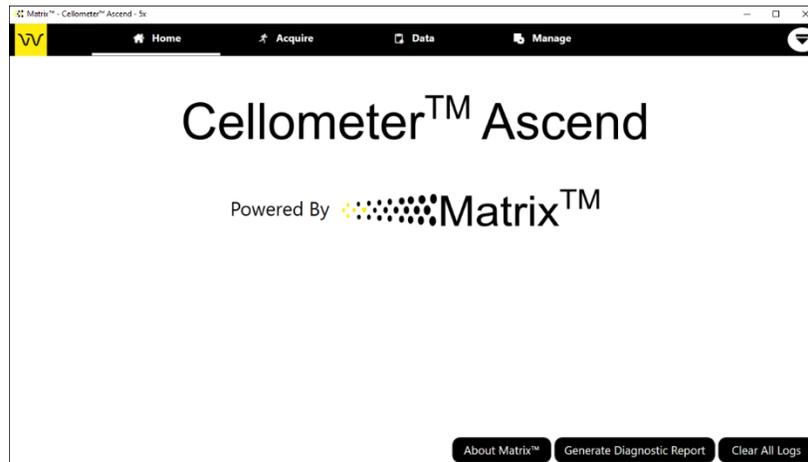
SLIDE-BASED AUTO FOCUS

Available for the Cellometer Ascend for the 3-Chamber Slide Only. The Slide-based auto focusing operation allows the software/instrument to focus on the consumable (without needing to focus directly on the sample) by using a target manufactured on the slide.

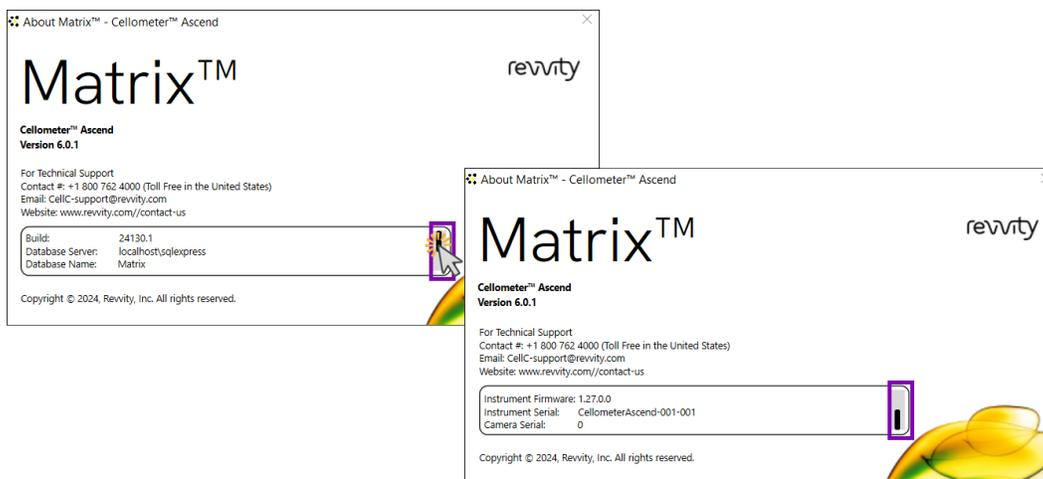
VIEWING SOFTWARE VERSION

The **Home** tab displays the logo for the instrument to which you are connected (e.g., Ascend as shown below).

For instruments that can be run using multiple versions of the software, the Powered By  Matrix™ logo distinguishes the Matrix software from previous versions available for the instrument.



The **Home** tab also contains the **About Matrix** button which displays product information (e.g., instrument type and software version), Support contact information, software details (e.g., build number, database server/name), and instrument details (e.g., instrument firmware version and instrument/camera serial numbers). *Users may need to click and drag the scroll bar to view all information.*



For details on using the **Generate Diagnostic Report** and **Clear All Logs** buttons to assist the Support team with troubleshooting technical issues, see *Generating Diagnostic Reports* on page 104.



WARNING: Logs may be required to maintain a historical archive. As using the Clear All Logs button will *permanently* remove accumulated logs for the installed Matrix software version, it is recommended that you contact IT *before* clearing logs from your system.

This page intentionally left blank.

Chapter 3. Acquiring Sample Data

This chapter describes functionality available in the two screens used to acquire sample data – *Setup* and *Preview*. After loading samples into the instrument, clicking the **Acquire** tab to launch the data acquisition workflow. Users can select a favorite (with all settings pre-defined), or enter *General* and *Consumable Details* information to select an assay, before advancing to preview the sample and confirm focus/fluorescent exposure for each channel.

- In the **Favorite Selection** area, you can choose a favorite and click **Preview** (if enabled) or **Count** to proceed with data acquisition.



- Upon entering setup details, selecting at least one well and then clicking the **Preview** button located at the bottom of the Setup screen, the instrument engages its camera for viewing live channel images of the sample. *If it is necessary to modify setup details while you are in the Preview screen, click the **Back** button to return to the Setup screen.*
- After adjusting instrument focus/fluorescent exposure for each channel, click the **Count** button at the bottom of the Preview screen. The instrument camera acquires sample images as specified by the selected assay which are then processed by the Matrix imaging and pattern-recognition software to decluster, identify and count individual cells according to defined cell type parameters.

Once a count has been performed, the scan result created as a result of the count is displayed at the top of the Results List. See *Chapter 4. Analyzing Scan Results* starting on page 21 for details.

ENTERING SETUP DETAILS

As the Matrix software is shared by multiple product families, it must account for physical differences between instruments such as the format of consumables. To accommodate for these differences, a variation exists in the *Well Details* area of the Setup screen as indicated by the instrument-specific sections below. Although the visual presentation of these differences varies, the core functionality driving these tasks remains the same.

Upon launch of the Matrix software, the **Acquire** tab *Setup* option is displayed by default. Personalize the details of your experiment by entering a consumable ID, selecting an assay, indicating specific wells for which to capture images (if applicable for your instrument type) and defining how the results data are to be exported.

Note: After initial setup, an instrument must be calibrated using the Matrix software prior to first use. If you are prompted with a message that the instrument has not been calibrated, contact Support by visiting <https://www.revvy.com/contact-us> or send email to: CellC-support@revvity.com

Entering Parameter Settings

The *Setup Details* area allows users to enter a customized plate name, select an assay and view/edit its details, select a tag, and enter a dilution factor.

The screenshot shows a 'General' section with the following elements:

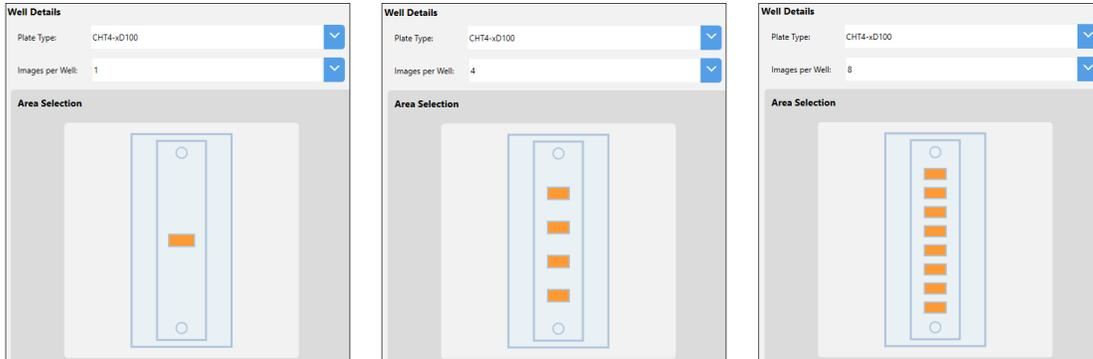
- Consumable ID:** A dropdown menu with 'New Sample' selected. To its right is a button with a checkmark icon and the text 'Add Timestamp'.
- Select Assay:** A dropdown menu with 'Select Assay...' selected. To its right is a button with a document icon and the text 'View'.
- Assay Description:** A text input field.
- Tag:** A dropdown menu with 'Select Tag...' selected.
- Dilution Factor:** A text input field containing the value '0.1'.
- Skip Preview:** Two buttons labeled 'No' and 'Yes'.

Users can enter the following information in the *Setup Details* area.

- Consumable ID** Customize the default plate name of “New Sample” to identify contents of the consumable. If a Consumable ID is *not* entered, a date/time stamp will be appended to the “New Sample” default (e.g., *New Sample 2023/08/25-10:58:09*). The **Add Timestamp** button toggles the timestamp append feature on and off.
- Select Assay** Select an assay from the dropdown and verify the assay description displayed. To view and/or edit assay details, click the **View** button. Click the **Back** or **Save and Back** (after making any changes) button to return to the Setup screen.
- Note:** If an assay is locked for editing (i.e., the **Category** field displays System or *Locked* instead of *Unlocked*), users must first click **Save As** to save it as a new assay before they can edit parameters. See *Editing an Assay* on page 51 for details.
- Tag** Select a tag from the dropdown or enter a new one. Tags are used to logically group scan results together for the purpose of custom reporting (e.g., Time Course Analysis). Tags have a maximum length of 32 characters and can be applied to any number of results, but can only be applied to one result derived from each scan. See *Creating a Time Course Series* on page 35 for details on how tags can be used to create custom reports.
- Dilution Factor** Enter the final dilution factor for the sample.
- Skip Preview** In the **Skip Preview** field, select **No** to enable the **Preview** button or **Yes** to skip previewing the sample and proceed to performing the count

Entering Well Details

The *Well Details* area for the Cellometer Ascend/K2 identifies the consumable **Type** used for the sample and allows you to select the loaded wells and number of **Images Per Well** (i.e., 1, 4, or 8) to be captured during the count, as shown below. *For the Cellometer Ascend, the default well selection has all wells selected.*



The visualization of *Images Per Well* for the slide depicts the selected number of images (i.e., cross-sections of the area in the counting chamber or “well”) to be captured during the scan. A comparison of *Total Volume Measured* for the Cellometer Ascend and K2 in relation to hemocytometer quadrants is provided below.

Total Volume Measured For Ascend Images Per Well:

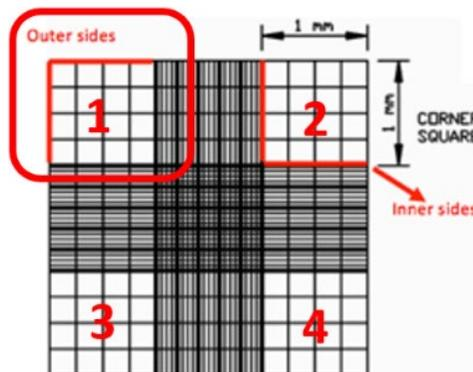
- Count from 2 images per well = 1.7 μL (or 17 quadrants of a hemocytometer)
- Count from 4 images per well = 3.4 μL (or 34 quadrants of a hemocytometer)
- Count from 8 images per well = 6.9 μL (or 69 quadrants of a hemocytometer)

Total Volume Measured For K2 Images Per Well:

- Count from 1 image per well = 0.15 μL
- Count from 4 images per well = 0.6 μL (or 6 quadrants of a hemocytometer)
- Count from 8 images per well = 1.2 μL (or 12 quadrants of a hemocytometer)

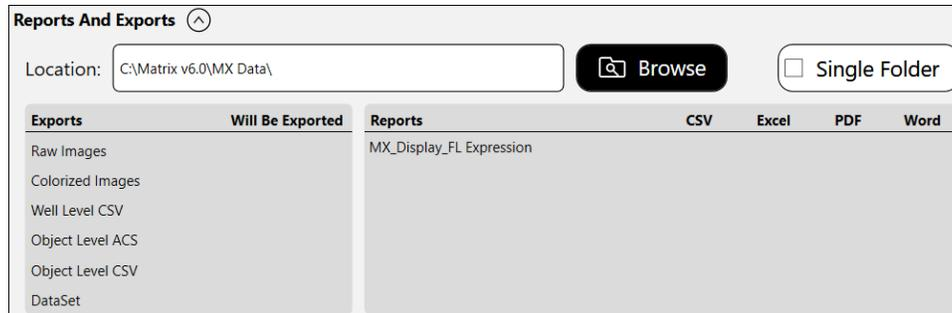
Total Volume Measured For Hemocytometer Quadrants:

- Count for 1 quadrant = 0.1 μL *Volume Per Quadrant* using an average of 4 quadrants result in 0.4 μL *Total Volume Measured*



Setting Auto Export Location

Expand the *Reports and Exports* area to identify a **Location** for automatic exports of images/data and generated output files for reports. Click the **Browse** button and navigate to a folder on your Operating Computer or network to define the default export path. *This path will remain as the default in the software until it is manually changed.*



In addition, *Exports* (see descriptions below) and output files for *Reports* (e.g., *CSV*, *Excel*, *PDF*, and *Word*) defined for the current assay are displayed.

Sample *Exports* include:

- Raw Images** Represents *Black and White* high-resolution PNG images for each channel in the assay.
- Colorized Images** Represents colorized, high-resolution PNG images of all acquired channels. *A single colorized image represents individual channel images superimposed one on top of another.*
- Well Level CSV** Represents well-level data including well name, calculation run, channel, count, mean size, etc. for each selected well in Comma Separated Values (CSV) format.
- Object Level ACS** Represents object-level data for each well including type, classification, size, circularity, area, perimeter, etc. in Image Cytometry Experiment (ICE) format.
Exported data may be associated with an ACS template and automatically opened for viewing in De Novo Software FCS Express (if specified in auto export options for the assay).
- Object Level CSV** Represents object-level data for each well including type, classification, size, circularity, area, perimeter, etc. in Comma Separated Values (CSV) format.
- Data Set** Represents a database file containing all images, results, assays, cell types and report templates associated with the scan result. *A data set is required to reload data for analysis or if you need help from Support to optimize assay and/or cell type parameters.*

To manage *Exports* and *Reports* defined for the selected assay, click the **View** button and expand the *Reports and Exports* option. Select *Exports* by clicking on file type buttons and manage *Reports* as necessary. Changes to the assay can either be saved to the current assay (i.e., any changes made will also be applied to other scan results that use the assay) or saved as a copy with a new name. Edited assays can be used for data acquisition. See *Managing Assay Reports and Exports* on page 57 for details.

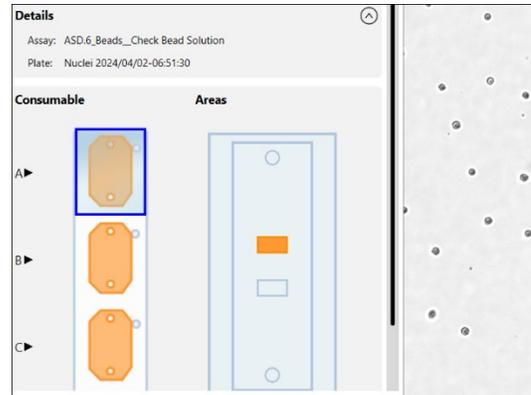
PREVIEWING THE SAMPLE

Once you have completed entering setup details for your sample, click the **Preview** button located at the bottom of the Setup screen (if enabled). The instrument engages its camera for viewing the sample and displays the Preview screen. Users can view live images of samples in selected wells, preview available channels for *Imaging Mode* associated with the assay, adjust instrument focus, and confirm fluorescence exposure for each channel.

Previewing Live Images

In the *Consumable* area, click on highlighted image sections in the slide visualization (i.e., representing cross-sections of the counting chamber or “well”). As you move from section to section, the live image changes per your selection.

To zoom in/out of an image, move the mouse to hover cursor over the viewing pane and turn the scroll wheel or, if using the touchscreen, apply universal gestures (e.g., touch center of image with two fingers and then spread them apart to zoom in and reverse this action to zoom out.) Current *Zoom* magnification is displayed in bottom right corner of the viewing pane.



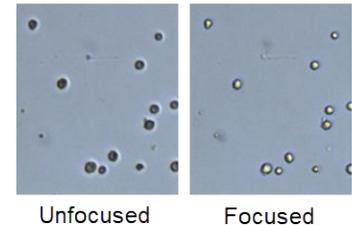
To move a zoomed image around, click and drag the image to a new location as needed.

Adjusting Focus

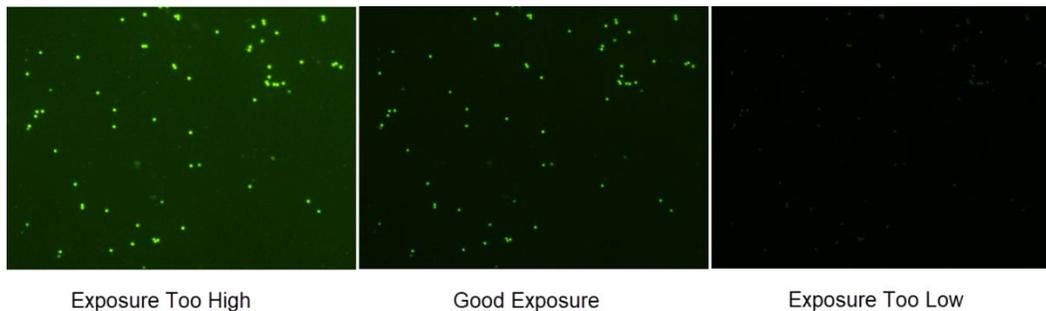
To adjust the focus of the live image being previewed, slowly turn the Focus Knob located on the right side of the instrument until cells/beads are in focus.

Obtaining good focus is key to ensuring accurate cell counts. Once good focus has been achieved, the instrument should perform most counting operations with only minor adjustments.

Once focus has been properly adjusted, continue by reviewing channel images.



When previewing fluorescent images, confirm that FL signal is strong but has a low, dark background. Modify the **FL Exposure** default to increase (prolong) or decrease (shorten) the exposure time accordingly.



PERFORMING A COUNT

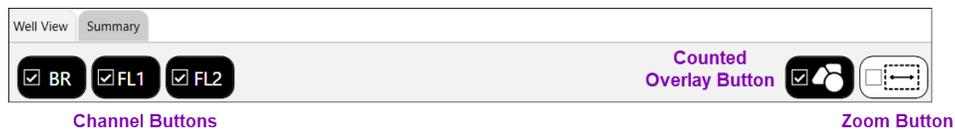
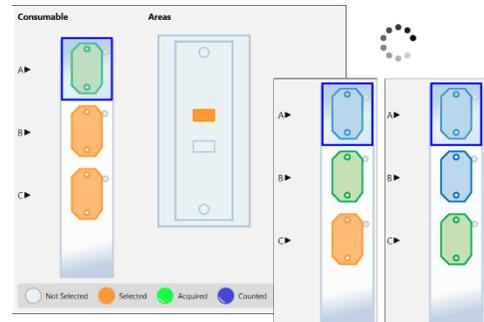
Once you have completed previewing the live image for the sample, click the **Count** button located at the bottom of the Preview screen. The instrument camera acquires sample images as specified by the selected assay which are then used by the Matrix software to calculate count results according to defined cell type parameters.

Note: Depending on the number of selected wells/images per well and defined assay parameters, the counting process can take a few seconds to up to a few minutes.

As the system acquires sample images and calculates count results, the colors used to mark selected wells will change to indicate status (i.e., from *Selected* to *Acquired* to *Counted*) as shown in the legend appearing below the Well Map.

As soon as a well is *Counted*, you can click on it to display count results below the viewing pane. *Count results will be displayed, printed and exported based on templates defined for the assay.*

Well images displayed can be varied by toggling on/off available channel buttons (displayed across the top of the viewing area of the **Well View** tab) and enhancing the *Zoom* magnification.



To zoom in and out of sample images, move the mouse to hover the cursor over the viewing pane and turn the scroll wheel or, if using the touchscreen, apply universal gestures (e.g., touch the center of the image with two fingers and then slowly spread them apart to zoom in and reverse this action to zoom out). Current magnification is displayed in bottom right corner of viewing pane.

When counting of all wells is complete the scan result will be added to the top of the Results List displayed on the Select screen. In addition, *Reports and Exports* output files defined for the current assay are automatically stored in the specified location. See *Chapter 4. Analyzing Scan Results* starting on page 21 for details.

When viewing count results, a single image is displayed representing the sample in the outlined chamber. Click on other highlighted chambers to view their sample images. If the consumable format allowed you to select the number of *Images Per Well*, you can click on any area (in the Consumable Areas visualization) to view the image of that area.

To move an image around, click once inside the viewing pane and drag the image to a new location as needed.

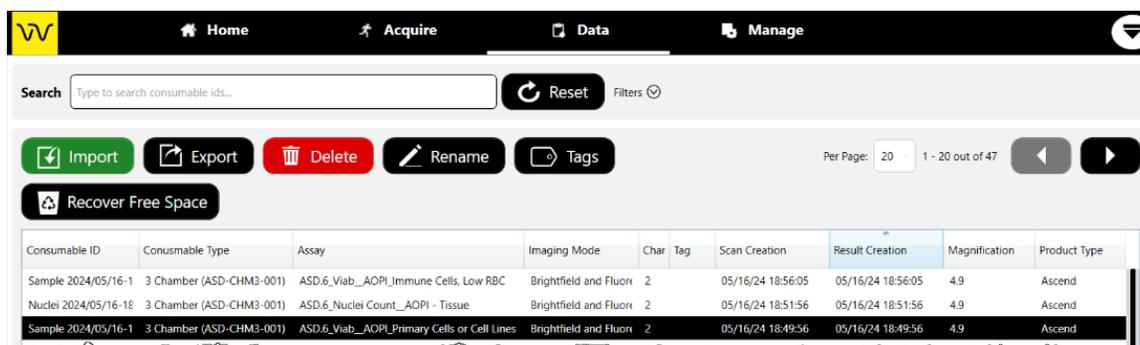
Chapter 4. Analyzing Scan Results

This chapter describes functionality available in the three screens used to analyze scan result data – *Select*, *Results*, and *Recount*. To view scan results, click the **Data** tab to launch the data analysis workflow. Users must first *Select* a scan result before you can view its count *Results*. You can then perform a *Recount* by fine-tuning assay parameter settings to be used in the analysis to create a new scan result.

- To open a scan result contained in the *Results List* of the Select screen you can either double-click the result, or click it once (to highlight it in the list) and then click the **View** button. See *Selecting Scan Results*, below.
- Once a scan result is displayed in the Results screen you can analyze it by selecting/de-selecting channels to vary the image and click on highlighted wells to review the data. *If you find it necessary to choose another scan result, click the **Back** button to return to the Results List.* See *Viewing Count Results* on page 24 for details.
- To perform a *Recount* you can select either a new assay or click the **View** button for the current assay to edit its parameter settings, and then click the **Recount** button. See *Performing a Recount* on page 28 for details.

SELECTING SCAN RESULTS

The Select screen displays all *Scan Results* currently generated or loaded into the database for your system. From this screen you can search for, select and view scan results. In addition, scan results can be imported/exported, deleted or renamed. If scan results are deleted, users can recover free space available in the database.

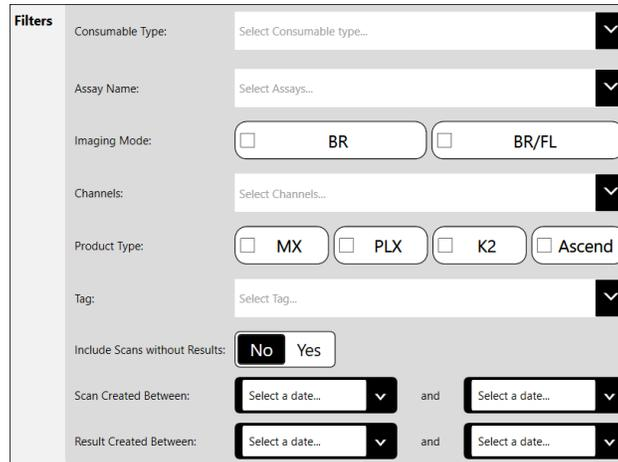


Use the **Per Page** control dropdown (contains values from 5 to 500) located in the upper right corner of this screen to select the number of scan results to be displayed per page, and the arrows to move back and forth between list pages.

See *Managing the Results List* on page 23 for more information on managing scan results contained in this list.

Searching for a Scan Result

To search for a scan result, enter a few key characters from a plate name into the **Search** field and/or expand the **Filters** area by clicking the down arrow  to specify search criteria.



Filters

Consumable Type: Select Consumable type... ▼

Assay Name: Select Assays... ▼

Imaging Mode: BR BR/FL

Channels: Select Channels... ▼

Product Type: MX PLX K2 Ascend

Tag: Select Tag... ▼

Include Scans without Results:

Scan Created Between: Select a date... ▼ and Select a date... ▼

Result Created Between: Select a date... ▼ and Select a date... ▼

Refine your search by selecting a plate type or assay name, clicking on an imaging mode or product type, selecting a tag, indicating whether to include scans without results, or entering a range of dates between which a scan and/or result was created. As you enter search criteria, the Results List is updated with matching entries.

If you find it necessary to clear all selected filters and begin the search again, click the **Reset** button.

Note: The Matrix software is designed to support multiple product families (e.g., Cellometer Ascend/K2 and Cellaca MX/PLX) and displays *all* available scan results in the Results List regardless of whether the Operating Computer is physically connected to that instrument.

Viewing a Scan Result

To open a scan result contained in the *Results List* you can either double-click the result or click it once to highlight it in the list and then click the **View** button. The Results screen is displayed.



Note: As *all* available scan results are contained in the Results List, you may be able to view scan results taken by an instrument that is different from the one to which your system is currently connected.

MANAGING THE RESULTS LIST

The following functionality is available when viewing the Results List.

- Import** Allows user to import previously stored scan results (stored as *.SCANRESULT* files) from an external location.
- Export** Allows users to export selected scan results (to be stored as *.SCANRESULT* files) to an external location.
- Delete** Allows users to delete the selected scan result. *If the Matrix 21 CFR Part 11 module is enabled, users will be prompted to enter a reason prior to deletion.*
After data has been deleted, click the **Recover Free Space** button (if displayed) to recover space that the data previously occupied in the database.
- Rename** Allows users to rename the selected scan result. *New plate name entered must be unique.*
- Tags** Allows users to select a tag from the dropdown or enter a new one to apply it to the currently selected scan result. *Tags can be used to create custom reports.*

Recover Free Space Button	<p>Allows users to recover free space that continues to be occupied in the database even <i>after</i> scan result data previously stored in that space has been deleted.</p> <p>Note: When scan results are deleted, the database does not immediately return space previously used by deleted data to the operating system. Instead, this process occurs gradually over time. The Recover Free Space button allows users to return unoccupied database space to the operating system at the time the operation is performed.</p> <p><i>Users must have System Admin rights for the database for this button to be displayed. In addition, as this process requires high CPU utilization during operation it is recommended that you not click perform other functions until this operation is complete.</i></p>
View	<p>Allows users to open the selected scan result for viewing.</p>

VIEWING COUNT RESULTS

Once you have opened a scan result for viewing, you can view count results presented in the Results screen.

Understanding Default Report Tabs

Report tabs initially displayed across the top of the viewing pane are associated with report templates that have been assigned to the current assay and are enabled by default to format the presentation of scan result data.



Although the **Well View** tab will always be enabled for display, you can change the report template assigned to this tab. In addition, you can disable any other default report tabs (e.g., **Summary** and **.csv** tabs for Matrix v4.0+; **All Wells** tab for Matrix v3.0), change the report templates used for display of these tabs, and/or add new tabs. See *Chapter 5. Customizing Scan Result Reports* on page 31 for details.

Note: As of the Matrix v4.0 release, users can customize report tabs displayed when viewing count results. Tabs enabled for display by default have been changed as noted below:

- The **Well View** report tab available in earlier releases will always be enabled. However, users can change the *Display* report template assigned to this tab or edit the template directly. *Any changes made to a report template in use by other assays will also be applied to other scan results the next time a recount is performed.*
- The **All Wells** tab available in earlier releases has been replaced with the **Summary** tab. In addition, a new **.csv** tab may be enabled. Users can change report templates assigned to these tabs, edit the templates directly (i.e., any changes made will also be applied to other scan results using these templates), disable them from displaying as tabs or delete them from the assay.
- Tabs enabled for display by default are identified by version numbers included in an assay name (e.g., **MX405.0_AOPI_Cell Lines** will include default tabs enabled in the Matrix v4.0 release).

CHANGING WELL VIEW IMAGE DISPLAY

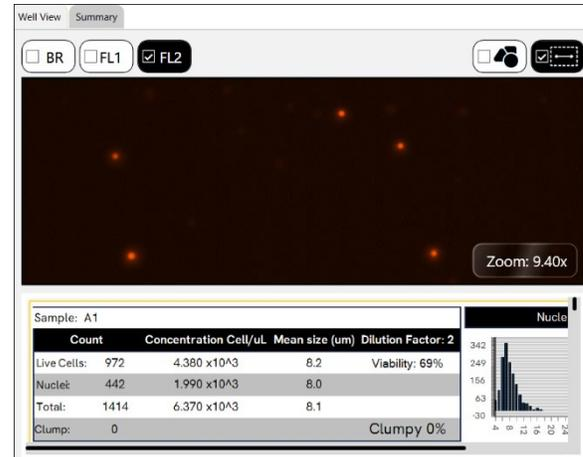
The **Well View** tab will always be displayed in the Results screen, and includes both an image and the associated count results for the selected well.

To move the image around, click and drag to a new location.

You can zoom in/out of the well view image by clicking once in the viewing pane and turning the mouse scroll ball or, if using a touchscreen, by applying universal gestures (e.g., touching center of image with two fingers and then spreading them apart to zoom in and reverse this action to zoom out).

Zoom magnification is indicated in bottom right corner of the viewing pane and can be increased up to 10.00x.

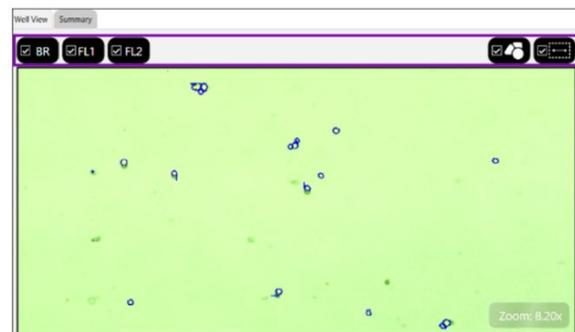
At the bottom of the **Well View** tab is a report containing well-level data for the sample in the selected well.



VARYING WELL VIEW CHANNELS/COUNTED OVERLAY DISPLAY

In the **Well View** report tab displayed by default, the following buttons may be available across the top of the viewing pane based on the Imaging Mode defined in the current assay.

- Click **Brightfield (BR)** or **Fluorescence** (e.g., **FL1**, **FL2**, etc.) buttons to select/de-select channels used in the image display. *Channel views are overlaid on top of each other.*
- Click the **Counted Overlay** button to show/hide the graphic overlay that identifies *Counted* cells by surrounding them with color-coded outlines. *For 2-channel Viability assays, Green is used for counted/live cells, Red for dead cells, and Yellow for cells not counted (e.g., if larger than the specified cell diameter). For Expression assays, Blue is used for outlining total cells in the masked channel.*



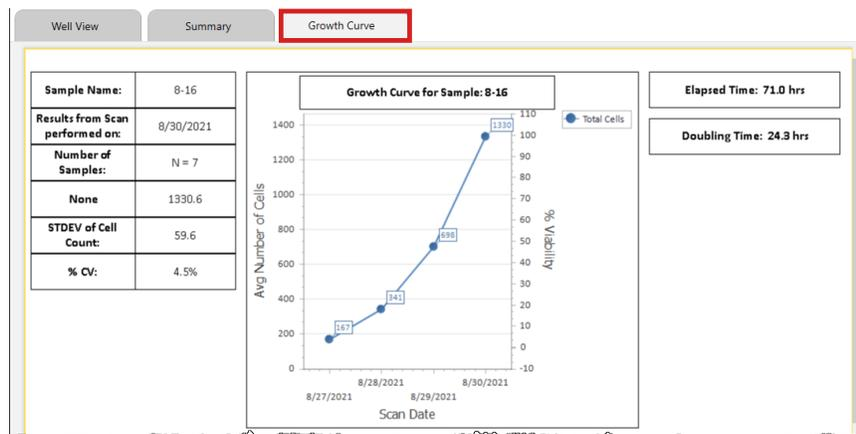
- Click the **Zoom** button to enable/disable display of current *Zoom* magnification in bottom right corner of viewing pane. *Zoom feature will still be functional even if not displayed.*

Understanding Custom Reporting

The Matrix custom reporting feature allows you to assign and/or modify report templates to be used by assays when generating scan results. You can change the report template used for display of the **Well View** report (which is always displayed), and add report templates to be enabled as tabs or output as specific file types (e.g., *CSV*, *Excel*, *PDF*, and *Word*) that can be opened/printed upon report generation.

Note: Any changes made to *Reporting* options for an assay will be saved with the assay and applied to all other scan results using that assay the next time a new scan or recount is performed.

For example, a custom **Growth Curve** report tab has been enabled in the scan result shown below. Clicking this tab displays data in the format defined by the associated report template (i.e., illustrating calculation of a growth curve for a time course series of scan results).



See *Chapter 5. Customizing Scan Result Reports* on page 31 for details on customizing *Reporting* options.

Verifying Auto Exports

Expand the *Exports* area in the Results screen to verify that automatic exports were completed and click the location link to open folder where the exported scan result files are stored.

In addition, if you defined any output file types (e.g., *CSV*, *Excel*, *PDF*, and *Word*) to be generated for assay report templates, you can verify that selected file type exports were also completed.

Exports C:\Matrix v6.0\MX Data\New Sample_20240331-001158\20240331...				
Exports	Status			
Raw Images	✓			
Colorized Images	✓			
Well Level CSV	✓			
Object Level ACS	✓			
Object Level CSV	✓			
DataSet	✓			
Reports	CSV	Excel	PDF	Word
PLX_Annexin V-FITC Hoechst P		✓		

Exporting and Printing Scan Results (Manual Settings)

To manually select additional export options (images/data) for scan result files as well as generated output files for reports, click the **Export** button located at the bottom of the Results screen. The Export dialog is displayed.

Confirm the export **Location** and if necessary update this path by clicking the **Browse** button, navigating to a folder on your Operating Computer or network, and clicking **OK**.

Select scan result **Images** (e.g., *Raw Images* or *Colorized Images*), **Data** (e.g., *Well Level CSV*, *Object Level CSV*, and *Object Level ACS*), and **Archive** (e.g., *Data Set*) file options to be exported to the specified location. *Data Sets* are stored as *.SCANRESULT* files.

In addition, for reports associated with the scan result you can select generated output file types (e.g., *CSV*, *Excel*, *PDF*, and *Word*) to be exported and indicate if files are to be opened automatically and/or printed upon export.

Note: If the *Object Level ACS* option is selected, the screen expands to display if an ACS template has been assigned to the assay and offers an *Auto Open* option after the report is generated. *If no ACS template is displayed, users can still export object level ACS data assuming they have the export privilege.* To change the ACS template assigned to an assay, users will need to edit the assay.

For report templates associated with the scan result you can select generated output file types (e.g., *CSV*, *Excel*, *PDF*, and *Word*) to be exported, and indicate if files are to be opened automatically and/or printed upon export.

Note: Any changes to export options are applied only when manually exporting scan results (i.e., by clicking the **Export and Print** button) and will *not* be saved with the assay or scan result.

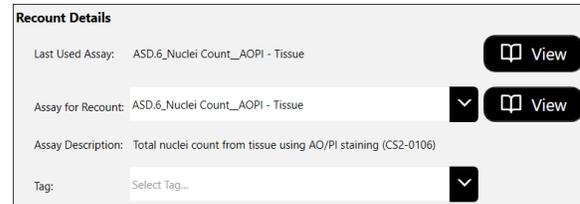
PERFORMING A RECOUNT

If you find it necessary to fine-tune assay parameters after reviewing your data results, click the **Recount** button located at the bottom of the Results screen. The Recount screen is displayed.

Once you have made any necessary changes (per the options described below), click the **Recount** button located at the bottom of the screen. After a recount is performed, the Navigation Bar returns to the Results screen.

Refining Assay Details/Selecting New Assay

In the *Recount Details* area you can view parameter settings for the *Last Used* assay, select a new assay from the dropdown or view the current assay to edit parameter settings to be used in the recount. See *Editing an Assay* on page 51 for details.



The screenshot shows the 'Recount Details' interface with the following fields and controls:

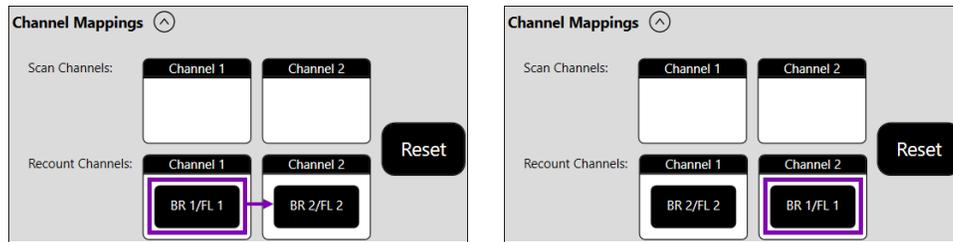
- Last Used Assay:** ASD.6_Nuclei Count_AOPI - Tissue (with a 'View' button)
- Assay for Recount:** ASD.6_Nuclei Count_AOPI - Tissue (with a dropdown arrow and a 'View' button)
- Assay Description:** Total nuclei count from tissue using AO/PI staining (CS2-0106)
- Tag:** Select Tag... (with a dropdown arrow)

Managing Channel Mappings

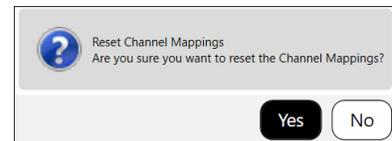
Expand the *Channel Mappings* area to identify *Scan Channels* and *Recount Channels* for the Imaging Mode associated with the current assay. Performing a recount uses channel mappings as defined by default (i.e., *Recount Channels* will remain in same positions used for the original *Scan Channels*).

Note: Each channel will have separate brightfield and fluorescent images as depicted in the mapping indicators (i.e., *BR1/FL1* and *BR2/FL2*).

To manage channel mappings for use in a recount, click on a mapping indicator and drag it to a new channel. The mapping indicator in that channel will swap positions with the one you are dragging automatically.

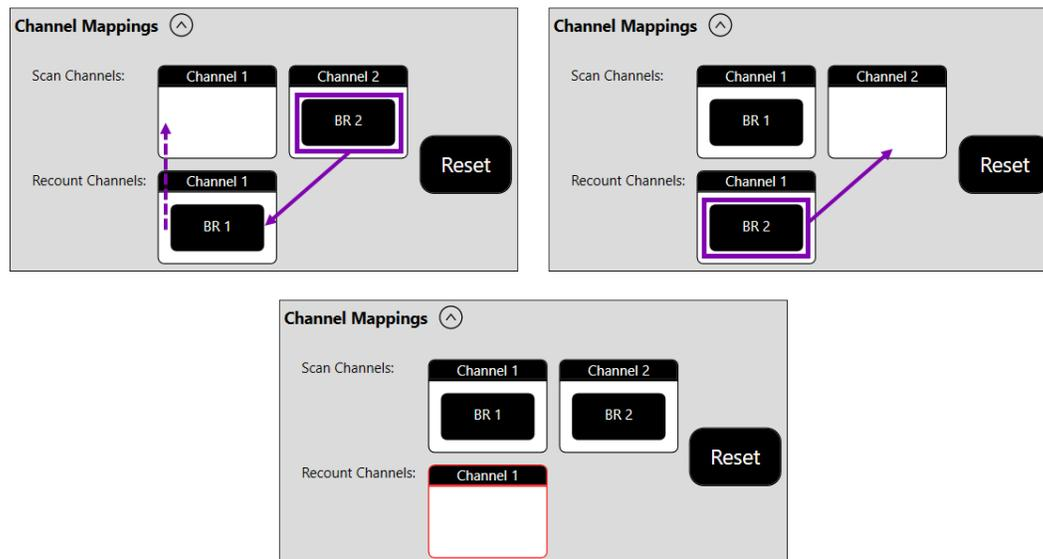


To return channel mappings to the positions they were in when you first expanded the *Channel Mappings* area, click the **Reset** button followed by **Yes** to confirm the action.



If a new assay selected for recount has fewer channels or you edit the assay previously used to reduce the number of channels, mappings displayed are updated to reflect available channels. To manage channel mappings for use in the recount, click on mapping indicators and drag them to an available position in the *Recount Channels* area.

2 Channel Assay Reduced to 1 Channel for Recount



If you drag indicators from *Recount Channels* to their home *Scan Channels* locations, a red outline is used to highlight empty *Recount Channels* locations. All *Recount Channels* must contain an indicator prior to recount.

Modifying Auto Export Options

Expand the *Reports and Exports* area to identify a **Location** for automatic exports of images/data and generated output files for reports. Click the **Browse** button and navigate to a folder on your Operating Computer or network to define the default export path. *This path will remain as the default in the software until it is manually changed.*

Exports	Will Be Exported	Reports	CSV	Excel	PDF	Word
Raw Images	<input checked="" type="checkbox"/>	MX_Display_FL Expression	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
Colorized Images	<input checked="" type="checkbox"/>		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
Well Level CSV	<input checked="" type="checkbox"/>		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
Object Level ACS	<input checked="" type="checkbox"/>		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
Object Level CSV	<input checked="" type="checkbox"/>		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>
DataSet	<input checked="" type="checkbox"/>		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>

In addition, *Exports* (e.g., *Raw Images*, *Colorized Images*, *Well Level CSV*, *Object Level ACS*, *Object Level CSV*, and *Dataset*) and output file types for *Reports* (e.g., *CSV*, *Excel*, *PDF*, and *Word*) defined for the assay will be indicated with a blue checkmark.

Note: To modify *Report and Export* settings displayed, click the **View** button for the selected assay in *Setup Details* area and expand the *Reports and Exports* section to update *Exports* and *Reports* file types selected. See *Managing Assay Reports and Exports* on page 57 for details.

Clicking the Recount Button

Once you have completed entering recount details for the samples, click the **Recount** button located at the bottom of the Recount screen. The Matrix software performs a recount of the scan using the modified parameters and displays the new scan result.

Chapter 5. Customizing Scan Result Reports

The Matrix custom reporting feature allows you to assign and/or modify report templates to be used by assays when generating scan results in the following ways:

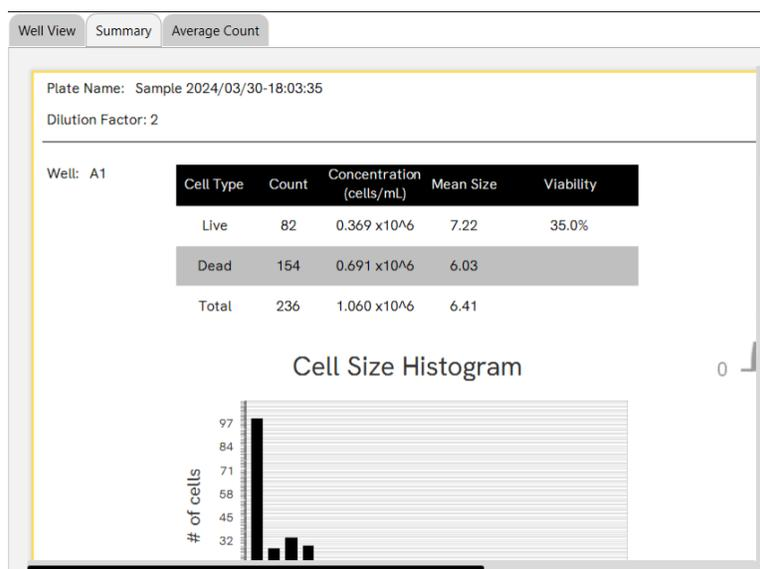
- The display template used for the **Well View** report tab may be changed. *This tab will always be displayed.*
- Report templates can be enabled/disabled as tabs and arranged to appear in sequence as desired.
- Report templates can be used to output specific file types (e.g., *CSV, Excel, PDF, and Word*) and open/print upon report generation.

Note: Any changes made to *Reporting* options for an assay will be saved with the assay and applied to all other scan results using that assay the next time a new scan or recount is performed.

VIEWING CUSTOM REPORTS

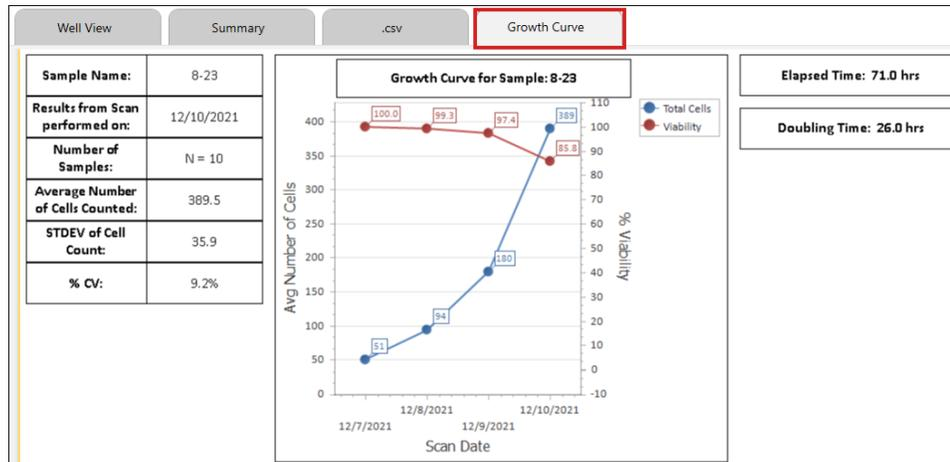
Although the **Well View** tab will always be enabled for display, you can change the report template assigned to this tab. In addition, you can disable any other default report tabs (e.g., **Summary** and **.csv** tabs for Matrix v4.0+; **All Wells** tab for v3.0), change the report templates used for display of these tabs, add new tabs and/or view a report template to edit its format directly.

For example, clicking the **Summary** tab for a scan result displays a full page view containing report data for all selected wells. *The Summary tab is enabled for Matrix v4.0+ assays by default and the assigned report template contains information similar to what was displayed in the All Wells tab available in earlier releases.*



You can either change the report template used for display of this tab or view it to modify the overall format of how report objects and data selections are presented. Any changes made to report templates will be saved with the assay and applied to all other scan results using that assay the next time a new scan or recount is performed.

As another example, a custom **Growth Curve** tab has been enabled for display in the scan result shown below. Clicking this tab displays consolidated sample data and line graphs for each group of wells illustrating the calculation of a growth curve for a time course series of scan results. *This report was generated on a Cellaca MX.*



Note: The Growth Curve shown above was created by assigning the *MX5_Growth Curve* report template (included as a custom report in the Matrix Report Templates Library) to an assay that was run repeatedly on the same sample in a time course series with a single tag applied to multiple scan results. See *Creating a Time Course Series* on page 35 for details on how create this custom report.

SETTING REPORTS AND EXPORTS FOR AN ASSAY

You can manage *Reports and Exports* options for an assay in the following areas of the Matrix software:

- When entering *Setup Details* in preparation for a count (i.e., **Acquire** tab Setup screen), select an assay and click the **View** button to edit assay details.
- When entering *Recount Details* for a selected scan result in preparation for a recount (i.e., **Data** tab Recount screen), select an assay to be used for the recount and click the **View** button to edit assay details.
- When managing details for a selected assay (**Manage** tab > Assays option) in the Assays List. Double-click on an assay in the list or highlight it once and then click the **View** button to edit assay details.

Once you are viewing details for an assay, expand the *Reports and Exports* area by clicking the  down arrow to view available options. When editing of the assay is complete, click **Save and Back** to save your changes to the assay or **Save As** to save a copy of the assay that includes your changes with a new name.

Note: As of the Matrix v4.0 release, report templates are no longer limited to the *Display*, *Export*, or *Print* template types used in earlier releases. A report template can be created for all these purposes and multiple templates assigned to an assay as needed. In addition, report tabs enabled by default will vary based on the Matrix software version associated with an assay (e.g., **Well View**, **Summary**, and **.csv** tabs are displayed for v4.0+ assays; **Well View** and **All Wells** tabs for v3.0 assays). Users can change report templates assigned to these tabs, edit the templates directly (i.e., any changes made will also be applied to other scan results using these templates), disable them from displaying as tabs or delete them from the assay.

Changing Template used in Well View Report Display

The **Well View** tab is enabled by default for all assays and cannot be disabled. However, you can change the report template used for the display of this tab or view the assigned report template and edit its format as necessary.

To change the report template currently assigned for the display of this tab, select a new template from the **Display** dropdown. Click **View** to confirm the selected template suits display needs for the report and edit it if necessary.

The screenshot shows a panel titled "Reports and Exports" with a dropdown menu labeled "Display" containing the text "MX5_Display_FL Expression_Well View". To the right of the dropdown is a blue button labeled "View".

Editing Assay Exports

Current assay *Exports* for scan result **Images** (e.g., *Raw Images* or *Colorized Images*), **Data** (e.g., *Well Level CSV*, *Object Level CSV*, and *Object Level ACS*), and **Archive** (e.g., *Data Set*) file options to be exported can be edited as necessary. *Data Sets are stored as .SCANRESULT files.*

If *Object Level ACS* is selected, the screen is expanded to display if an ACS template has been assigned and offers an *Auto Open* option. *If no template is displayed, users can still export object level ACS data assuming they have the export privilege.* To change the ACS template assigned to the assay, select a new one from the dropdown.

The screenshot shows an "Exports" panel with three sections: "Images" (with checkboxes for "Raw Images" and "Colorized Images"), "Data" (with checkboxes for "Well Level CSV", "Object Level CSV", and "Object Level ACS"), and "Archive" (with a checkbox for "Data Set").

The screenshot shows a "Data" panel with checkboxes for "Well Level CSV", "Object Level CSV", and "Object Level ACS". Below these is an "Object Level ACS Options" section with a checked "Use Template" checkbox and a dropdown menu labeled "Select An ACS Template To Use With the Export...". To the right is an "Auto Open" checkbox.

Managing the Assay Reports List

The Reports List for an assay may initially contain two report templates enabled by default used in the display of the **Summary** and **.csv** tabs. *Note that the .csv tab may not be available for assays using Expression imaging mode.*

The screenshot shows a "Reports" table with columns for Report Template, Display (Tab Name), CSV (Export, Auto Open, Print), Excel (Export, Auto Open, Print), PDF (Export, Auto Open, Print), and Word (Export, Auto Open, Print). Two rows are visible: "MX5_Display_BR and FL Viability" with Summary tab and "Default_Export_BR and FL Viability" with .csv tab. The "View" button above the table is highlighted with a red box.

Report Template	Display Tab Name	CSV			Excel			PDF			Word		
		Export	Auto Open	Print									
MX5_Display_BR and FL Viability	Summary	✓	✓										
Default_Export_BR and FL Viability	.csv	✓	✓										

To select a report template and edit how it is being applied to this assay, you can either double-click it in this list or click on the report once to highlight it and then click the **View** button.

You can use *Report Tab Options* to disable/enable the display of report templates as tabs and edit/enter tab names. *A total of five (5) custom tabs may be enabled for display at any one time.* In addition, you can select *Output File Options* for report templates and indicate if files are to be opened automatically or printed upon report generation.

When you have finished defining options for a report, click the **Update Report** button to save your changes and view the updated Reports List. Repeat this process for each report template to be modified.

To delete a report template from the Reports List for an assay, click on the report once to highlight it and then click the **Delete** button. You can also add a new report template by clicking the **Create** button. See *Adding Report Templates*, below for details.

Report Template	Display	CSV			Excel			PDF			Word			
		Tab Name	Export	Auto Open	Print									
MX4_Display_2FL Viability	Summary		✓			✓								
Default_Export_2FL Viability	Data					✓								
MX4_Growth Curve	Growth							✓	✓		✓	✓		
Default_Print_2FL Viability								✓		✓				

Note: If the report template deleted was from the Report Template Library provided with the Matrix software, it can be re-imported if necessary. If the report template was from a custom library or created using the Matrix software, it may be permanently deleted unless it was exported to an external location and saved prior to deletion.

As report tabs enabled for display are presented in the order in which they appear in the Reports List, you can change the sequence of tabs by highlighting a report and clicking the **Move Up** or **Move Down** buttons.

Adding Report Templates

To add a new report template, click the **Create** button. Select a template in the **Report Template** field, choose if the report should be enabled as a tab for display (i.e., **Yes**) and enter a tab name.

Select any output file types for the report and indicate if they are to be opened automatically (*Auto Open*) or printed upon report generation whenever the assay is used to perform a count/recount. *Output of files is independent of whether a template is also enabled as a tab.* Click **Add Report** to add the report to the Reports List.

CREATING A TIME COURSE SERIES

Creating a time course series of scan count results is useful when the intent of an experiment is to analyze the evolution of cell samples over time. The Matrix software custom reporting feature uses tags to logically group scan results in the order they are generated to identify them as part of a sequence. Tags can later be modified when viewing scan results in the Results List.

In this example, a time series of count results was acquired over the course of four days. For the first scan, a plate name of *Day 1* was entered for cell samples, an appropriate assay (e.g., *MX505.0_AOPI_Cell Lines*) was selected and the tag *Growth* added. *Once a tag is added you will be able to select it from the **Tag** dropdown in future scans.*

The screenshot shows the 'Setup Details' and 'Reports and Exports' sections of the Matrix software interface. In the 'Setup Details' section, the 'Plate Name' is 'Day 1', 'Select Assay' is 'MX505.0_AOPI_Cell Lines', 'Assay Description' is 'Total cell count and % viability using AOPI staining', 'Tag' is 'Growth', and 'Dilution Factor' is '2'. The 'Reports and Exports' section shows the 'Display' dropdown set to 'MX5_Display_2FL Viability_Well View'. The 'Reports' section is expanded, showing 'Report Template' as 'MX5_Growth Curve', 'Display in Tab' as 'No', and 'Tab Name' as 'Growth Curve'. A 'Reports' table is also visible, listing report templates and their associated display and export options.

Report Template	Display	CSV
	Tab Name	Export
MX5_Display_2FL Viability	Summary	
Default_Export_2FL Viability	.csv	
MX5_Growth Curve	Growth Curve	✓

The **View** button for the assay was selected to present assay details and the *Reporting* area expanded to create a **Growth Curve** tab (i.e., by assigning the *MX5_Growth Curve* report template) enabled for display.

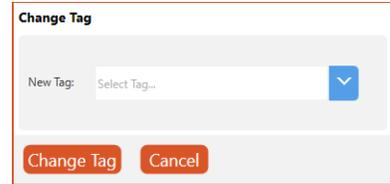
As the experiment progressed and subsequent counts were performed at determined intervals, new plate names were entered to reflect the stage of the experiment (e.g., *Day 2*, *Day 3*, *Day 4*) with users continuing to select the same tag (e.g., *Growth*) from the dropdown.

When the fourth day of the experiment was complete, scan results appeared in the Results List as shown below.

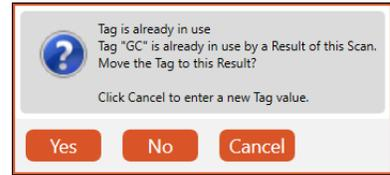
The screenshot shows the 'Results' list in the Matrix software interface. The 'Results' section includes buttons for Import, Export, Delete, Rename, and Tags. The 'Results' table shows four rows of scan results for 'Day 1' through 'Day 4', all with the 'Growth' tag. The 'Tag' column is highlighted with a red box.

Plate Name	Plate Type	Assay	Imaging Mode	Tag	Scan Creation	Result Creation	Magnification	Product Type
Day 4	12x2 Nexcelom Plate (CHM24-A100)	MX405.0_AOPI_Cell Lines	2 Brightfield and 2 Fluorescent	Growth	12/10/21 12:51:30	12/16/21 20:09:15	2.4	Mx
Day 3	12x2 Nexcelom Plate (CHM24-A100)	MX405.0_AOPI_Cell Lines	2 Brightfield and 2 Fluorescent	Growth	12/09/21 12:57:51	12/16/21 20:03:26	2.4	Mx
Day 2	12x2 Nexcelom Plate (CHM24-A100)	MX405.0_AOPI_Cell Lines	2 Brightfield and 2 Fluorescent	Growth	12/08/21 13:03:46	12/16/21 19:53:46	2.4	Mx
Day 1	12x2 Nexcelom Plate (CHM24-A100)	MX405.0_AOPI_Cell Lines	2 Brightfield and 2 Fluorescent	Growth	12/07/21 13:09:29	12/16/21 19:51:35	2.4	Mx

Once a scan result is available in the Results List, users can highlight it and click the **Tags** button to change the tag (i.e., add a new tag, edit/replace an existing tag or remove the tag by clearing the field). *Scan results associated with the same tag will be used collectively to create a time series.*



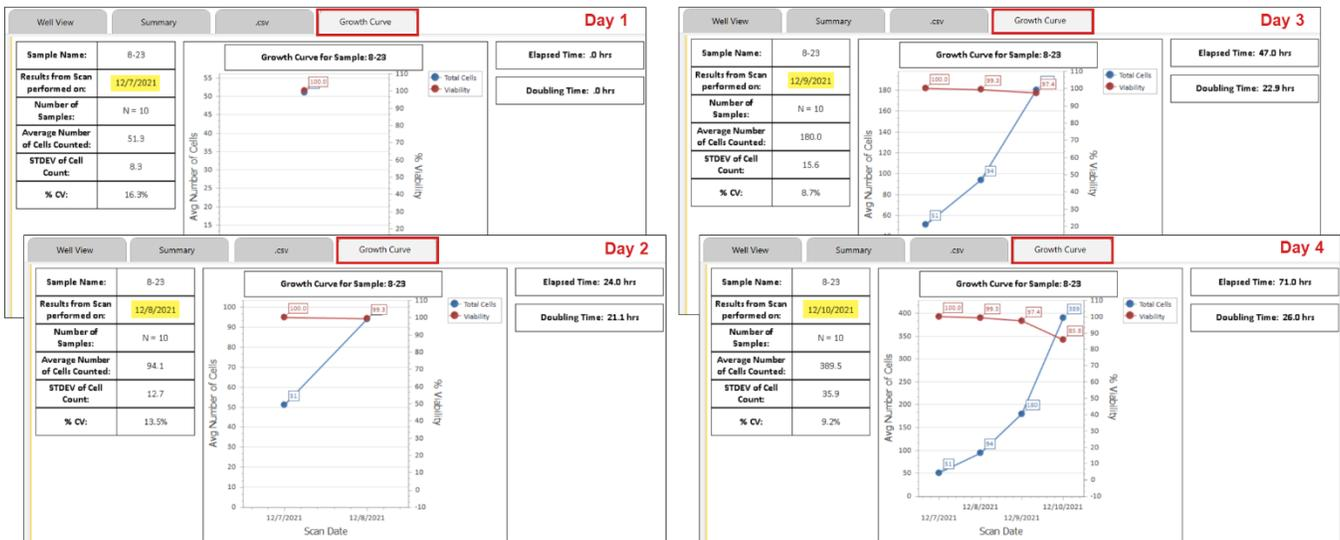
If users find it necessary to refine assay parameters and perform a recount for a scan result associated with a tag, they will be presented with a notification prompt that the tag is already in use by a result of the scan. Users can choose to move the tag to the new scan result (**Yes**), create the new scan result without the tag (**No**), or cancel the recount to select a new tag.



Note: Although a single tag may be applied to multiple scan results, only one result derived from each scan may be associated with the tag.

Another key concept when creating a time course series is that well names used in scan results must be consistent throughout the entire series and if multiple wells on the same plate use the same name, data from those wells will be averaged together before generation of the report. *Wells can be renamed while viewing a scan result.*

To create a **Growth Curve** report for this example containing all four scan results, users must select the scan result representing the *end* of the time series to include all data points (i.e., if users viewed the **Growth Curve** report generated for the scan result with plate names *Day 1*, *Day 2*, or *Day 3* as shown in the series as below, only the data points available at each specific point in time for the series will be included in the line chart).



Details for creating a simple custom time course report template using the DevExpress *Report Designer for WPF* plugin are provided in *Creating a Time Course Series* on page 35.

For assistance with creating or editing more complex time course time course reporting templates, contact Support by visiting <https://www.revvy.com/contact-us> or send email to: CellC-support@revvy.com

Best Practices and Workflow Tips

- Time course data is accumulated based on the well names used in scan results. If wells are *not* consistently named throughout the entire series, well data presented in the report may exceed the number of wells on the plate and/or include gaps. In addition, if the same well name applies to multiple wells on a plate, data in those wells will be averaged together before generation of the report. *Wells can easily be renamed after a scan and a recount performed to update the scan result.*
- When performing a recount on a scan result currently tagged to be part of a time course series, you will be prompted with a *Tag Already in Use* notification pop-up since the updated result will be derived from the same initial scan. Click **Yes** to move the tag to the updated result.
- Tags may consist of up to 32 characters in length and are displayed in a column in the Results List. *Click on the Tag column header to sort scan results in the list by tags.* In addition, tags may be used to filter the Results List to limit the display of scan results.
- If a tag is associated with a time course series, selecting the scan result with the most recent *Scan Creation* timestamp will include all data available for the report. *Selecting a scan result in the middle of the series will display only data available at that specific point in time.*
- To add a tag to be associated with multiple scan results, highlight all results before clicking the **Tags** button. *Click once to highlight the first result, and then hold either the **Shift** key (to select a block of results) or **Ctrl** key (to select non-contiguous results) down while clicking additional results.* Any changes you make to the tag will be applied to all selected scan results.
- To change a tag associated with multiple scan results that are assigned the same tag, highlight all results before clicking the **Tags** button. *Click once to highlight the first result, and then hold either the **Shift** key (to select a block of results) or **Ctrl** key (to select non-contiguous results) down while clicking additional results.* Any changes you make to the tag will be applied to all selected scan results.
- To delete tags associated with one or more scan results, highlight all results containing the tags to be deleted before clicking the **Tags** button. *Click once to highlight the first result, and then hold either the **Shift** key (to select a block of results) or **Ctrl** key (to select non-contiguous results) down while clicking additional results.* Ensure the **New Tag** field is blank and click **Change Tags** to remove the tags from all selected scan results.

CREATING A TIME COURSE REPORT TEMPLATE

Users can create a time course report template to illustrate data collected from a time course series by selecting the **Manage** tab > *Report Templates* option and clicking the **Create** button located at bottom of the screen. See *Editing a Report Template* on page 82 for details on using the *Report Template Designer*.

Follow these steps to create a simple time course report template containing a chart:

1. From the Control Toolbox displayed along the left side of viewing pane, select the **Chart**  object and drag it onto the template. The *Chart Designer* window opens automatically.
2. Click the **Change Chart Type**  icon located above the navigation panel on the left. Use the scroll bar to select the desired chart type (e.g., *Line* in the *Line Series* section).
3. In the navigation panel, click “Series 1” once to highlight it.
4. In the panel displayed on the right, click the **Data** tab. From the *objectDataSource1* list displayed, expand the *TimeCourseData* list (must be expanded for both levels).
5. Select and drag *ScanCreationTimeStamp* into the **Argument** cell below the list.
6. Select and drag *LiveCount* into the **Value** cell below the list.
7. Click **OK** followed by the **Select Result** button (located at bottom of the screen).
8. Choose a result that has a value in the *Tag* column and click **Select**. *If a tag is associated with multiple results, selecting the result with the most recent Scan Creation timestamp will include all the results in the chart while selecting a scan result in the middle of the series will display only data available at that point in time.*
9. Click the **Preview** tab (located in upper right corner of screen) to preview the chart with selected data.
10. Click **Save and Back** to enter a name for the time course report template.

Once you’ve created a time course report template, assign it to an assay being used for the collection of time series data and enable it for display as a report tab (e.g., **Growth Curve**). Ensure that tags have been added to scan results identifying key data points in the time course series and select a scan result associated with that tag to be used as the endpoint for the report.

To create a more complex time course report template or for assistance using *Report Template Designer*, contact Support by visiting <https://www.revivity.com/contact-us> or send email to: CellC-support@revivity.com

Chapter 6. Managing Favorites

This chapter describes how favorites are displayed in the *Favorites List* and how to manage the list.

A *Favorite* is a collection of count parameters such as an assay, consumable name/type, and reports/exports settings that can be saved for quick re-use. Users can assign images to be associated with favorites for quick reference, specify a prefix to be used for a series of consumables within an experiment, and enable the option of skipping the Preview screen (in favor of proceeding directly to the Count screen).

VIEWING THE FAVORITES LIST

You can manage favorites available to your instrument system by clicking the **Manage** tab > *Favorites* option to view the *Favorites List*. From this screen you can import/export, rename and delete favorites appearing in the list, as well as show/hide favorites displayed in the **Favorites Selection** panel in the **Acquire** tab Setup screen.

Name	Description	Assay Name	Product	Consumable Type	Skip Preview	Tag	State	Shown
AOPI Immune_MX	Concentration and viability measurement using AOPI on primar	MX_6_Viab_AOPI_Primary Cells	MX	12x2 Plate (CHM24-A100)	✓		🔒	✓
GFP Percent_MX	Concentration and GFP population percentages	MX_6_FL_Proteins_GFP Transfection Rate_	MX	12x2 Plate (CHM24-A100)	✓		🔒	✓
Nuclei AOPI_MX	Concentration of nuclei using AOPI for single-cell seq on tissue	MX_6_Nuclei Count_AOPI - Tissue	MX	12x2 Plate (CHM24-A100)	✓		🔒	✓

Use the **Per Page** control to change the number of favorites displayed per page, and the arrows to move back and forth between pages in the list.

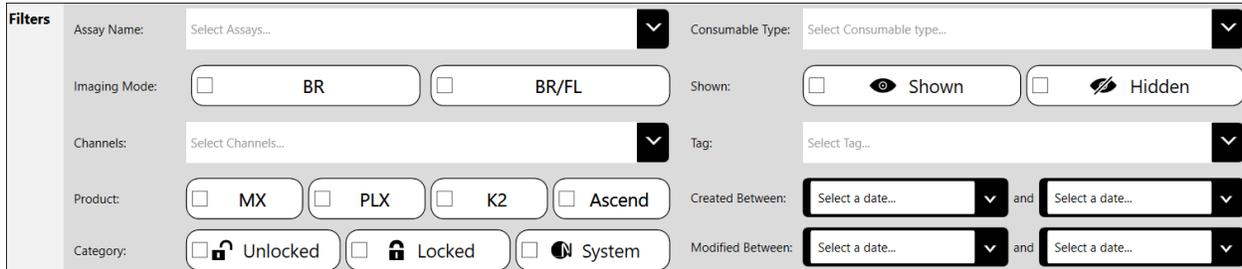
Favorites can be sorted by clicking column headings (i.e., ascending/descending indicators will be displayed), and information presented includes template name, description, assay name, product, consumable type, and tag.

Icons displayed in the *State* column indicate whether a favorite was provided as a Revvity *System* standard (which cannot be edited), or if a favorite is currently *Locked* or *Unlocked*. When viewing *System* and *Locked* favorites, a **Save As** button allows users to copy defined parameter settings as a source for creating a new favorite.

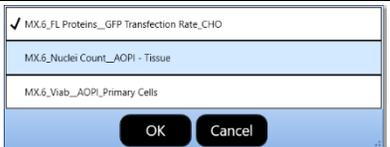
Checkmarks displayed in the *Skip Preview* column indicates the feature has been enabled and in the *Shown* column indicates the favorite is displayed in the **Favorite Selection** panel.

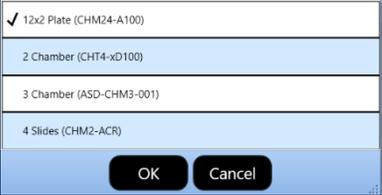
SEARCHING FOR FAVORITES

To search for a favorite, enter a few key characters of a favorite name or description in the **Search** field and/or expand the **Filters** area by clicking the down arrow  to specify search criteria. *Depending on monitor display size, you may need to collapse the Filters area to view search results.*



As you choose filters (e.g., by selecting *Assay Name*, *Imaging Mode*, *Channels*, *Product*, *Category*, *Consumable Type*, *Shown*, or *Tag* options) or enter a *Created Between/Modified Between* range of dates, the *Favorites List* is updated automatically to display matching entries.

Assay Name	Enter the first few characters of a favorite and then click the dropdown to select from matching entries. Click OK .	
Imaging Mode	Choose from the following imaging modes: BR – Filters for assays with Brightfield BR/FL – Filters for assays with Brightfield and Fluorescence	
Channels	Use the dropdown to select one or more channels to be included in the search, then click OK .	
Product	Choose from the following products: MX – Filters for favorites selected to be used for the Cellaca MX PLX – Filters for favorites selected to be used for the Cellaca PLX K2 – Filters for favorites selected to be used for the Cellometer K2 Ascend – Filters for favorites selected to be used for the Cellometer Ascend	
Category	Choose from the following categories: Unlocked – Filters for favorites that are unlocked and can be edited Locked – Filters for favorites that are locked and cannot be edited System – Filters for favorites provided by Revvity (locked and cannot be edited)	

Consumable Type	Enter the first few characters of a consumable type for an instrument and then click OK .	
Shown	<p>Choose from the following <i>Shown</i> states:</p> <p>Shown – Filters for favorites selected to appear in Favorites Selection panel</p> <p>Hidden – Filters for favorites that do <i>not</i> appear in Favorites Selection panel</p> <p>The Favorites Selection panel is available in Acquire tab Setup screen.</p>	
Tag	Enter the first few characters of favorites tag and then click the dropdown to select from matching entries. Click OK .	
Created Between	Enter a <i>Created Between</i> range indicating the start/end dates between which to filter for favorites created in that time frame.	
Modified Between	Enter a <i>Modified Between</i> range indicating the start/end dates between which to filter for favorites modified in that time frame.	

If you find it necessary to clear all selected filters and begin the search again, click the **Reset** button.

MAINTAINING THE FAVORITES LIST

When you select a favorite, buttons at the top of the *Favorites List* become available to perform the following functions. *The **Import** button will always be enabled as it does not require the selection of a favorite.*

Note: If the *Matrix 21 CFR Part 11* module is enabled, buttons will only be available for users who have been granted permission to perform specific functions.

Importing Favorites

1. While viewing the *Favorites List*, click the **Import** button.
2. Navigate to a folder where a favorite was previously saved.
3. Select one or more *.favorite* files to be imported.
4. Click **Open**. If selected favorites already exist in your database, respond to the confirmation prompt by clicking **Overwrite** to overwrite the file in your system, **Auto-Rename** to automatically add “(#)” to the end of the file name indicating the imported file is a copy (where # represents a value of 1, 2, 3, etc.) or **Cancel** to abort the import.
5. Click **OK** to acknowledge the successful import and confirm that imported favorites are displayed in the *Favorites List*.

Exporting Favorites

1. Select one or more favorites from the Favorites *List* to be exported by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of favorites) or **Ctrl** key (to select non-contiguous favorites) down while clicking additional favorites.
2. Click the **Export** button.
3. Navigate to a folder where favorites are to be saved.
4. Click **OK** to save *.favorite* files in the export location.

Renaming Favorites

1. Select the favorite to be renamed and click the **Rename** button.
2. Edit the favorite name and click **Rename** to save your changes.

Deleting Favorites

1. Select one or more favorites from the Favorites to be deleted by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of favorites) or **Ctrl** key (to select non-contiguous favorites) down while clicking on additional favorites.

Click the **Delete** button followed by **Yes** to confirm the action.

If the Matrix 21 CFR Part 11 module is enabled, users will be prompted to enter a reason prior to deletion.

Note: If favorite deleted was from the Favorites Library provided with the Matrix software, it can be re-imported if necessary. If favorite was from a custom library, it may be permanently deleted unless favorite was exported to an external location and saved prior to deletion.

Showing/Hiding Favorites in Selection Panel

*The **Favorites Selection** panel is available in the Acquire tab Setup screen.*

1. Select one or more favorites from the Favorites *List* to be shown or hidden in the **Favorites Selection** panel by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of favorites) or **Ctrl** key (to select non-contiguous favorites) down while clicking on additional favorites.
2. Click the **Show** or **Hide** button depending on whether you want to show/hide the selected favorite in the **Favorites Selection** panel.

EDITING A FAVORITE

Editing a favorite allows users to change basic *Favorite Details* (such as name, description, icon, category, consumable ID, assay, etc. as described below). *Not all options shown in this section are available for all products and instrument configurations.*

To select a favorite and view its details, double-click it in the *Favorites List*, or click the assay once to highlight it and then click the **View** button located at the bottom of the screen.

Note: The favorite displayed contains many sections in which the content of each section may vary depending on the product type and the channels chosen to be associated with the assay.

Edit the assay as described in each of the sections presented below:

- *Modifying Favorite Details* on page 44
- *Choosing a Consumable Type* on page 45
- *Managing Favorite Reports and Exports* on page 45

When viewing/editing of a favorite is complete, click one of the following buttons to return to the *Favorites List*.



Back Button: Click **Back** to return to previous screen without saving any changes. *If favorite has unsaved changes, click **OK** in response to confirmation prompt.*



Save Button: Click **Save** to save your changes and return to previous screen. Click **Auto Back** to toggle functionality of the **Save** button between **Save** (to save and continue editing) and **Save and Back** (to save and return to previous screen).



Save As Button: Click **Save As** to save your changes as a copy with a new name or select another favorite from the dropdown to override it and return to the previous screen. *If the favorite is locked, this also allows you to copy the favorite to use as a source for creating a new favorite.*

Modifying Favorite Details

In the *Favorite Details* area, you can edit the favorite *Name*, *Description*, *Icon*, *Category*, *Consumable ID*, *Assay*, *Tag*, or *Skip Preview* feature as indicated below.

The screenshot shows the 'Favorite Details' form with the following fields and controls:

- Name:** Text input field containing 'AOPI Immune_ASD'.
- Description:** Text input field containing 'AOPI Immune cell assay using 3-chamber slide with slide autofocus'.
- Icon:** Image preview showing a yellow 3D model of a cell, with 'Browse...' and 'Clear' buttons.
- Category:** Radio buttons for 'Unlocked' (checked) and 'Locked'.
- Consumable ID:** Text input field containing 'Sample', with an 'Add Timestamp' checkbox.
- Assay:** Dropdown menu showing 'ASD.6_Viab_AOPI_Immune Cells, Low RBC' and a 'View' button.
- Assay Description:** Text input field containing 'Total cell count and % viability using AO/PI staining (CS2-0106)'.
- Tag:** Text input field with placeholder 'Type here to enter a Tag...'.
- Skip Preview:** Radio buttons for 'No' and 'Yes'.

Note: You can only edit a favorite if the category displayed is *Unlocked*. If the category displayed is *Locked*, you must first copy the favorite by clicking the **Save As** button and entering a new name to save defined parameters as a source for creating a new assay. Once saved, you can edit copied parameters in the new assay as necessary.

To edit favorite details, modify information contained in the following fields.

- Name** Displays the name of the favorite. *Must be unique*.
- Description** Displays a brief description which can be used to identify the purpose of the favorite or include defined parameters to help users distinguish it from others presented in the Favorites *List*.
- Icon** Displays an image which can be used to quickly identify the favorite and help users distinguish it from others presented in the Favorite Selection panel. Click the **Browse** button to select a new image or the **Clear** button to remove the current image.
- Category** Indicates if the current status of a favorite is *Unlocked* or *Locked*. When viewing details for a *Locked* or *System* assay, use the **Save As** button to enter a new name and copy defined parameters to be used as a source for creating a new assay.

The locked state of an assay is displayed in the *State* column of the *Favorites List* using the Revvity *System*, *Unlocked*, or *Locked* icons.

If an assay is unlocked, you can select the **Locked**  button to prevent the assay from being edited by other users.

Note: It is recommended that you do *not* select the **Locked** button until *after* editing is complete as once the assay is locked, you will no longer be able to make any changes.

- Consumable ID** Displays the consumable ID of the favorite. Click the Add Timestamp button to enable/disable automatically appending a timestamp to the consumable ID.
- Assay** Displays the assay associated with the favorite. Click the dropdown to change the selected assay.
- Tag** Displays any tags assigned to the favorite.
- Skip Preview** Toggles the display of the **Preview** (if set to **No**) and **Count** (if set to **Yes**) buttons.

Choosing a Consumable Type

In the *Consumable* area, click the Plate Type dropdown to display consumable types available for the instrument.

If the **Images Per Well** option is available, click the dropdown and select an option.

A preview of the consumable type is displayed.

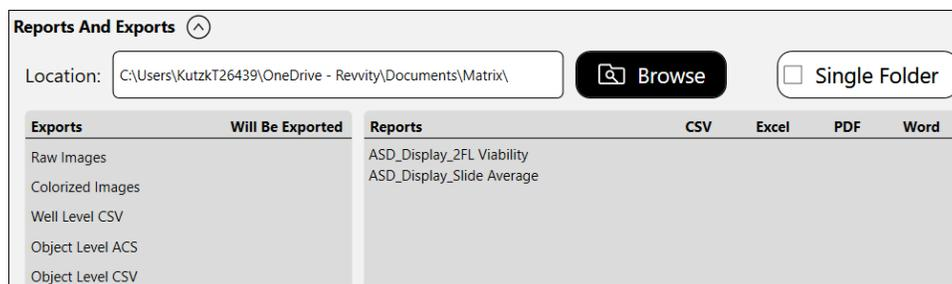


Managing Favorite Reports and Exports

In the *Reports and Exports* area you can assign report templates to control how data is displayed, exported and printed. In addition to the **Well View** tab which will always be displayed, a total of five (5) custom reporting tabs may be enabled for display at any one time.

Note: As of the Matrix v4.0 release, report templates are no longer limited to the *Display, Export, or Print* template types used in earlier releases. A report template can be created for all these purposes and multiple templates assigned to an assay as needed. In addition, report tabs enabled by default will vary based on the Matrix software version associated with an assay (e.g., **Well View, Summary** and **.csv** tabs are displayed for v4.0 assays; **Well View** and **All Wells** tabs for v3.0 assays). Users can change report templates assigned to these tabs, edit the templates directly (i.e., any changes made will also be applied to other scan results using these templates), disable them from displaying as tabs, or delete them from the assay.

If *Reports and Exports* parameters are not displayed, expand the area by clicking the down arrow to view available options.



Reports and Exports options are associated with the defined assay. See *Managing Assay Reports and Exports* on page 57 to manage these options for assay assigned to the favorite.

Creating a New Favorite

You can create a new favorite by clicking the **Manage** tab > *Favorites* option to display the *Favorites List* and then clicking the **Create** button located at the bottom of the screen.

The screenshot shows the 'Favorite Details' form within the 'Manage' tab. The form includes the following fields and controls:

- Name:** A text input field with the placeholder text 'Type here to enter a favorite name...'.
- Description:** A text input field with the placeholder text 'Type here to enter a favorite description...'.
- Icon:** A square icon placeholder, a 'Browse...' button, and a 'Clear' button.
- Category:** Two radio buttons labeled 'Unlocked' (checked) and 'Locked'.
- Consumable ID:** A text input field with the placeholder text 'Type here to enter a Consumable ID...' and a checked 'Add Timestamp' checkbox.
- Assay:** A dropdown menu with the text 'Select Assay...' and a 'View' button.
- Assay Description:** A text input field.

At the bottom of the screen, there is a navigation bar with a 'Back' button, an 'Auto Back' toggle (checked), and 'Save' and 'Save As' buttons.

Enter basic *Favorite Details* (such as name, description, icon, category, consumable ID, assay, and tag), choose if the **Skip Preview** feature is to be enabled, select a consumable type, specify a *Reports and Exports location*, and then click the **Save and Back** button to add the assay to the *Assays List*.

To copy defined parameter settings for an existing favorite to be used as a source for creating a new favorite, select a favorite and view its details by double-clicking it in the *Favorites List* or clicking it once to highlight it and then clicking the **View** button located at the bottom of the screen. Click the **Save As** button and enter a new name to save a copy of the favorite, and then click the **Save** button.

Edit favorite parameters as described in *Editing a Favorite* on page 43 and click the **Save** button (with **Auto Back** enabled) to confirm the new favorite has been saved to the *Favorites List*.

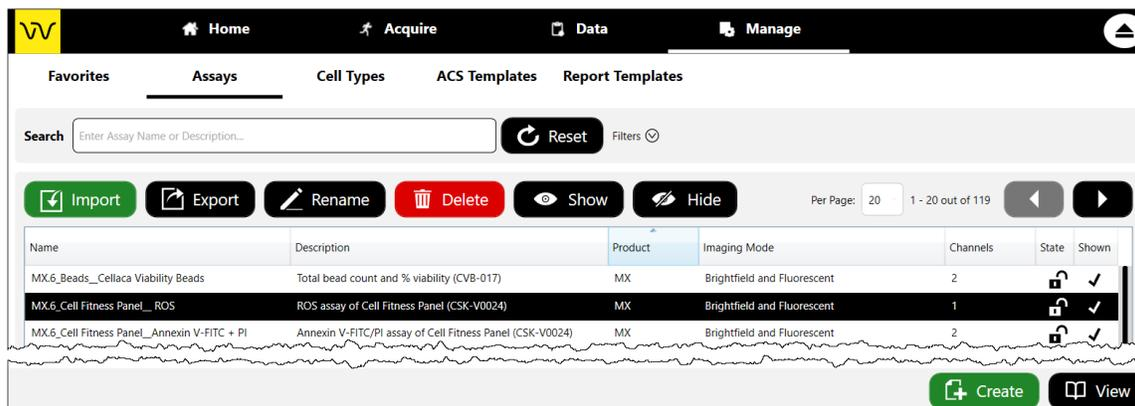
Chapter 7. Managing Assays

This chapter describes how assays are displayed in the *Assays List* and how to manage the list to keep it current. Details on how to edit and create assays are also provided. Assay parameters can be optimized before each count /recount to meet your analysis needs.

Note: When migrating to the Matrix database from an earlier version, updated assays can be imported via the software *after* migration is complete. For a list of assays provided with the Matrix software or for assistance in defining a custom assay, contact Support by visiting <https://www.revvy.com/contact-us> or send email to: CellC-support@revvity.com

VIEWING THE ASSAYS LIST

You can manage assays available for your instrument system by clicking the **Manage** tab > *Assays* option to view the *Assays List*. From this screen you can import/export, rename, and delete assays appearing in the *Assays List*, as well as show/hide assays displayed in the *Assays* dropdown appearing in other screens.



In addition, you can select an assay from the list to view/edit its details, create a new assay by defining custom parameter settings or use a locked assay as a source for creating a new assay based on its parameter settings.

Note: If the Matrix 21 *CFR Part 11* module is enabled, the **Create** button (and **Save and Back** button when viewing assay details) will only be available for users who have been granted permission to perform these functions.

Use the **Per Page** control to change the number of assays displayed per page, and the arrows to move back and forth between pages in the list.

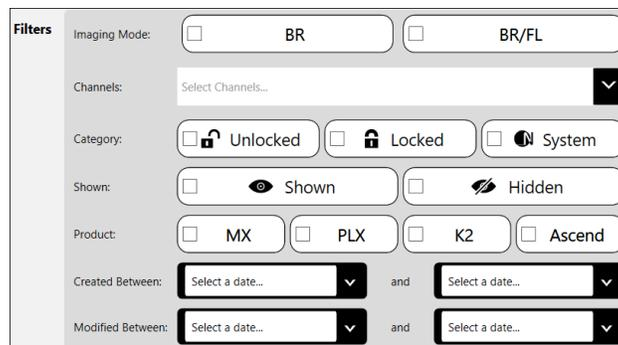
Assays can be sorted by clicking on column headings (i.e., ascending/descending indicators will be displayed), and information presented for each assay includes name, description, the product for which it is intended, imaging mode defined for the assay, and created/last modified dates.

Icons displayed in the *State* column indicate whether an assay was provided as a Revvity *System* standard (which cannot be edited), or if an assay is currently *Locked* or *Unlocked*. When viewing *System* and *Locked* assays, a **Save As** button allows users to copy defined parameter settings as a source for creating a new assay.

A checkmark ✓ displayed in the *Shown* column indicates the assay is to be included in the *Assays* dropdown available in other screens (e.g., **Select Assay** field in the **Acquire** tab Setup screen and **Assay for Recount** field in the **Acquire** tab *Recount* screen).

SEARCHING FOR ASSAYS

To search for an assay, enter a few key characters of an assay name or description in the **Search** field and/or expand the *Filters* area by clicking the down arrow  to specify search criteria. *Depending on monitor display size, you may need to collapse the Filters area to view search results.*



As you choose filters (e.g., by selecting *Imaging Mode*, *Category*, *Shown*, and *Product* options) or enter a *Created Between/Modified Between* range of dates, the *Assays List* is updated automatically to display matching entries.

Imaging Mode

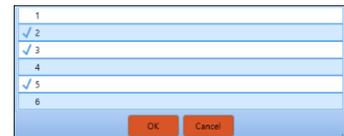
Choose from the following imaging modes:

BR – Filters for assays with Brightfield

BR/FL – Filters for assays with Brightfield and Fluorescence

Channels

Use the dropdown to select one or more channels to be included in the search, then click **OK**.



Category

Choose from the following categories:

Unlocked – Filters for assays that are unlocked and can be edited

Locked – Filters for assays that are locked and cannot be edited

System – Filters for assays provided by Revvity (locked and cannot be edited)

System assays are no longer available in the Assay Library provided with the Matrix software but may exist on your system if imported from an earlier release.

- Shown** Choose from the following *Shown* states:
- Shown** – Filters for assays selected to appear in the *Assays* dropdown
 - Hidden** – Filters for assays that do *not* appear in the *Assays* dropdown
- Note:** The *Assays* dropdown is available in other screens (e.g., **Select Assays** field in the Setup screen and **Assay for Recount** field in the Recount screen).
- Product** Choose from the following products:
- MX** – Filters for assays selected to be used for the Cellaca MX
 - PLX** – Filters for assays selected to be used for the Cellaca PLX
 - K2** – Filters for assays selected to be used for the Cellometer K2
 - Ascend** – Filters for assays selected to be used for the Cellometer Ascend
- Created Between** Enter a *Created Between* range by selecting start/end dates representing a time period during which to filter for assays that were created.
- Modified Between** Enter a *Modified Between* range by selecting start/end dates representing a time period during which to filter for assays that were modified.

If you find it necessary to clear all selected filters and begin the search again, click the **Reset** button.

To select an assay and view its details, double-click it in the *Assays List*, or click the assay once to highlight it and then click the **View** button located at the bottom of the screen.

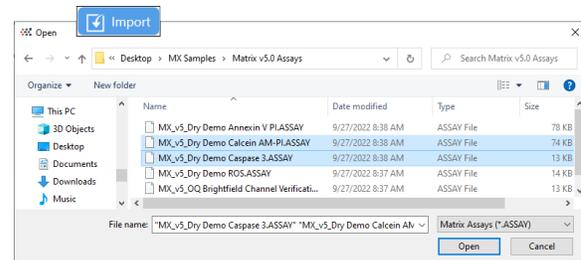
MAINTAINING THE ASSAYS LIST

When you select an assay, buttons at the top of the *Assays List* become available to perform the following functions. *The **Import** button will always be available as it does not require the selection of an assay in your library.*

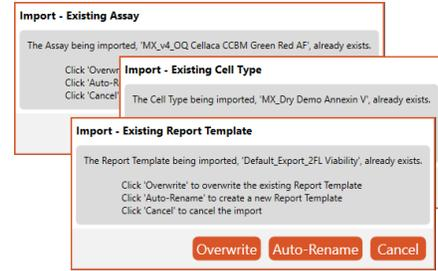
Note: If the *Matrix 21 CFR Part 11* module is enabled, buttons will only be available for users who have been granted permission to perform specific functions.

Importing Assays

1. While viewing the *Assays List*, click the **Import** button.
2. Navigate to a folder where an external assay library is available or an assay was previously saved.
3. Select one or more *.ASSAY* files to be imported. *Keep in mind that cell types and report templates associated with selected assays will also be included in the import.*

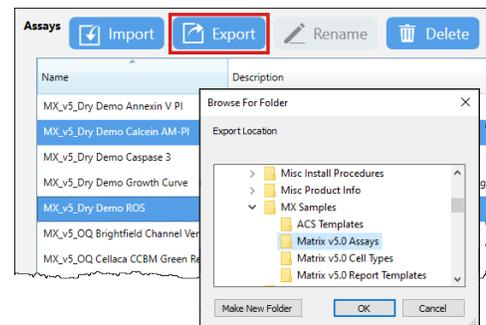


- Click **Open**. If selected assays (or cell types and report templates associated with the assays) already exist in your database, respond to confirmation prompts by overwriting the files in your system, auto renaming the files on import or canceling import of the file.
- Click **OK** to acknowledge the successful import and confirm that imported assays are displayed in the *Assays List*.



Exporting Assays

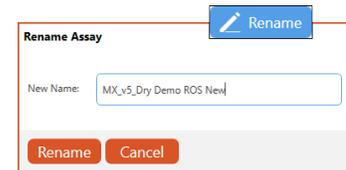
- Select one or more assays from the *Assays List* to be exported by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of assays) or **Ctrl** key (to select non-contiguous assays) down while clicking additional assays.
- Click the **Export** button.
- Navigate to a folder where assays are to be saved.
- Click **OK** to save .ASSAY files in the export location.



Renaming Assays

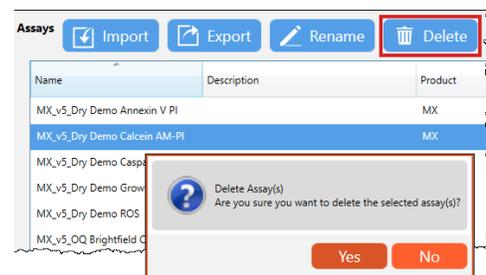
The **Rename** button will only be enabled for unlocked assays.

- Select the assay to be renamed and click the **Rename** button.
- Edit the assay name and click **Rename** to save your changes.



Deleting Assays

- Select one or more assays from the *Assays List* to be deleted by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of assays) or **Ctrl** key (to select non-contiguous assays) down while clicking on additional assays.
- Click the **Delete** button followed by **Yes** to confirm the action. *If the Matrix 21 CFR Part 11 module is enabled, users will be prompted to enter a reason prior to deletion.*



Note: If assays deleted were from the Assay Library provided with the Matrix software, they can be re-imported if necessary. If assays were from a custom library or created using the Matrix software, they may be permanently deleted unless assays were exported to an external location and saved prior to deletion.

Showing/Hiding Assays in Dropdown

The Assays dropdown is available in other screens (e.g., **Select Assay** field in the **Acquire** tab Setup screen and **Assay for Recount** field in the **Acquire** tab Recount screen).

1. Select one or more assays from the Assays List to be shown or hidden in the Assays dropdown by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of assays) or **Ctrl** key (to select non-contiguous assays) down while clicking on additional assays.
2. Click the **Show** or **Hide** button depending on whether you want to show/hide the selected assays in the Assays dropdown available in other screens.

Name	Description	Product	Imaging Mode	Channels	Created	Modified	State	Shown
MX_v5_Dry Demo Annexin V PI		MX	Brightfield and Fluorescer	2	8/28/2020 10:48:19 AM	9/27/2022 8:44:38 AM		<input checked="" type="checkbox"/>
MX_v5_Dry Demo Calcein AM-PI		MX	Brightfield and Fluorescer	2	8/28/2020 9:54:34 AM	9/22/2022 11:26:01 AM		<input checked="" type="checkbox"/>
MX_v5_Dry Demo Caspase 3		MX	Brightfield and Fluorescer	1	9/9/2020 10:14:49 AM	9/27/2022 8:44:31 AM		<input checked="" type="checkbox"/>
MX_v5_Dry Demo Growth Curve	Calculation of Doubling Time using scans with	MX	Brightfield and Fluorescer	2	9/17/2021 1:47:13 PM	9/27/2022 8:44:57 AM		<input checked="" type="checkbox"/>
MX_v5_Dry Demo ROS		MX	Brightfield and Fluorescer	1	8/28/2020 9:54:05 AM	9/27/2022 8:44:54 AM		<input checked="" type="checkbox"/>

EDITING AN ASSAY

Editing an assay allows users to change basic *Assay Details* (such as name, description, category, and product type), defined *Imaging and Analysis* parameter settings, and *Reporting* templates currently assigned to the assay. *Not all options shown in this section are available for all products and instrument configurations.*

To select an assay and view its details, double-click it in the Assays List, or click the assay once to highlight it and then click the **View** button located at the bottom of the screen.

Note: The assay displayed contains many sections in which the content of each section may vary depending on the product type and the channels chosen to be associated with the assay.

Edit the assay as described in each of the sections presented below:

- *Modifying Basic Assay Details* on page 52
- *Defining Assay Imaging and Analysis Parameters* on page 53
- *Managing Assay Reports and Exports* on page 57

When viewing/editing of an assay is complete, click one of the following buttons to return to the Assays List.



Back Button: Click **Back** to return to previous screen without saving any changes. *If favorite has unsaved changes, click **OK** in response to confirmation prompt.*



Save Button: Click **Save** to save your changes and return to previous screen. Click **Auto Back** to toggle functionality of the **Save** button between **Save** (to save and continue editing) and **Save and Back** (to save and return to previous screen).



Save As Button: Click **Save As** to save your changes as a copy with a new name or select another favorite from the dropdown to override it and return to the previous screen. *If the favorite is locked, this also allows you to copy the favorite to use as a source for creating a new favorite.* Click  to toggle functionality of the **Save As** button between **Save As** (to save as and continue editing) and **Save As and Back** (to save as and return to previous screen).

Modifying Basic Assay Details

In the *Assay Details* area, you can edit the assay *Name*, *Description*, *Category*, or *Product* as indicated below.

Note: You can only edit an assay if the category displayed is *Unlocked*. If the category displayed is *Locked* or if the assay is a Revvity *System* default, you must first copy the assay by clicking the **Save As** button and entering a new name to save defined parameters as a source for creating a new assay. Once saved, you can edit copied parameters in the new assay as necessary.

To edit basic assay details, modify information contained in the following fields.

Name Displays the name of the assay. *Must be unique.*

Description Displays a brief description which can be used to identify the purpose of the assay or include defined parameters to help users distinguish it from others presented in the *Assays List*.

Category Indicates if an assay was provided as a Revvity default   *System* (i.e., *System* assays cannot be edited) or if its current status is *Unlocked* or *Locked*. When viewing details for a *Locked* or *System* assay, use the **Save As** button to enter a new name and copy defined parameters to be used as a source for creating a new assay.

The locked state of an assay is displayed in the *State* column of the *Assays List* using the Revvity *System*, *Unlocked*, or *Locked* icons.

If an assay is unlocked, you can select the **Locked**  button to prevent the assay from being edited by other users.

Note: It is recommended that you do *not* select the **Locked** button until *after* editing is complete as once the assay is locked, you will no longer be able to make any changes.

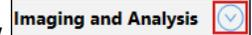
Product Indicates product type for which the assay was created (e.g., *MX*, *PLX*, *K2*).

Note: As the Matrix software is shared by multiple product families, it must account for physical differences between instruments. To accommodate for these differences, assay details available for editing may vary from what is presented in this guide based on the selected product.

Defining Assay Imaging and Analysis Parameters

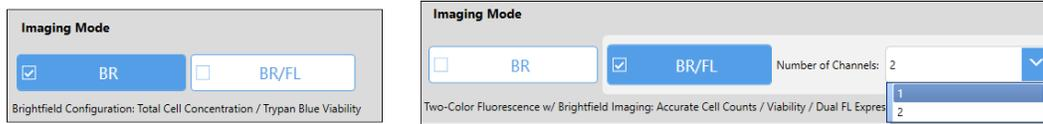
In the *Imaging and Analysis* area, you can edit defined assay parameter settings for *Imaging Mode*, *Analysis Mode*, *Focusing Mode* (if available based on product), *Dilution* and *Channel Imaging Parameters* as indicated below.

If *Imaging and Analysis* parameters are not displayed, expand the area by clicking the down arrow to view available options.



IMAGING MODE OPTIONS

Edit *Imaging Mode* selected for the assay by clicking an imaging mode option.



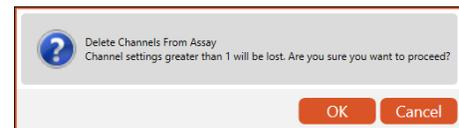
BR **Brightfield Only:** Indicates assay will use a brightfield image to determine *Total Cell Concentration* and *Trypan Blue Cell Viability*.

BR/FL **Brightfield/Fluorescence:** Indicates assay will use single or dual channel imaging by analyzing fluorescent images in conjunction with their associated brightfield images.

Number of Channels: Dropdown lists available channels for the assay (i.e., 1-2 for Cellometer Ascend/K2).

- *BR/FL 1* – Brightfield Image, Single Fluorescent Image Analysis
- *BR/FL 2* – Brightfield Image, Two Fluorescent Image Analysis

Note: If channels previously defined for assay will be lost as a result of a new imaging mode selected, you will be prompted with a confirmation message. Click **OK** to proceed or **Cancel** to retain current imaging mode.



ANALYSIS MODE OPTIONS

In the *Analysis Mode* area you can edit the analysis mode selected for the assay by clicking available options (as shown in the screens below) and entering values for any additional fields presented.

Note: Analysis modes displayed will vary based on the selected *Imaging Mode* for the assay. In addition, *Channel Imaging and Cell Type Parameters* displayed on page 55 will also vary as a result of the selected *Analysis Mode*.

FOR BR ONLY IMAGING MODE

Cell Count – Perform brightfield image analysis to determine *Total Cell Concentration*.

Trypan Viability – Perform brightfield image analysis of Trypan Blue exclusion to determine *Cell Viability*.

FOR BR/FL 1 CHANNEL IMAGING MODE

Cell Count – Count cells in brightfield/fluorescent images independently. Allows you to select either a **BR** or **FL** mask to assist in finding FL positive cells. *Use a fluorescent mask if cells are difficult to detect in brightfield images.*

Viability – Calculate cell concentration and % viability using a nuclear staining dye (single-fluorescence). Allows you to select a **BR**, **FL**, or **Hybrid** mask to assist in finding FL positive cells. In addition, you can select **FL Classification (Live or Dead)**.

Use a fluorescent mask if cells are difficult to detect in brightfield images OR a Hybrid mask when counting FL images to exclude corresponding cells found in BR images (e.g., dead cells with faint walls that only show up using FL).

Expression – Perform analysis using a single fluorophore (e.g., GFP, RFP, etc.). Allows you to select either a **BR** or **FL** mask to assist in finding FL positive cells. *Use a fluorescent mask if cells are difficult to detect in brightfield images.*

You can also expand (*increase **Expand** value*) or contract (*decrease **Expand** value*) the mask around objects found to accurately collect fluorescent intensity measurements.

Expression assays use a single mask for segmentation while pulling fluorescent values from the same area across all channels. By enabling the software to read FL values from cells without requiring those cells to be segmented in each FL channel using traditional methods, weaker FL intensities can be identified.

FOR BR/FL 2 CHANNEL IMAGING MODE

Viability – Calculate cell concentrations and % viability using two nuclear staining dyes (dual-fluorescence). Allows you to select either no mask (**None**) or a brightfield (**BR**) mask to assist in finding FL positive cells. You can also select a **FL Classification** (**Total**, **Live**, or **Dead populations**) for each channel.

Total classification can only be assigned to one FL channel. If currently selected in one channel, selecting it in the other channel automatically swaps the first channel options.

Analysis Mode
 Viability Expression
 Dual-Fluorescence Cell Viability: Report Concentrations And Viability Using Nuclear Staining Dyes
 Mask: None BR
 FL1 Classification: Total Live
 FL2 Classification: Total Dead

Expression – Perform analysis using dual fluorescence for samples with two FL stains. Allows you to select either a **BR** or **FL** mask to assist in finding FL positive cells. *Use a fluorescent mask if cells are difficult to detect in brightfield images.*

You can also expand (*increase **Expand** value*) or contract (*decrease **Expand** value*) the mask around objects found to accurately collect fluorescent intensity measurements.

Analysis Mode
 Viability Expression
 Dual Fluorescence Analysis For Samples Containing Two FL Stains
 Mask: BR FL
 Uses the Brightfield image to aid in the finding of FL positive Cells
 Expand (µm):
 Amount, in microns, to expand or contract the found mask object which is used to collect FL intensity measurements in all channels

Expression assays use a single mask for segmentation while pulling fluorescent values from the same area across all channels. By enabling the software to read FL values from cells without requiring those cells to be segmented in each FL channel using traditional methods, weaker FL intensities can be identified.

DILUTION FACTOR

In the *Dilution* area, you can edit the final dilution factor by clicking in the field and modifying the displayed value.

Dilution
 Dilution Factor For General Assay As Indicated By Sample Preparation Protocol

CHANNEL IMAGING AND CELL TYPE PARAMETERS

Channel Imaging Parameters are grouped using tabs and display will vary based on the selected *Imaging Mode* (i.e., *Channel 1* tab is displayed for **BR** or **BR/FL 1** imaging modes, and *Channel 2* tab is displayed for **BR/FL 2**).

Note: *Channel 1* refers to **BR** only or **BR1|FL1** images taken by the instrument's camera during a scan through the first filter set and *Channel 2* refers to the **BR2|FL2** images taken through the second filter set.

You can edit *Channel Imaging Parameters* (e.g., *Brightfield*, *Fluorescence* and *Filters*) for the assay. If **Channel 1** and **Channel 2** tabs are available, toggle between the display for each channel by clicking on the applicable tab. *Channel 1 Imaging Parameters are displayed by default.*

The screenshot displays two panels for channel configuration. The top panel is for Channel 1, with 'Channel 1' selected in the tab bar. It contains three sections: 'Brightfield' with 'Use Custom Exposure' set to 'No' and 'Custom Exposure Factor' at 1.0; 'Fluorescence' with 'Fluorophore Name' as 'AO' and 'Exposure (ms)' as 700; and 'Filters' with '535' selected and '660' unselected. The bottom panel is for Channel 2, with 'Channel 2' selected. It contains similar sections: 'Brightfield' with 'Use Custom Exposure' set to 'No' and 'Custom Exposure Factor' at 1.0; 'Fluorescence' with 'Fluorophore Name' as 'PI' and 'Exposure (ms)' as 2500; and 'Filters' with '535' unselected and '660' selected. Below the Channel 2 parameters is a dropdown menu showing 'K2_PI Stained Cells' and a 'View' button.

Brightfield **Use Custom Exposure** – Indicates if custom exposure factors are being used for the channel. Click **Yes** to edit value in the **Custom Exposure Factor** field (below).

Custom Exposure Factor – Value prolongs or shortens exposure time for brightfield image.

Fluorescence **Fluorophore Name** – Displays fluorophore name being used for the channel.

Exposure (ms) – Value indicates exposure time for fluorescent image.

Filters **535** – Excitation 470 nm/Emission 535 nm

660 – Excitation 540 nm/Emission 660 nm

In addition, you can edit *Cell Type Parameters* used for the channel by changing the currently selected cell type (i.e., use the dropdown to choose another cell type) and clicking the **View** button to modify cell type parameters.

Note: When using the **Viability** Analysis Mode, cell type parameters can be set independently for each channel. However, when using the **Expression** Analysis Mode with more than one channel, cell type parameters are only required for *Channel 1*.

To manage cell types available on your instrument system or to edit/create cell types, see *Chapter 8. Managing Cell Types* on starting on page 61.

If you click the **View** button to modify cell type parameters, whether you can edit parameter settings displayed will depend on the cell type's *Locked State* indicated in the *Cell Type Details* section.

Note: You can only edit parameter settings for cell types that are *Unlocked*. However, for *Revvity System* or *Locked* cell types you can click the **Save As** button located at the bottom of the screen to save cell type parameters by entering a new name and clicking **Save**, then edit its parameter settings accordingly. When you click **Save and Back**, the cell type currently selected has been automatically updated to the cell type you just saved/modified.

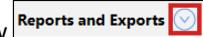
If the cell type is *Unlocked* you can edit parameter settings and then click **Save and Back** to save your changes. See *Editing a Cell Type* on page 65 for more information on editing cell type parameter settings.

Managing Assay Reports and Exports

In the *Reports and Exports* area you can assign report templates to control how data is displayed, exported and printed. In addition to the **Well View** tab which will always be displayed, a total of five (5) custom reporting tabs may be enabled for display at any one time.

Note: As of the Matrix v4.0 release, report templates are no longer limited to the *Display*, *Export*, or *Print* template types used in earlier releases. A report template can be created for all these purposes and multiple templates assigned to an assay as needed. In addition, report tabs enabled by default will vary based on the Matrix software version associated with an assay (e.g., **Well View**, **Summary** and **.csv** tabs are displayed for v4.0 assays; **Well View** and **All Wells** tabs for v3.0 assays). Users can change report templates assigned to these tabs, edit the templates directly (i.e., any changes made will also be applied to other scan results using these templates), disable them from displaying as tabs, or delete them from the assay.

If *Imaging and Analysis* parameters are not displayed, expand the area by clicking the down arrow to view available options.



Reports and Exports ⌵

Display: MX5_Display_2FL Viability_Well View ⌵ View **Report template used for Well View tab**

Exports

Images: Raw Images Colorized Images

Data: Well Level CSV Object Level CSV Object Level ACS

Archive: Data Set

Reports Create Delete View Move Up Move Down

Report Template	Display	CSV	Excel
	Tab Name	Export	Export
		Auto Open	Auto Open
		Print	Print
MX5_Display_2FL Viability	Summary		
Default_Export_2FL Viability	.csv		

Default report templates enabled as tabs

To change the report template currently assigned to the **Well View** tab, use the **Display** dropdown to select a new template. Click **View** to confirm the selected template suits your display needs and edit it if necessary.

To define assay *Exports* such as *Images*, *Data*, and *Archive* output files (i.e., exported automatically after analysis is complete), select any of the following options.

- Raw Images** Represents *Black and White* high-resolution PNG images.
- Colorized Images** Represents *Fluoresced* high-resolution PNG images.
- Well Level CSV** Represents well-level data including well name, calculation run, channel, count, mean size, etc. for each selected well in Comma Separated Values (CSV) format.
- Object Level CSV** Represents object-level data including type, classification, size, circularity, area, perimeter, etc. in Comma Separated Values (CSV) format.
- Object Level ACS** Represents object-level data including type, classification, size, circularity, area, perimeter, etc. in Image Cytometry Experiment (ICE) format. *Data may be associated with an ACS template for import into the De Novo Software FCS Express application. See Chapter 9. Managing ACS Templates starting on page 71.*
- Data Set** Represents a database file containing all images, results, assays, cell types and report templates associated with the scan result.

To select a report template and edit how it is being applied to an assay, double-click it in the *Reports* area, or click on a report in the list once to highlight it and then click the **View** button.

Reports										
Create Delete View Move Up Move Down										
Report Template	Display	CSV	Excel	PDF	Auto Open	Print	Auto Open	Print	Auto Open	Print
	Tab Name	Export	Export	Export						
MX5_Display_2FL Viability	Summary									
Default_Export_2FL Viability	.csv									

You can use report tab options to disable/enable the display of report templates as tabs and edit/enter tab names. A total of five (5) custom tabs may be enabled for display at any one time. In addition, you can select output file types for report templates, and indicate if files are to be opened automatically or printed upon report generation. Click the **Update Report** button to save your changes and view the updated Reports List.

Reports																									
Create Delete View Move Up Move Down																									
Report Template:	Default_Export_2FL Viability	File Type <table border="1"> <thead> <tr> <th></th> <th>Auto Open</th> <th>Print</th> </tr> </thead> <tbody> <tr> <td><input type="checkbox"/> CSV</td> <td>No Yes</td> <td>No Yes</td> </tr> <tr> <td><input type="checkbox"/> Excel</td> <td>No Yes</td> <td>No Yes</td> </tr> <tr> <td><input type="checkbox"/> PDF</td> <td>No Yes</td> <td>No Yes</td> </tr> <tr> <td><input type="checkbox"/> Word</td> <td>No Yes</td> <td>No Yes</td> </tr> </tbody> </table>										Auto Open	Print	<input type="checkbox"/> CSV	No Yes	No Yes	<input type="checkbox"/> Excel	No Yes	No Yes	<input type="checkbox"/> PDF	No Yes	No Yes	<input type="checkbox"/> Word	No Yes	No Yes
	Auto Open	Print																							
<input type="checkbox"/> CSV	No Yes	No Yes																							
<input type="checkbox"/> Excel	No Yes	No Yes																							
<input type="checkbox"/> PDF	No Yes	No Yes																							
<input type="checkbox"/> Word	No Yes	No Yes																							
Display in Tab:	No Yes	Report Tab Options																							
Tab Name:	.csv	Output File Options																							
										Update Report Cancel															

To manage this list for an assay, select a template and click the **Create/Delete** or **Move Up/Move Down** buttons.

Reports										
Create Delete View Move Up Move Down										
Report Template	Display	CSV	Excel	PDF	Auto Open	Print	Auto Open	Print	Auto Open	Print
	Tab Name	Export	Export	Export						
MX5_Display_2FL Viability	Summary									
Default_Export_2FL Viability	.csv	✓								
MX5_Growth Curve	Growth Curve		✓							✓
MX5_Display_2FL Viability Plate View 12x2										

- Add a report template by clicking the **Create** button. Select a template in the **Report Template** field, choose if the report should be enabled as a tab for display (i.e., *Yes*) and enter a tab name. Select any output file types for the report, and indicate if they are to be opened automatically (*Auto Open*) or printed upon generation whenever the assay is used to perform a count/recount. *Output of files is independent of whether a template is also enabled as a tab.* Click the **Add Report** button to add the report to the Reports List.
- Delete a template by clicking on the report once to highlight it and then clicking the **Delete** button. *If report template deleted was from the Report Template Library provided with the Matrix software, it can be re-imported if necessary. If report template was from a custom library or created using the Matrix software, it may be permanently deleted unless it was exported to an external location and saved prior to deletion.*
- Change the order in which report tabs are displayed by changing the sequence of templates in the Report List. Highlight a report template to select it and click the **Move Up** or **Move Down** buttons accordingly.

For complete step-by-step instructions on setting reporting options for an assay such as changing the display template in the **Well View** report, managing an assay Report List or adding report templates, see *Chapter 5. Customizing Scan Result Reports* starting on page 31 for details.

CREATING A NEW ASSAY

You can create a new assay by clicking the **Manage** tab > *Assays* option to display the *Assays List* and then clicking the **Create** button located at the bottom of the screen.

Enter basic *Assay Details* (such as name, description, category, and product type), *Imaging and Analysis* parameter settings, and *Reporting* options, and then click the **Save and Back** button to add the assay to the *Assays List*.

To copy defined parameter settings for an existing assay to be used as a source for creating a new assay, select an assay and view its details by double-clicking it in the *Assays List*, or clicking it once to highlight it and then clicking the **View** button located at the bottom of the screen. Click the **Save As** button and enter a new name to save a copy of the assay, and then click the **Save** button.

Edit assay parameters as described in *Editing an Assay* on page 51 and click the **Save and Back** button to confirm that the new assay has been saved to the *Assays List*.

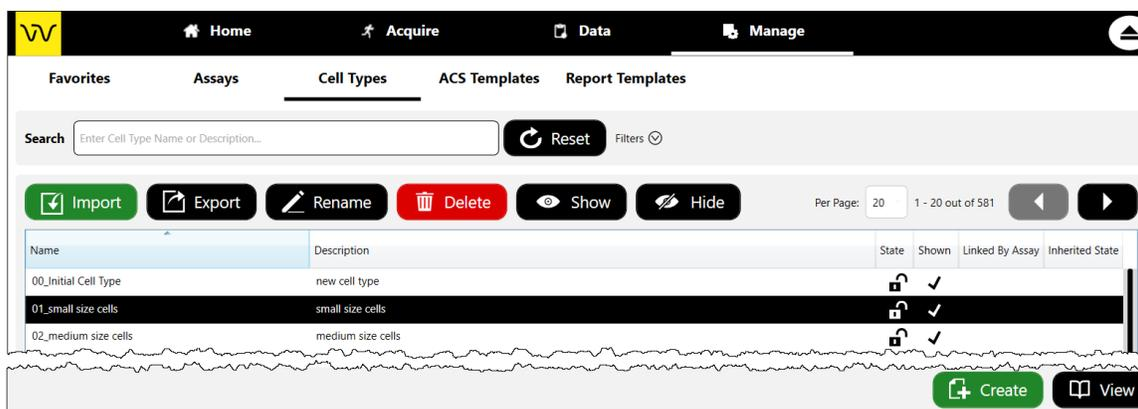
Chapter 8. Managing Cell Types

This chapter describes how cell types are displayed in the *Cell Types List* and how to manage the list to keep it current. Details on how to edit and create cell types are also provided. Cell type parameters can be optimized to meet your data acquisition needs.

Note: When migrating to the Matrix database from an earlier version, updated cell types can be imported via the software *after* migration is complete. For a list of cell types provided with the Matrix software or for assistance in defining a custom cell type, contact Support by visiting <https://www.revivity.com/contact-us> or send email to: CellC-support@revivity.com

VIEWING THE CELL TYPES LIST

You can manage cell types for your instrument system by clicking the **Manage** tab > *Cell Types* option to view the *Cell Types List*. From this screen you can import/export, rename, and delete cell types appearing in the *Cell Types List*, as well as show/hide cell types displayed in the *Cell Types* dropdown appearing in other screens.



In addition, you can select a cell type from the list to view/edit its details, create a new cell type by defining custom parameter settings or use a locked cell type as a source for creating a new cell type based on its parameter settings.

Note: If the Matrix 21 *CFR Part 11* module is enabled, the **Create** button (and **Save and Back** button when viewing cell type details) will only be available for users who have been granted permission to perform these functions.

Use the **Per Page** control to change the number of cell types displayed per page, and the arrows to move back and forth between pages in the list.

Cell types can be sorted by clicking on column headings (i.e., ascending /descending indicators will be displayed), and information presented for each cell type includes name, description, and created/last modified dates.

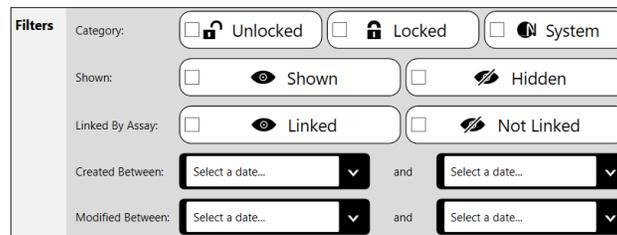
Icons displayed in the *State* column indicate whether a cell type was provided as a Revvity *System* standard (which cannot be edited), or if a cell type is currently *Locked* or *Unlocked*. When viewing *System* and *Locked* cell types, a **Save As** button allows users to copy defined parameter settings as a source for creating a new cell type.

The remaining columns display checkmarks ✓ for the cell type based on the conditions noted below:

- A checkmark displayed in the *Shown* column indicates the cell type is to be included in the *Cell Type* dropdown available in other screens (e.g., **Cell Type Parameters** field in the **Manage** tab > Assays option Details screen).
- A checkmark displayed in the *Linked by Assay* column indicates the cell type is associated with an assay. *Cell types must be unlinked from all assays before they can be deleted.*
- A checkmark displayed in the *Inherited State* column indicates the cell type was locked prior to importing it into your instrument library (i.e., its *Locked State* was inherited during import). *Keep in mind that cell types may be imported directly via the **Manage** tab > Cell Types option or as components when importing assays/datasets.*

SEARCHING FOR CELL TYPES

To search for a cell type, enter a few key characters of a cell type name or description in the **Search** field and/or expand the *Filters* area by clicking the down arrow  to specify search criteria. *Depending on monitor display size, you may need to collapse the Filters area to view search results.*



As you choose filters (e.g., by selecting *Category*, *Shown*, or *Linked By Assay* options) or enter a *Created Between/Modified Between* range of dates, the *Cell Types List* is updated automatically to display matching entries.

Category

Choose from the following categories:

Unlocked – Filters for cell types that are unlocked and can be edited

Locked – Filters for cell types that are locked and cannot be edited

System – Filters for cell types provided by Revvity (locked and cannot be edited)

System cell types are no longer available in the Cell Type Library provided with the Matrix software but may exist on your system if imported from an earlier release.

Shown

Choose from the following *Shown* states:

Shown – Filters for cell types selected to appear in the *Cell Types* dropdown

Hidden – Filters for cell types that do *not* appear in the *Cell Types* dropdown

Note: The *Cell Types* dropdown is available in other screens (e.g., **Cell Type Parameters** field in the Assay Details screen).

Linked By Assay

Choose from the following *Linked by Assay* states:

Linked – Filters for cell types that are currently associated with an assay

Not Linked – Filters for cell types that are *not* currently associated with an assay

Cell types must be unlinked from all assays before they can be deleted.

- Created Between** Enter a *Created Between* range indicating the start/end dates between which to filter for cell types created in that time frame.
- Modified Between** Enter a *Modified Between* range indicating the start/end dates between which to filter for cell types modified in that time frame.

If you find it necessary to clear all selected filters and begin the search again, click the **Reset** button.

To select a cell type and view its details, double-click it in the *Cell Types List*, or click the cell type once to highlight it and then click the **View** button located at the bottom of the screen.

MAINTAINING THE CELL TYPES LIST

When you select a cell type, buttons at the top of the *Cell Types List* become available to perform the following functions. *The **Import** button will always be enabled as it does not require the selection of a cell type in your library.*

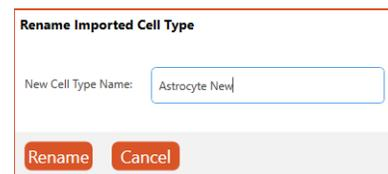
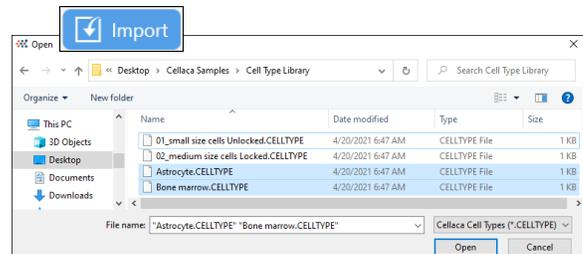
Note: If the Matrix 21 CFR Part 11 module is enabled, buttons will only be available for users who have been granted permission to perform specific functions.

Importing Cell Types

1. While viewing the *Cell Types List*, click the **Import** button.
2. Navigate to a folder where an external cell type library is available or a cell type was previously saved.
3. Select one or more *.CELLTYPE* files to be imported.
4. Click **Open**. If selected cell types already exist in your database, respond to the confirmation prompt by clicking **Yes** to overwrite the file in your system or **No** to abort the import. If you choose **No** you will be prompted to enter a new name under which to import the cell type (and then click **Rename**) or click **Cancel** to abort the import.

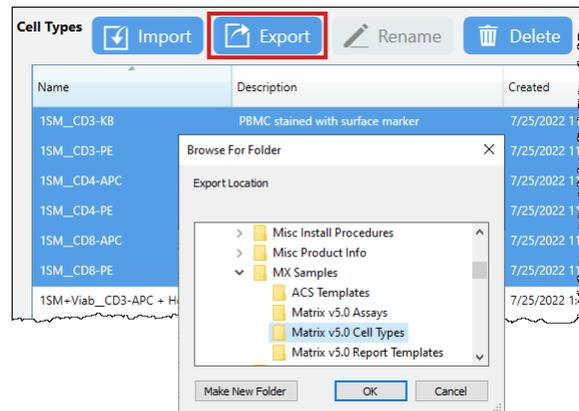
Note: When importing a *System* or *Locked* cell type, you will be prompted to import it under a different name. Click **Yes**, enter a new name, and then click **Rename** to complete the import.

5. Click **OK** to acknowledge the successful import and confirm that imported cell types are displayed in the *Cell Types List*.



Exporting Cell Types

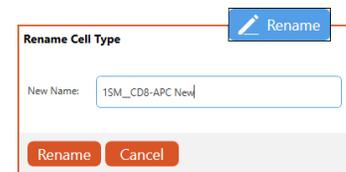
1. Select one or more cell types from the *Cell Types List* to be exported by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of cell types) or **Ctrl** key (to select non-contiguous cell types) down while clicking additional cell types.
2. Click the **Export** button.
3. Navigate to a folder where cell types are to be saved.
4. Click **OK** to save *.CELLTYPE* files in the export location.



Renaming Cell Types

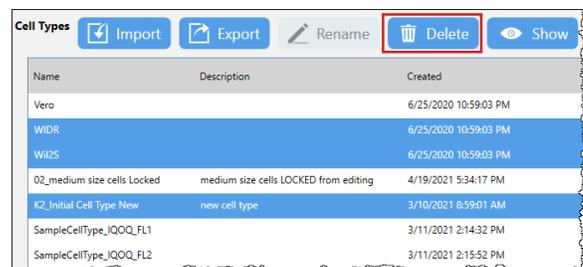
The **Rename** button will only be available for unlocked cell types.

1. Select the cell type to be renamed and click the **Rename** button.
2. Edit the cell type name and click **Rename** to save your changes.



Deleting Cell Types

1. Select one or more cell types from the *Cell Types List* to be deleted by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of cell types) or **Ctrl** key (to select non-contiguous cell types) down while clicking on additional cell types.
2. Click the **Delete** button. *If the Matrix 21 CFR Part 11 module is enabled, users will be prompted to enter a reason prior to deletion.*



Note: If cell types to be deleted were from the Cell Type Library provided with the Matrix software, they can be re-imported if necessary. If cell types were from a custom library or created using the Matrix software, they may be permanently deleted unless cell types were exported to an external location and saved prior to deletion.

Cell types must be unlinked from all assays before they can be deleted.

Showing/Hiding Cell Types in Dropdown

The *Cell Types* dropdown is available in other screens (e.g., **Cell Type Parameters** field in the Assay Details screen).

1. Select one or more cell types from the *Cell Types List* to be shown or hidden in the *Cell Types* dropdown by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of cell types) or **Ctrl** key (to select non-contiguous cell types) down while clicking on additional cell types.

- Click the **Show** or **Hide** button depending on whether you want to show/hide the selected cell types in the *Cell Types* dropdown available in other screens.

Name	Description	Created	Modified	State	Shown	Linked By Assay	Inherited State
1SM+Viab_CD3-APC + Hoechst + DeadGreen	PBMC stained with surface marker and dyes	7/25/2022 1:45:25 PM	8/26/2022 11:34:26 AM		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	
1SM+Viab_CD3-KB + Hoechst + RubyDead	PBMC stained with surface marker and dyes	7/25/2022 1:47:39 PM	8/26/2022 11:34:32 AM		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	
1SM+Viab_CD3-PE + Hoechst + RubyDead	PBMC stained with surface marker and dyes	7/25/2022 1:49:39 PM	9/26/2022 11:29:39 AM		<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	

EDITING A CELL TYPE

Editing a cell type allows users to change basic *Cell Type Details* (such as name, description and locked state), defined *Brightfield Parameters* settings and *Fluorescence Parameters* settings.

To select a cell type and view its details, double-click it in the *Cell Types List* or click the cell type once to highlight it and then click the **View** button located at the bottom of the screen.

Cell Type Details

Name:

Description:

Locked State: Unlocked Locked Last Modified On: 06/26/2020

Brightfield Parameters

Cell Attributes

Cell Diameter (µm): to

Roundness:

Contrast Enhancement:

Declustering No Yes

Edge Factor:

Threshold Factor:

Background Adjustment:

Trypan Blue

Dead Cell Diameter (µm): to

Sensitivity:

Uniformity:

Very Dim Dead Cells: No Yes

Contrast Enhancement:

Fluorescence Parameters

Cell Attributes

Thresholding Manual Auto

Auto Back

Edit the cell type as described in each of the sections presented below:

- *Basic Cell Type Details* on page 66
- *Brightfield Parameters* on page 68
- *Fluorescence Parameters* on page 69

When viewing/editing of a cell type is complete, click one of the following buttons to return to the *Cell Types List*.



Back Button: Click **Back** to return to previous screen without saving any changes. *If cell type has unsaved changes, click **OK** in response to confirmation prompt.*



Save Button: Click **Save** to save your changes and return to previous screen. Click **Auto Back** to toggle functionality of the **Save** button between **Save** (to save and continue editing) and **Save and Back** (to save and return to previous screen).



Save As Button: Click **Save As** to save your changes as a copy with a new name or select another cell type from the dropdown to override it and return to the previous screen. *If the cell type is locked, this also allows you to copy the cell type to use as a source for creating a new cell type.* Click **Auto Back** to toggle functionality of the **Save As** button between **Save As** (to save as and continue editing) and **Save As and Back** (to save as and return to previous screen).

Basic Cell Type Details

In the *Cell Type Details* area, you can edit the cell type *Name*, *Description* and *Locked State* as indicated below.

Note: You can only edit a cell type if the category displayed is *Unlocked*. If the category displayed is *Locked* or if the cell type is a Revvity *System* default, you must first copy the cell type by clicking the **Save As** button and entering a new name to save defined parameters as a source for creating a new cell type. Once saved, you can edit copied parameters for the new cell type as necessary.

To edit basic cell type details, modify information contained in the following fields.

- | | |
|---------------------|---|
| Name | Displays the name of the cell type. <i>Must be unique.</i> |
| Description | Displays a brief description which can be used to identify the cell type or include defined parameters to help users distinguish it from others presented in the <i>Cell Types List</i> . |
| Locked State | <p>Indicates if a cell type was provided as a Revvity default <input checked="" type="checkbox"/> System (i.e., <i>System</i> cell types cannot be edited) or if its current status is <i>Unlocked</i> or <i>Locked</i>. When viewing details for a <i>Locked</i> or <i>System</i> cell type, use the Save As button to copy defined parameters as a source to be used for creating a new cell type.</p> <p>The locked state of a cell type is displayed in the <i>State</i> column of the <i>Cell Types List</i> using the Revvity <i>System</i>, <i>Unlocked</i>, or <i>Locked</i> icons.</p> <p>If a cell type is unlocked, you can select the Locked <input type="checkbox"/> Locked button to prevent the cell type from being edited by other users.</p> <p>Note: It is recommended that you do <i>not</i> select the Locked button until <i>after</i> editing is complete as once the cell type is locked, you will no longer be able to make any changes.</p> |

Brightfield Parameters

In the *Brightfield Parameters* area, you can edit defined settings for *Cell Attributes*, *Declustering*, and *Trypan Blue* parameters as indicated below.

Cell Attributes

Defines basic brightfield parameter attributes for cells to be counted.

Cell Diameter (µm) Indicates range of cell diameter *minimum* to *maximum* values to be counted. *Cells with diameters that fall outside of this range will not be counted.*

Roundness Indicates minimum cell shape *roundness* factor to be counted. *Values range from 0.00 (includes all cell shapes) to 1.00 (includes only perfectly round cells).*

Contrast Enhancement Defines contrast enhancement value for cells in relation to the background. *Values range from 0.00 (cells with high contrast to background) to 0.80 (cells with low contrast to background); recommended value is 0.4.*

The screenshot shows the 'Brightfield Parameters' dialog box with the 'Cell Attributes' section active. It contains three input fields: 'Cell Diameter (µm)' with a range from 2.0 to 22.0, 'Roundness' set to 0.05, and 'Contrast Enhancement' set to 0.80.

Declustering

Defines whether individual cells within a clump are to be counted. Turning this feature off allows clumps to be counted as one unit if its diameter falls within the minimum/maximum cell diameter range. *The **Yes** button indicates declustering parameters are to be used (enabled by default). Click **No** to turn OFF declustering.*

Edge Factor Indicates degree to which cell edges must be enhanced for optimal declustering. *Values range from 0.0 (clearly defined edges) to 1.0 (edges difficult to distinguish from the background).*

Threshold Factor Indicates threshold ratio between cell signal and background. *Values range from 0.0 (cell signal to background is very low) to 1.0 (cell signal to background is high). Lower values help to normalize 'noise' from fluorescent signals.*

Background Adjustment Indicates the adjustment ratio between cell signal and background. *Values range from 0.0 (cell signal to background is very low) to 1.0 (cell signal to background is high). Lower values help to normalize 'noise' from brightfield signals.*

The screenshot shows the 'Declustering' dialog box. It features a 'Declustering' section with 'No' and 'Yes' buttons, where 'Yes' is selected. Below are three input fields: 'Edge Factor' set to 0.7, 'Threshold Factor' set to 1.0, and 'Background Adjustment' set to 1.0.

Trypan Blue

Defines trypan blue viability cell detection parameter settings for cells to be counted.

Dead Cell Diameter (µm)	Indicates range of dead cell diameter <i>minimum</i> to <i>maximum</i> values to be counted. <i>Dead cells with diameters that fall outside of this range will not be counted.</i>
Sensitivity	Adjusts sensitivity of the camera to darkness level of trypan blue stained cells to be counted. <i>Values range from 0.00 (detects very dark stained cells) to 10.0 (detects more mixed staining populations).</i>
Uniformity	Indicates trypan blue staining uniformity to be counted. <i>Values range from 100 (stained cells are all uniform in color) to 255 (stained cells have non-uniform dark and light areas).</i>
Very Dim Dead Cells	When feature is enabled (<i>Yes</i>), helps to detect very dim stained dead cells.
Contrast Enhancement	<i>Very Dim Dead Cells</i> feature must be selected for this field to be enabled. Refines the division between the background and cells with low contrast (i.e., no defined edges). <i>Suggested enhancement value is 0.60; values range from 0.00 (stained cells have medium contrast to background) to 0.80 (stained cells have very low contrast to background).</i>

Trypan Blue

Dead Cell Diameter (µm): to

Sensitivity:

Uniformity:

Very Dim Dead Cells: No Yes

Contrast Enhancement:

Fluorescence Parameters

Defines fluorescent parameter attributes for cells to be counted.

Cell Diameter (μm) Indicates range of cell diameter *minimum* to *maximum* values to be counted. *Cells with diameters that fall outside of this range will not be counted.*

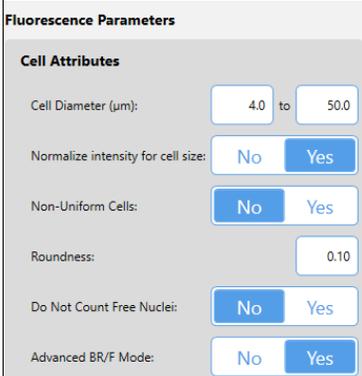
Normalize Intensity for Cell Size When feature is enabled (*Yes*), adjusts fluorescent intensity to the size of the cell. For example, if cells don't uptake the staining all at once and both light and dark areas exist within cells, enabling this feature normalizes FL intensity throughout the entire cell rather than having multiple intensity readings.

Non-Uniform Cells When feature is enabled (*Yes*), identifies a cell as positive even if only a portion of the cell has fluorescence intensity above the defined threshold. Enabling this feature is useful if cells in the sample have a mixed staining population that contains both light and dark areas.

Roundness Indicates minimum cell shape *roundness* factor to be counted. *Values range from 0.00 (includes all cell shapes) to 1.0 (includes only perfectly round cells).*

Do Not Count Free Nuclei When feature is enabled (*Yes*), uses a proprietary image analysis algorithm that excludes free floating nuclei from being counted as cells.

Advanced BR/F Mode When feature is enabled (*Yes*), uses a proprietary counting method combining brightfield and fluorescence imaging to perform enhanced declustering.



Fluorescence Parameters

Cell Attributes

Cell Diameter (μm): 4.0 to 50.0

Normalize intensity for cell size: No Yes

Non-Uniform Cells: No Yes

Roundness: 0.10

Do Not Count Free Nuclei: No Yes

Advanced BR/F Mode: No Yes

Thresholding

Indicates fluorescence intensity required for cells to be counted.

% of Image Range to Count Defines percentage of FL threshold to be counted. For *Manual* thresholding, entering a value in this field defines the percentage of pixels to be counted thus excluding cells that fall below that value. *Values range from 1-100.*



Thresholding Manual Auto

% of Image Range to Count: 10

Threshold Factor: 1.0

For *Auto* thresholding, this feature establishes maximum intensity of all images taken for the sample to be 100%. Entering a value in this field defines a percentage of that maximum (e.g., 10 %) thus excluding cells with an intensity that fall below that value.

A general rule is to lower the threshold value to include dim cells or increase threshold value to exclude cells.

Threshold Factor Indicates threshold ratio between cell signal and background. *Values range from 0.0 (cell signal to background is very low) to 1.0 (cell signal to background is high). Lower values help to normalize 'noise' from fluorescent signals.*

CREATING A NEW CELL TYPE

You can create a new cell type by clicking the **Manage** tab > *Cell Types* option to display the *Cell Types List* and then clicking the **Create** button located at the bottom of the screen.

Enter basic *Cell Type Details* (such as name, description, and locked state), define *Brightfield Parameters* and *Fluorescence Parameters* settings, and then click the **Save and Back** button to add the cell type to the *Cell Types List*.

To copy defined parameter settings for an existing cell type to be used as a source for creating a new cell type, select a cell type and view its details by double-clicking it in the *Cell Types List*, or clicking it once to highlight it and then clicking the **View** button located at the bottom of the screen. Click the **Save As** button and enter a new name to save a copy of the cell type, and then click the **Save** button.

Edit cell type parameters as described in *Editing a Cell Type* on page 65 and click the **Save and Back** button to confirm that the new cell type has been saved to the *Cell Types List*.

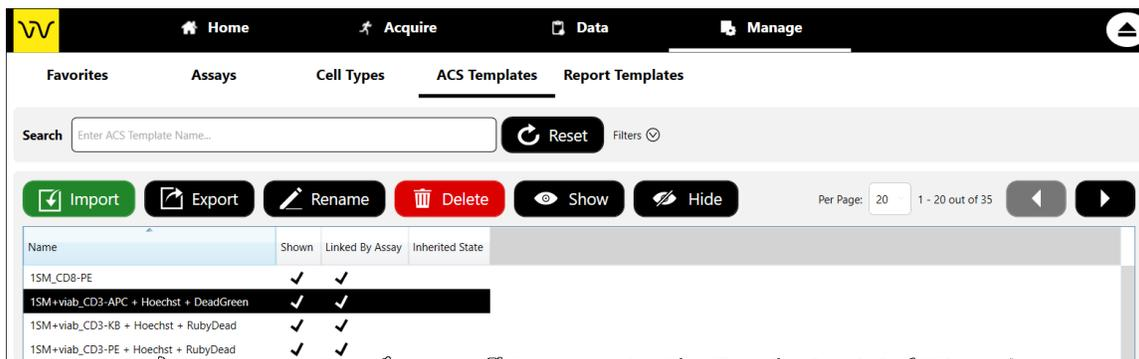
Chapter 9. Managing ACS Templates

This chapter describes how ACS templates are displayed in the *ACS Templates List* and how to manage the list to keep it current. ACS templates can be assigned to assays if the *Object Level ACS* export option is selected.

ACS templates can be assigned to assays as an output file type when exported data in Image Cytometry Experiment (ICE) format is to be imported into the De Novo Software FCS Express application.

VIEWING THE ACS TEMPLATES LIST

You can manage ACS templates available to your instrument system by clicking the **Manage** tab > *ACS Templates* option to view the *ACS Templates List*. From this screen you can import/export, rename, and delete ACS templates appearing in the *ACS Templates List*, as well as show/hide ACS templates displayed in the *ACS Templates* dropdown appearing in other screens.



Note: ACS templates can only be created, edited, and viewed in the De Novo Software FCS Express application, and then imported via the Matrix software for your instrument system. In addition, no verification is performed on ACS templates other than checking for the .FEY file extension.

Use the **Per Page** control to change the number of ACS templates displayed per page, and the arrows to move back and forth between pages in the list.

ACS Templates can be sorted by clicking on column headings (i.e., ascending/descending indicators will be displayed), and information presented for each template includes name and created/last modified dates.

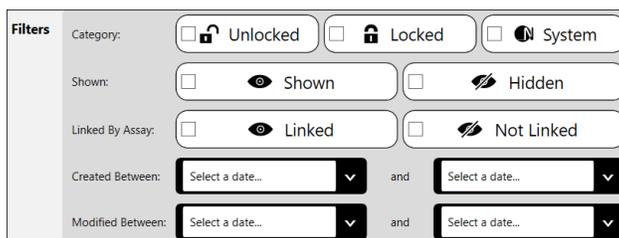
The remaining columns display checkmarks for the ACS template based on the conditions noted below:

- A checkmark displayed in the *Shown* column indicates the ACS template is to be included in the *ACS Templates* dropdown available in other screens (e.g., in *Reports and Exports/Data* section of Assay Details screen).
- A checkmark displayed in the *Linked by Assay* column indicates the ACS template is associated with an assay. *ACS templates must be unlinked from all assays before they can be deleted.*

- A checkmark displayed in the *Inherited State* column indicates the ACS template was locked prior to importing it into your instrument library (i.e., its *Locked State* was inherited during import). *Keep in mind that ACS templates may be imported directly via the **Manage** tab > ACS Templates option or as components when importing assays/datasets.*

SEARCHING FOR ACS TEMPLATES

To search for an ACS template, enter a few key characters of an ACS template name or description in the **Search** field and/or expand the *Filters* area by clicking the down arrow  to specify search criteria. *Depending on monitor display size, you may need to collapse the Filters area to view search results.*



As you choose filters (e.g., by selecting *Category*, *Shown*, or *Linked By Assay* options) or enter a *Created Between/Modified Between* range of dates, the *ACS Templates List* is updated automatically to display matching entries.

Category

Choose from the following categories:

Unlocked – Filters for ACS templates that are unlocked and can be edited

Locked – Filters for ACS templates that are locked and cannot be edited

System – Filters for ACS templates provided by Revvity (locked and cannot be edited)

System ACS templates may not be included in a release.

Shown

Choose from the following *Shown* states:

Shown – Filters for ACS templates selected to appear in *ACS Templates* dropdown

Hidden – Filters for ACS templates that do *not* appear in *ACS Templates* dropdown

The *ACS Templates* dropdown is available in other screens (e.g., in *Reports and Exports/ Data* section of the Assay Details screen).

Linked By Assay

Choose from the following *Linked by Assay* states:

Linked – Filters for ACS templates that are currently associated with an assay

Not Linked – Filters for ACS templates that are *not* currently associated with an assay

ACS Templates must be unlinked from all assays before they can be deleted.

Created Between

Enter a *Created Between* range indicating the start/end dates between which to filter for ACS templates created in that time frame.

Modified Between Enter a *Modified Between* range indicating the start/end dates between which to filter for ACS templates modified in that time frame. *As ACS Templates cannot be modified in the Matrix software, modification dates for templates file are updated only when you overwrite an existing file on import.*

If you find it necessary to clear all selected filters and begin the search again, click the **Reset** button.

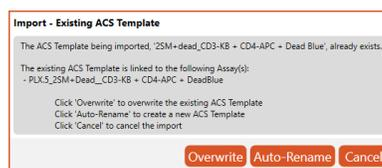
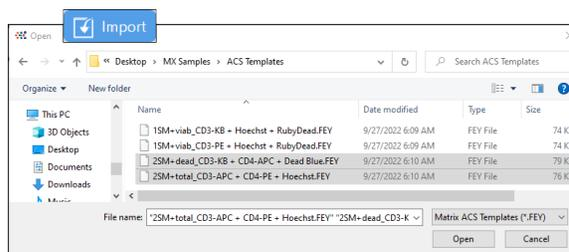
MAINTAINING THE ACS TEMPLATES LIST

When you select an ACS template, buttons at the top of the *ACS Templates List* become available to perform the following functions. *The **Import** button will always be enabled as it does not require selection of an ACS template in your library.*

Note: If the Matrix 21 *CFR Part 11* module is enabled, buttons will only be available for users who have been granted permission to perform specific functions.

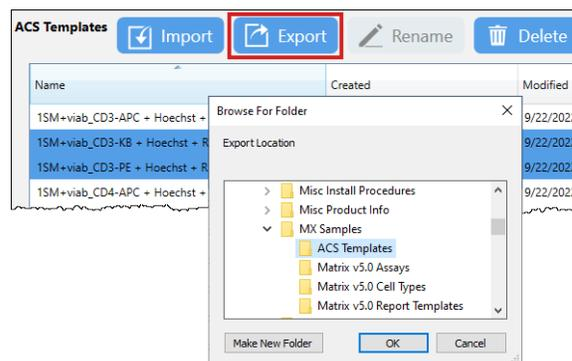
Importing ACS Templates

1. While viewing the *ACS Templates List*, click the **Import** button.
2. Navigate to a folder where an ACS template library is available or an ACS template was previously saved.
3. Select one or more *.FEY* files to be imported.
4. Click **Open**. If selected ACS templates already exist in your database, respond to the confirmation prompt by clicking **Overwrite** to overwrite the file in your system, **Auto-Rename** to automatically add “(#)” to the end of the file name indicating the imported file is a copy (where # represents a value of 1, 2, 3, etc.) or **Cancel** to abort the import.
5. Click **OK** to acknowledge the successful import and confirm that imported ACS templates are displayed in the *ACS Templates List*.



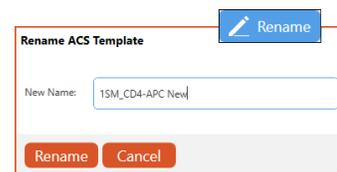
Exporting ACS Templates

1. Select one or more ACS templates from the *ACS Templates List* to be exported by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of templates) or **Ctrl** key (to select non-contiguous templates) down while clicking additional ACS templates.
2. Click the **Export** button.
3. Navigate to a folder where ACS templates are to be saved.
4. Click **OK** to save *.FEY* files in the export location.



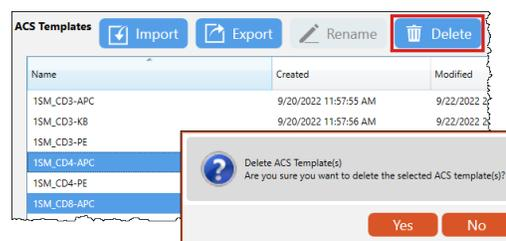
Renaming ACS Templates

1. Select the ACS template to be renamed and click the **Rename** button.
2. Edit the ACS template name and click **Rename** to save your changes.



Deleting ACS Templates

1. Select one or more ACS templates from the *ACS Template List* to be deleted by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of templates) or **Ctrl** key (to select non-contiguous templates) down while clicking on additional ACS templates.
- Click the **Delete** button followed by **Yes** to confirm the action.



If the Matrix 21 CFR Part 11 module is enabled, users will be prompted to enter a reason prior to deletion.

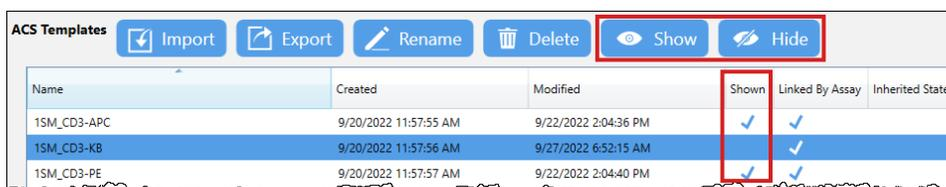
Note: If ACS template deleted was from the ACS Template Library provided with the Matrix software, it can be re-imported if necessary. If ACS template was from a custom library, it may be permanently deleted unless ACS template was exported to an external location and saved prior to deletion.

ACS Templates must be unlinked from all assays before they can be deleted.

Showing/Hiding ACS Templates in Dropdown

The ACS Templates dropdown is available in other screens (e.g., in Reports and Exports/Data section of Assay Details screen).

1. Select one or more ACS templates from the ACS Templates List to be shown or hidden in the ACS Templates dropdown by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of templates) or **Ctrl** key (to select non-contiguous templates) down while clicking on additional ACS templates.
2. Click the **Show** or **Hide** button depending on whether you want to show/hide the selected ACS templates in the ACS Templates dropdown available in other screens.



Name	Created	Modified	Shown	Linked By Assay	Inherited State
1SM_CD3-APC	9/20/2022 11:57:55 AM	9/22/2022 2:04:36 PM	✓	✓	
1SM_CD3-KB	9/20/2022 11:57:56 AM	9/27/2022 6:52:15 AM	✓	✓	
1SM_CD3-PE	9/20/2022 11:57:57 AM	9/22/2022 2:04:40 PM	✓	✓	

This page intentionally left blank.

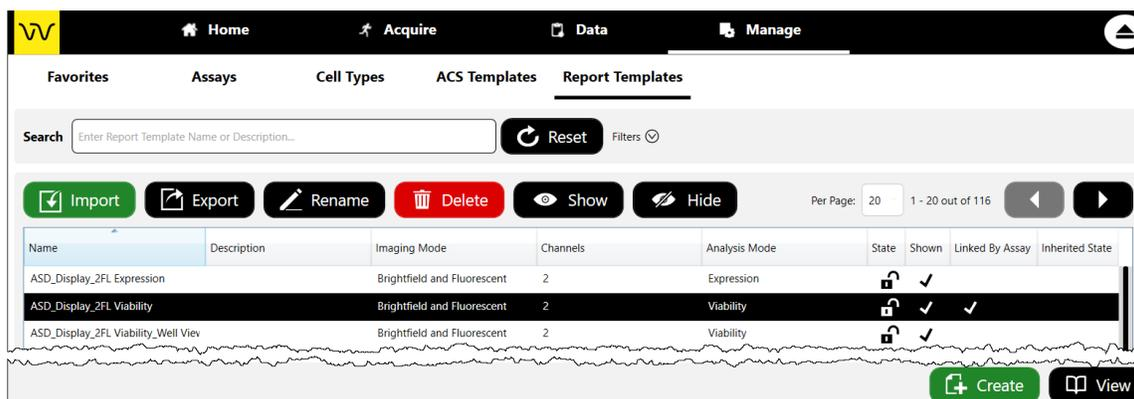
Chapter 10. Managing Report Templates

This chapter describes how report templates are displayed in the *Report Templates List* and how to manage the list to keep it current. An overview on how to edit and create report templates is also provided. Report templates are used to display, export, and print data, and can be customized to meet your reporting needs.

Note: When migrating to the Matrix database from an earlier version, updated report templates can be imported via the software *after* migration is complete. For list of report templates provided with Matrix software or for assistance in defining a custom report template, contact Support by visiting <https://www.revvy.com/contact-us> or send email to: CellC-support@revvity.com

VIEWING THE REPORT TEMPLATES LIST

You can manage report templates for your instrument system by clicking the **Manage** tab > *Report Templates* option to view the *Report Templates List*. From this screen you can import/export, rename, and delete report templates appearing in the *Report Templates List*, as well as show/hide report templates displayed in the *Report Templates* dropdown appearing in other screens.



In addition, you can select a report template from the list to view/edit its details, create a new report template by adding custom report elements/defining parameter settings, or use a locked report template as a source for creating a new report template based on its elements/parameter settings.

Note: If the Matrix 21 CFR Part 11 module is enabled, the **Create** button (and **Save and Back** button when viewing report templates) will only be available for users who have been granted permission to perform these functions.

Use the **Per Page** control to change the number of report templates displayed per page, and the arrows to move back and forth between pages in the list.

Report Templates can be sorted by clicking on column headings (i.e., ascending/descending indicators will be displayed), and information presented for each report template includes name, description, imaging mode/ number of channels, analysis mode defined for the report template, and created/last modified dates.

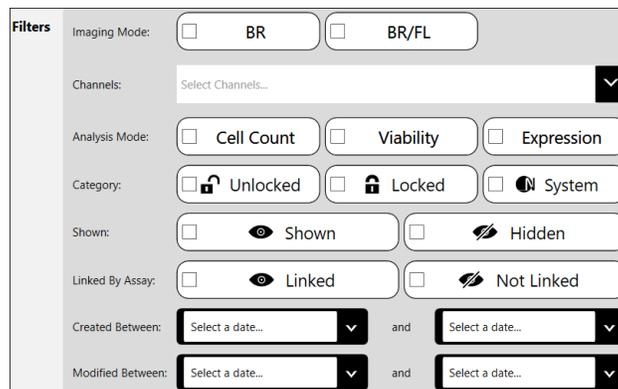
Icons displayed in the *State* column indicate whether a report template was provided as a Revvity *System* standard (which cannot be edited), or if a report template is currently *Locked* or *Unlocked*. When viewing *System* and *Locked* report templates, a **Save As** button allows users to copy report elements and defined parameter settings as a source for creating a new report template.

The remaining columns display checkmarks ✓ for the report template based on the conditions noted below:

- A checkmark displayed in the *Shown* column indicates the report template is to be included in the *Templates* dropdowns available in other screens (e.g., **Display Template**, **Print Template**, and **Export Template** fields in the Assay Details screen).
- A checkmark displayed in the *Linked by Assay* column indicates the report template is associated with an assay. *Report templates must be unlinked from all assays before they can be deleted.*
- A checkmark displayed in the *Inherited State* column indicates the report template was locked prior to importing it into your instrument library (i.e., its *Locked State* was inherited during import). *Keep in mind that report templates may be imported directly via the **Manage** tab > Report Templates option or as components when importing assays/datasets.*

SEARCHING FOR REPORT TEMPLATES

To search for a report template, enter a few key characters of a report template name or description in the **Search** field and/or expand the *Filters* area by clicking the down arrow  to specify search criteria. *Depending on monitor display size, you may need to collapse the Filters area to view search results.*



As you choose filters (e.g., by selecting *Imaging Mode*, *Channels*, *Analysis Mode*, *Category*, *Shown*, or *Linked By Assay* options) or enter a *Created Between/Modified Between* range of dates, the *Report Templates List* is updated automatically to display matching entries.

Imaging Mode

Choose from the following imaging modes:

BR – Filters for report templates with Brightfield

BR/FL – Filters for report templates with Brightfield and Fluorescence

Channels

Use the dropdown to select one or more channels to be included in the search, then click **OK**.

Report templates may include up to six fluorescent channels as defined by the associated assay.

**Analysis Mode**

Choose from analysis modes that can be associated with a template (e.g., *Cell Count, Viability, Expression*). *Analysis modes available for a report template will vary based on the imaging mode selected for the template.*

Category

Choose from the following categories:

Unlocked – Filters for report templates that are unlocked and can be edited

Locked – Filters for report templates that are locked and cannot be edited

System – Filters for report templates provided by Revvity (locked and cannot be edited)

System report templates are no longer available in the Report Template Library provided with the Matrix software but may exist on your system if imported from an earlier release.

Shown

Choose from the following *Shown* states:

Shown – Filters for report templates selected to appear in *Templates* dropdowns

Hidden – Filters for report templates that do *not* appear in *Templates* dropdowns

The *Templates* dropdowns are available in other screens (e.g., **Display Template**, **Export Template**, and **Print Template** fields in the Assay Details screen).

Linked By Assay

Choose from the following *Linked by Assay* states:

Linked – Filters for report templates that are currently associated with an assay

Not Linked – Filters for report templates that are *not* currently associated with an assay

When saving changes made to a report template linked by other assays, you will be prompted with a message indicating that all linked assays will also be updated.

Report Templates must be unlinked from all assays before they can be deleted.

Created Between

Enter a *Created Between* range indicating the start/end dates between which to filter for report templates created in that time frame.

Modified Between

Enter a *Modified Between* range indicating the start/end dates between which to filter for report templates modified in that time frame.

If you find it necessary to clear all selected filters and begin the search again, click the **Reset** button.

To select a report template and view its details, double-click it in the *Report Templates List* or click the report template once to highlight it and then click the **View** button located at the bottom of the screen.

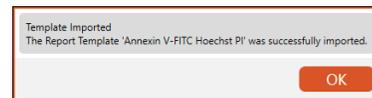
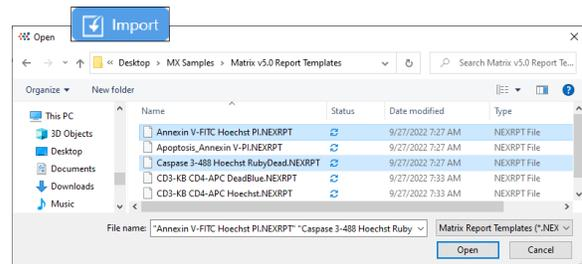
MAINTAINING THE REPORT TEMPLATES LIST

When you select a report template, buttons at the top of the *Report Templates List* become available to perform the following functions. The **Import** button will always be enabled as it does not require the selection of a report template in your library.

Note: If the Matrix 21 CFR Part 11 module is enabled, buttons will only be available for users who have been granted permission to perform specific functions.

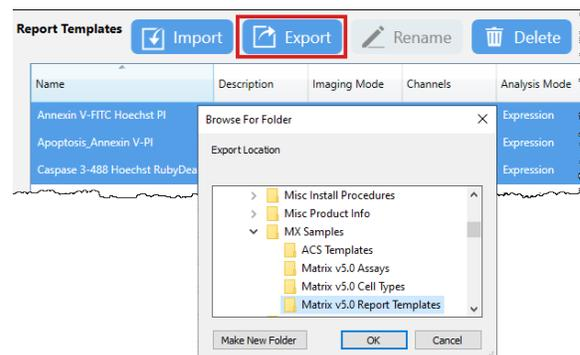
Importing Report Templates

1. While viewing the *Report Templates List*, click the **Import** button.
 2. Navigate to a folder where an external report template library is available or a report template was previously saved.
 3. Select one or more *.NEXRPT* files to be imported.
 4. Click **Open**. If selected report templates already exist in your database, respond to the confirmation prompt by clicking **Yes** to overwrite the file in your system or **No** to abort the import. If you choose **No** you will be prompted to enter a new name under which to import the report template (and then click **Rename**) or click **Cancel** to abort the import.
- Note:** When importing a *System* or *Locked* report template, you will be prompted to import it under a different name. Click **Yes**, enter a new name, and then click **Rename** to complete the import.
5. Click **OK** to acknowledge the successful import and confirm that imported report templates are displayed in the *Report Templates List*.



Exporting Report Templates

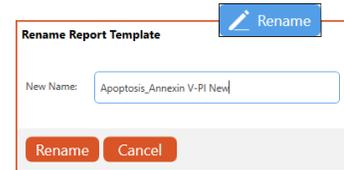
1. Select one or more report templates from the *Report Templates List* to be exported by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of templates) or **Ctrl** key (to select non-contiguous templates) down while clicking additional report templates.
2. Click the **Export** button.
3. Navigate to a folder where report templates are to be saved.
4. Click **OK** to save *.NEXRPT* files in the export location.



Renaming Report Templates

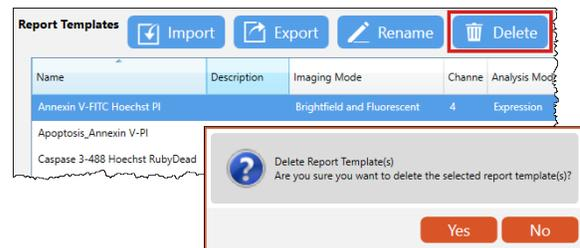
The **Rename** button will only be available for unlocked report templates.

1. Select the report template to be renamed and click the **Rename** button.
2. Edit the report template name and click **Rename** to save your changes.



Deleting Report Templates

1. Select one or more report templates from the *Report Template List* to be deleted by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of templates) or **Ctrl** key (to select non-contiguous templates) down while clicking on additional report templates.



Click the **Delete** button followed by **Yes** to confirm the action. *If the Matrix 21 CFR Part 11 module is enabled, users will be prompted to enter a reason prior to deletion.*

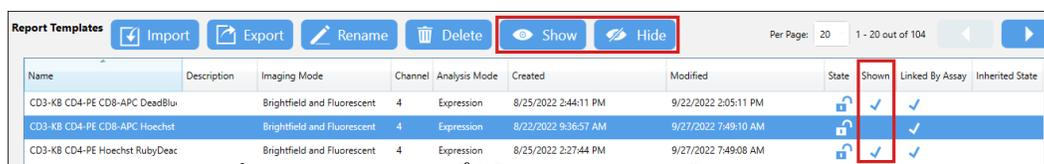
Note: If report template deleted was from the Report Template Library provided with the Matrix software, it can be re-imported if necessary. If report template was from a custom library or created using the Matrix software, it may be permanently deleted unless report template was exported to an external location and saved prior to deletion.

Report templates must be unlinked from all assays before they can be deleted.

Showing/Hiding Report Templates in Dropdown

The *Templates dropdown* is available in other screens (e.g., **Display Template**, **Export Template**, and **Print Template** fields in the Assay Details screen).

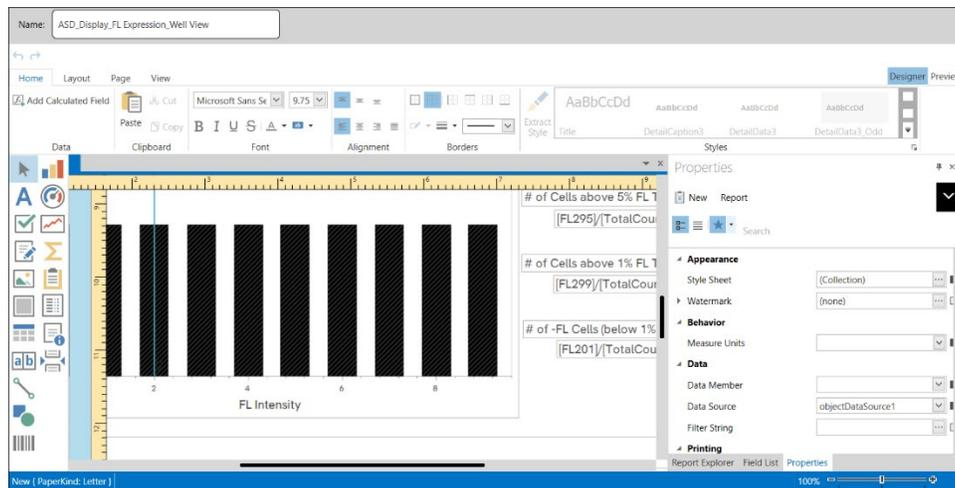
1. Select one or more report templates from the *Report Templates List* to be shown or hidden in the *Templates dropdown* by clicking once to highlight the first one, and then holding either the **Shift** key (to select a block of templates) or **Ctrl** key (to select non-contiguous templates) down while clicking on additional templates.
2. Click the **Show** or **Hide** button depending on whether you want to show/hide the selected report templates in the *Templates dropdown* available in other screens.



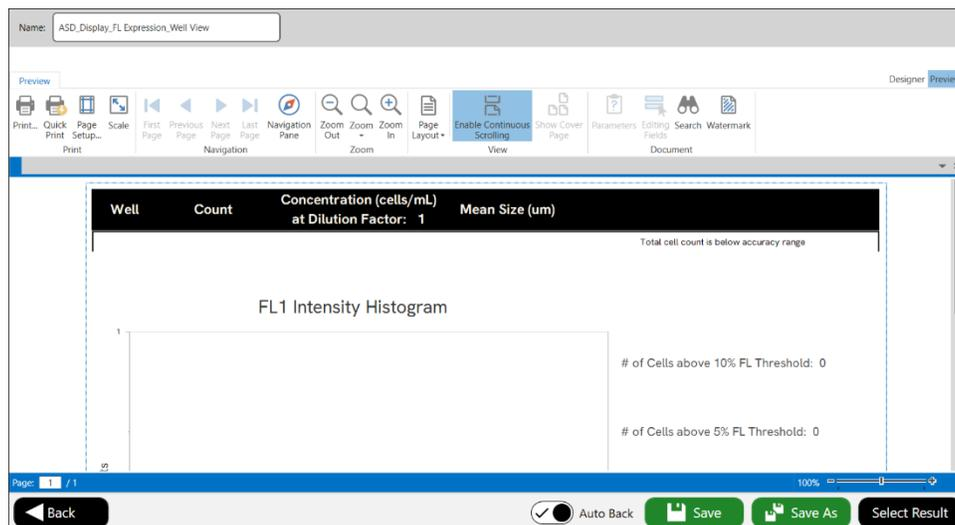
EDITING A REPORT TEMPLATE

Editing a report template allows users to customize the overall structure and individual elements used to comprise the report, as well as to modify basic *Report Template Details* (e.g., name, description, imaging mode/number of channels, analysis mode and locked state) available on the Save Report Template As screen.

To select a report template and view its details, double-click it in the *Report Templates List* or click the report template once to highlight it and then click the **View** button located at the bottom of the screen.



Report templates can be viewed in two modes – *Designer* (shown in sample screen above) and *Preview* (shown in sample screen below). *Designer* mode is displayed by default and allows you to edit template report elements while *Preview* mode displays the final layout of the template and allows you to populate it with sample data.



While viewing a report template, you can identify if it is locked (representing either a *System* or *Locked* template) based on whether the **Save and Back** button is available.

Edit the report template as described in each of the sections presented below:

- *Editing Basic Report Template Details*, below
- *Editing Report Elements* on page 85

When viewing/editing of a report template is complete, click one of the following buttons to return to the *Report Templates List*, edit defined parameters for the report template or view the template using a selected dataset.

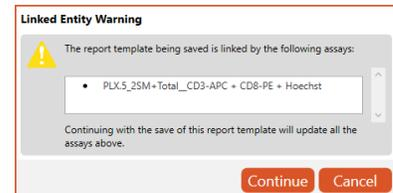


Back Button: Click **Back** to return to previous screen without saving any changes. *If report template has unsaved changes, click **OK** in response to confirmation prompt.*



Save Button: Click **Save** to save your changes and return to previous screen. Click **Auto Back** to toggle functionality of the **Save** button between **Save** (to save and continue editing) and **Save and Back** (to save and return to previous screen).

*If a report template is linked by an assay, you will be prompted to click **Continue** as changes saved will update all linked assays.*



Save As Button: Click **Save As** to save your changes as a copy with a new name or select another report template from the dropdown to override it and return to the previous screen. *If the report template is locked, this also allows you to copy the template to use as a source for creating a new report template.*

See *Editing Basic Report Template Details*, below for details regarding the Save Template As screen.



Select Result: Click **Select Result** to view the report template using a selected dataset. Choose a dataset from the displayed list and click **Select**. Click **Preview** to view the report with the selected data.

Editing Basic Report Template Details

When saving a report template (by clicking either the **Save and Back** or **Save As|Save As and Back** buttons), you can edit the *Name, Description, Imaging Mode/Number of Channels, Analysis Mode* and *Locked* state of the template.

Note: You can only edit a report template if it is unlocked. If the report template is locked or if the report template is a Revvity System default, you must first copy the template by clicking the **Save As** button and entering a new name. You can then edit defined parameters for the report template as necessary and click **Save** to save them as a source for a new report template.

To edit basic report template details, modify information contained in the following fields.

- Name** Displays the name of the report template. Enter a new name or use the dropdown to select a template to override. *Must be unique.*
- Description** Displays a brief description which can be used to identify the purpose of the report template.
- Imaging Mode** Indicates if Imaging Mode to be associated with the template is *Brightfield Only (BR)* or *Brightfield/Fluorescence (BR/FL)* including the number of channels. *Note that Cellometer Ascend/K2 supports up to 2 channels.*
- You can change the number of channels regardless of the previous value. In addition, the Imaging Mode selected will populate **Analysis Mode** field, below with available options.
- Analysis Mode** Indicates Analysis Mode to be used for the template. *Analysis modes displayed will vary based on the selected Imaging Mode.*
- See *Defining Assay Imaging and Analysis Parameters* on page 53 for options.
- Locked** Indicates if template is to be *Unlocked* (to allow editing) or *Locked* (to prevent editing).

Click **Save** to create the new report template and then edit its structure and elements accordingly. *Templates can only be edited if the Unlocked option was selected for the Locked field.*

Editing Report Elements

Report elements can be edited using the *Report Designer for WPF* application (a third-party plugin by *DevExpress*). See *Appendix A. Report Designer for WPF Reference* on page 105 for more information about using this application.

Key report elements are indicated below.

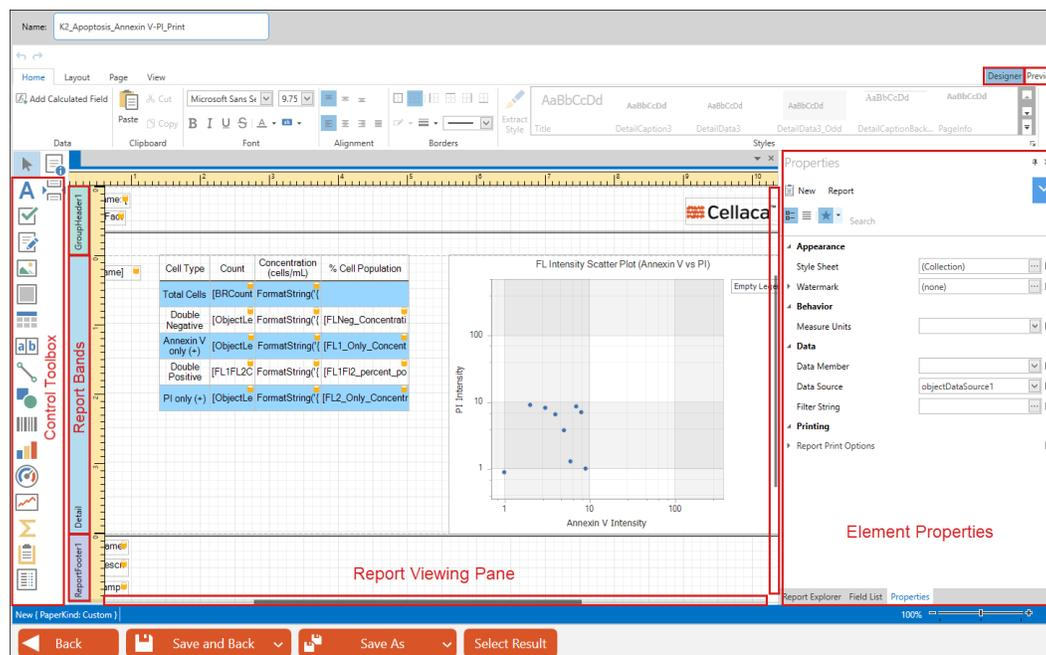
- The **Designer Tab** allows you to customize report elements using *Designer* mode. Features include:

Report Viewing Pane – Displays layout of elements in the report. Use the horizontal and vertical scroll bars to move around in the viewing pane.

Element Properties – Lists properties of elements that appear in the Report Viewing Pane. *Click on an element in the viewing page to show its properties.*

Control Toolbox – Allows you to add report controls to the template (e.g., text labels, check boxes, tables, etc. as indicated by toolbox icons). *Located on the left side of the screen.*

Report Bands – Allows you to customize the report element controlled by the band (i.e., colored boxes displayed on the left of the screen such as *GroupHeader1*, *Detail*, and *ReportFooter1* as shown in the sample below) by editing its element properties (displayed on the right of the screen). *Clicking on a specific report band varies the display of element properties.*



- The **Preview Tab** allows you to view final layout of the report template using *Preview* mode.

To add report elements to a template, click on an element in the *Control Toolbox* and drag it into the viewing pane or, click on an element in the viewing pane to reposition it and/or press **Delete** to remove it from the template.

To populate a report template with sample data, click the **Select Result** button located at the bottom of the screen to choose a scan result containing appropriate data for the template (i.e., click on the *Imaging Mode* column to sort data and choose a result using the imaging mode for which the template was designed) and click **Select**.

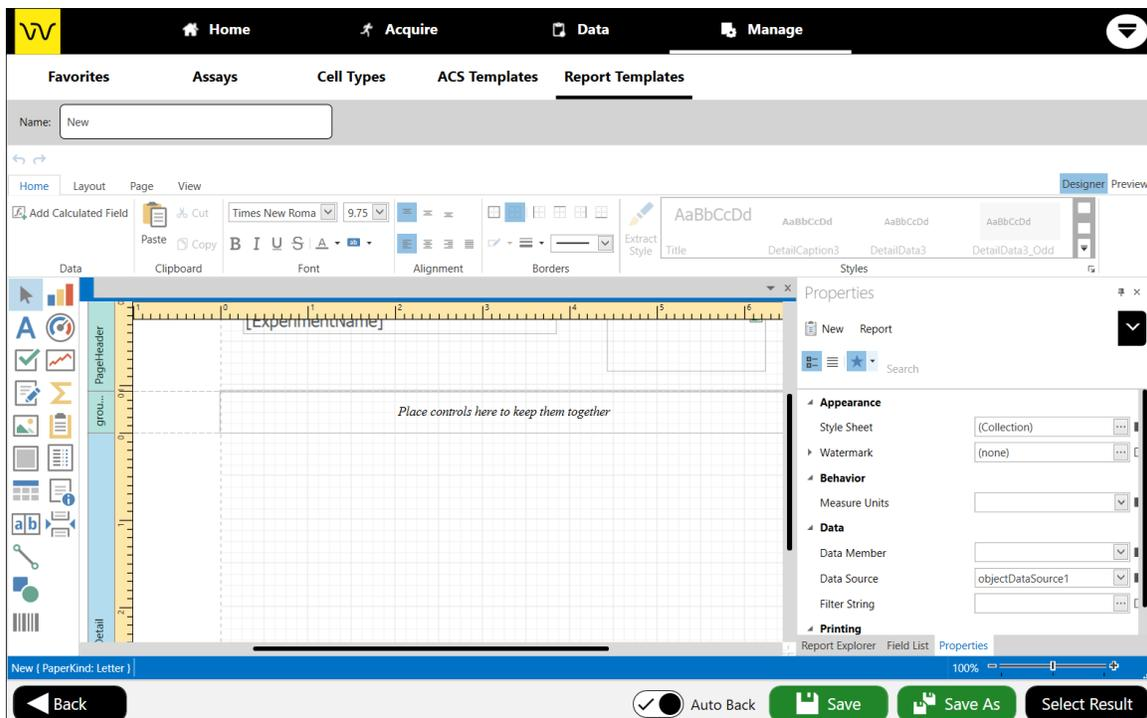
Note: Once scan result data is selected, you may need to toggle between the **Designer** and **Preview** tabs to refresh the screen and view the data.

CREATING A NEW REPORT TEMPLATE

You can create a new report template by clicking the **Manage** tab > *Report Templates* option to display the *Report Templates List* and then clicking the **Create** button located at the bottom of the screen.

Add report elements and then click the **Save and Back** button to enter basic *Report Template Details* such as the *Name*, *Description*, *Imaging Mode/Number of Channels*, *Analysis Mode*, and *Locked* state of the report template.

Note: Report elements are added using the *Report Designer for WPF* application by *DevExpress*. See *Appendix A. Report Designer for WPF Reference* on page 105 for more information about using this application.



To copy existing report elements to be used as a source for creating a new report template, select a report template and view its details by double-clicking it in the *Report Templates List*, or clicking it once to highlight it and then clicking the **View** button located at the bottom of the screen. Click the **Save As** button and enter a new name to save a copy of the report template, and then click the **Save** button.

Edit report elements as described in *Editing a Report Template* on page 82 and click the **Save and Back** button to confirm that the new report template has been saved to the *Report Template List*.

Chapter 11. Using the 21 CFR Part 11 Module

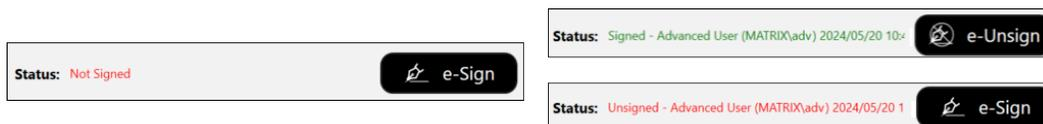
Functionality described in this chapter is available for users who have purchased a Matrix 21 CFR Part 11 module license and had the module enabled by Support. For more information about enabling this module, contact Support by visiting <https://www.revvy.com/contact-us> or send email to: CellC-support@revvy.com

FUNCTIONALITY OVERVIEW

The Matrix 21 CFR Part 11 module is compliant with the *Code of Federal Regulations (CFR) Title 21 Part 11 – Electronic Records: Electronic Signatures* published by the U. S. Food and Drug Administration (FDA). In summary, this module ensures that an organization’s use of electronic records and digital signatures in place of traditional paper-based documentation is in compliance with current FDA regulations.

Key elements in Matrix 21 CFR Part 11 module functionality are described below.

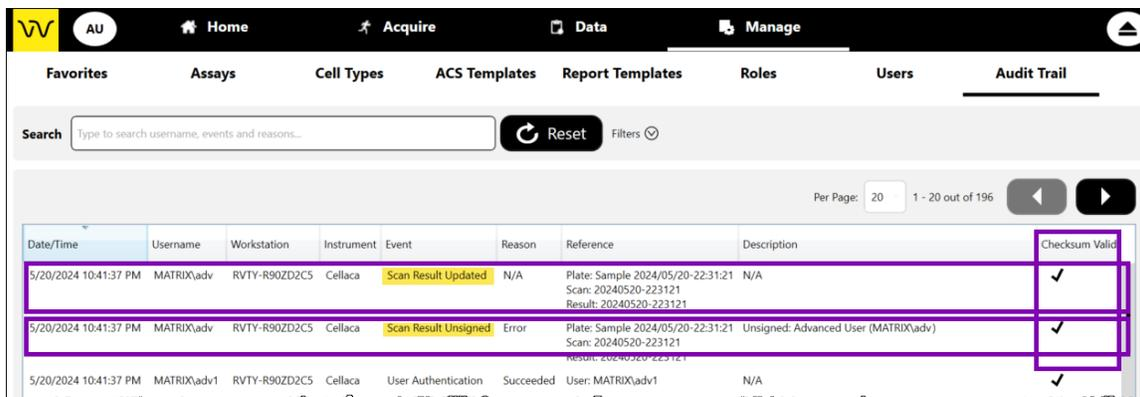
- *Electronic Signatures* are captured during the counting/analysis workflow using **e-Sign/e-Unsign** buttons and include user name, date/time when signature was executed, reason for the action, and meaning of signature (i.e., for *Review, Approval, or Rejection*). If a user’s assigned role does not have permissions to approve count results, a supervisor can log in to e-Sign the record on behalf of the user.



- *Electronic Records* are created as the result of linking electronic signatures to user actions ensuring that records have not been copied or falsified in any manner.

Consumable ID	Consumable Type	Assay	Imaging Mode	Channels	Tag	Scan Creation	Result Creation	Status	Signed/Unsigned On	User	User Name
Sample 2024/05/20-22:31:21	12x2 Plate (CHM24	MX.6_Viab_AOPL_Primary Cells	Brightfield and Fluor	2		05/20/24 22:31:21	05/20/24 22:31:21	Unsigned	05/20/24 22:41:37	MATRIX\adv	Advanced User

- An *Audit Trail* provides assurance regarding the integrity of an electronic record and continually monitors all users performing actions, the type of actions performed, and the date/time associated with user actions.



Audit trails can be retained either *Locally* on an instrument’s Operating Computer or on a *Windows Network* (i.e., using administrator-controlled access via user roles and privileges).

ESTABLISHING YOUR USER HIERARCHY

The Matrix *21 CFR Part 11* module requires that you establish a hierarchy of *Users* within your organization consisting of Administrator Users associated with full access to module functionality and other users with varying levels of access controlled through the assignment of *Roles* (i.e., sets of privileges that can be grouped and applied to multiple users sharing similar job responsibilities).

Monitoring User Actions

To create users and monitor their actions, options are added to the **Manage** tab for use by Administrator Users – *Roles*, *Users*, and *Audit Trail* – when the Matrix *21 CFR Part 11* module is enabled.

As users are created, they are added to the *Users List* which displays a summary of account details including role, personal information (*Title* and *First Name/Last Name*), and if the account is currently enabled.

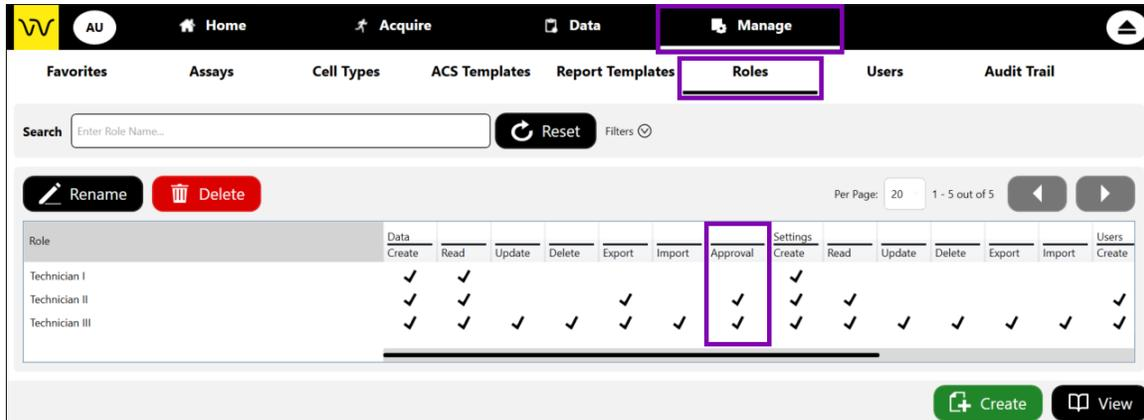
Name	Role	Title	First	Last	Enabled
MATRIX\ADM	Administrator	Administrator	Admin	User	✓
MATRIX\Tech I	Technician I	Technician I	Basic	User	✓
MATRIX\Tech II	Technician II	Technician II	Advanced	User	✓
MATRIX\Tech III	Technician III	Technician III	Supervisor	User	✓

When users log in to the Matrix software to perform daily tasks, the system automatically generates an *Audit Trail* capturing user actions and validates them against the database to ensure integrity of the data is maintained.

Date/Time	Username	Workstation	Instrument	Event	Reason	Reference	Description	Checksum Valid
5/20/2024 10:18:08 PM	MATRIX\ADM	RVTY-R90ZD2C5	Cellaca	User Added	N/A	User: Basic User	Added User: AccountName: MATRIX\Tech I FirstName: Basic LastName: User Title: Technician I IsActiveDirectoryUser: False IsEnabled: True IsSystemUser: False ForcePasswordReset: False Role Name: Technician I	✓
5/20/2024 10:17:00 PM	MATRIX\ADM	RVTY-R90ZD2C5	Cellaca	Role Added	N/A	Role: Technician I	Added Role: Role Name: Technician I PrivilegeFlags: Data, Create, Data_Read, Settings_Create IsAdministrator: False IsService: False IsSystemRole: False	✓
5/20/2024 10:16:17 PM	MATRIX\ADM	RVTY-R90ZD2C5	Cellaca	User Logged In	N/A	N/A	N/A	✓

Controlling User Access via Roles

Roles allow you to control varying levels of Matrix functionality to which specific groups of users will have access. When creating a role, select subsets of privileges that apply to users performing common tasks.



For example, you can create a basic *Technician I* role that is able to *Create* and *Read* count results but does *not* have access to add an electronic signature to approve those results. To support this basic role, you can create advanced roles (*Technicians II* and *III*) with the *Approval* privilege that are able to log in at technician workstations using the **e-Sign/e-Unsign** buttons to approve or reject displayed count results.

Understanding Role Privileges

Roles are used in the Matrix *21 CFR Part 11* module to control the following areas of responsibility:

- **Data** – Controls whether users can create, manage (i.e., read, update, and delete), export/import, and approve/reject data during the counting and analysis workflow.
- **Settings** – Controls whether users can access the **Manage** tab > *Assays, Cell Types, and Report Templates* options allowing them to create, manage (i.e., read, update, and delete), and export/import these entities.
- **Users, Roles, or Audits** – Controls whether users can access the **Manage** tab > *Users, Roles, or Audits* options (displayed only if the *21 CFR Part 11* module is enabled) allowing them to create or manage (i.e., read, update, and delete) Matrix software users and roles. In addition, controls whether users can monitor (i.e., read and export) the audit trail automatically generated by the system.

Privileges							
Data:	<input type="checkbox"/> Create	<input type="checkbox"/> Read	<input type="checkbox"/> Update	<input type="checkbox"/> Delete	<input type="checkbox"/> Export	<input type="checkbox"/> Import	<input type="checkbox"/> Approval
Settings:	<input type="checkbox"/> Create	<input type="checkbox"/> Read	<input type="checkbox"/> Update	<input type="checkbox"/> Delete	<input type="checkbox"/> Export	<input type="checkbox"/> Import	
Users:	<input type="checkbox"/> Create	<input type="checkbox"/> Read	<input type="checkbox"/> Update	<input type="checkbox"/> Delete			
Roles:	<input type="checkbox"/> Create	<input type="checkbox"/> Read	<input type="checkbox"/> Update	<input type="checkbox"/> Delete			
Audits:	<input type="checkbox"/> Read	<input type="checkbox"/> Export					

DATA PRIVILEGES

The following privileges control whether users can create, manage (i.e., read, update, and delete), export/import and approve/reject data during the counting and analysis workflow.

Note: The *Data: Read* privilege must be selected for the **Data** tab to be accessible to a user regardless of any other *Data* privileges assigned to that user.

Create Allows users to *create* a new scan result (i.e., **Count** and **Recount** buttons are enabled in the **Acquire** tab Setup screen) and count results will be displayed under the **Data** tab. *Users must have the Data: Read privilege to view scan results in the **Data** tab Select screen.*

Read Allows users to *view* all scan results available under the **Data** tab and to select a specific scan result (i.e., **View** button is enabled) to be displayed. *This privilege must be granted to all users needing to view the **Data** tab Select screen regardless of any other Data privileges assigned.*

Update *Feature not yet implemented.*

Delete Allows users to select a scan result available under the **Data** tab and delete it (i.e., **Delete** button is enabled) as necessary. *Users will be prompted to enter a reason prior to deletion.* In addition, this privilege enables the **Recover Free Space** button if users also have *sys_admin* rights to perform this action on the database server.

Export Allows users to select one or more scan results available under the **Data** tab and export them (i.e., **Export** button is enabled) to a user-defined location such as the Operating Computer desktop.

Import Allows users to select one or more scan results stored outside the Matrix software database and import them (i.e., **Import** button is enabled) to be displayed under the **Data** tab.

Approval Allows users to add an e-signature (i.e., use of **e-Sign** and **e-Unsign** buttons is authorized) for count results data currently displayed. If the logged-in user role performing the count does not include this privilege, another user with the privilege in their role can log in to e-Sign the record on their behalf.

SETTINGS PRIVILEGES

The following privileges control whether users can access the **Manage** tab > *Assays, Cell Types, ACS Templates* and *Report Templates* options allowing them to create, manage (i.e., read, update, and delete) and export/import these entities.

Note: The *Settings: Read* privilege must be selected for the **Manage** tab to be accessible to a user regardless of any other *Settings* privileges assigned to that user.

- Create** Allows users to *create* a new assay, cell type, or report template (i.e., **Create** button is enabled in the respective *Assays*, *Cell Types*, and *Report Templates* options available under the **Manage** tab) and save the entity (i.e., **Save and Back** and **Save As** buttons are enabled).
- Read** Allows users to *view* all assays, cell types, or report templates in the respective screens for these options available under the **Manage** tab and to select a specific entity (i.e., **View** button is enabled) to be displayed. *This privilege must be granted to all users needing to view the **Manage** tab > Assays, Cell Types, and Report Templates options regardless of any other Settings privileges assigned.*
- Update** Allows users to *modify* an assay, cell type, or report template in the respective screens for these options available under the **Manage** tab and save their changes (i.e., **Save and Back** and **Save As** buttons are enabled). *The selected entity must be unlocked to allow for editing.* Also enables the **Rename**, **Show**, and **Hide** buttons in the Assays, Cell Types, and Report Templates screens.
- Delete** Allows users to select an assay, cell type, ACS template, or report template in the respective screens for these options available under the **Manage** tab and delete it (i.e., **Delete** button is enabled) as necessary. *Users will be prompted to enter a reason prior to deletion.*
- Export** Allows users to select one or more assays, cell types, ACS templates, or report templates in the respective screens for these options available under the **Manage** tab and export them (i.e., **Export** button is enabled) to a user-defined location such as the Operating Computer desktop.
- Import** Allows users to select one or more assays, cell types, ACS templates, or report templates stored outside the Matrix software database and import them (i.e., **Import** button is enabled) to be displayed in the respective screens for these options to be available under the *Manage* tab.

USERS PRIVILEGES

The following privileges control whether users can access the **Manage** tab > *Users* option allowing them to create and manage Matrix software users.

Note: The *Users: Read* privilege must be selected for the **Manage** tab > *Users* option to be accessible to a user regardless of any other *Users* privileges assigned to that user.

Users: Create Read Update Delete

- Create** Allows users to *create* a new account (i.e., **Create** button is enabled) to be displayed in the Users List and save the account (i.e., **Save and Back** and **Save As** buttons are enabled).
- Read** Allows users to *view* all accounts available in the Users List including their assigned roles and to select a user (i.e., **View** button is enabled) to display account details. Passwords will *never* be displayed. *This privilege must be granted to all users needing to view the **Manage** tab > Users option regardless of any other Users privileges assigned.*
- Update** Allows users to modify an account by assigning a different role, disabling/enabling the account, editing personal information (*First Name*, *Last Name*, and *Title*) and to save their changes (i.e., **Save and Back** button is enabled). *Although the username for an account cannot be changed, users with*

*the User: Create, Delete privileges can easily delete accounts and create new ones as necessary. Also enables the **Enable**, **Disable**, and **Reset Password** buttons in the **Manage tab > Users** option.*

Delete Allows users to select an account available in the Users List and delete it (i.e., **Delete** button is enabled) as necessary.

ROLES PRIVILEGES

The following privileges control whether users can access the **Manage tab > Roles** option allowing them to create and manage Matrix software roles.

Note: The *Roles: Read* privilege must be selected for the **Manage tab > Roles** option to be accessible to a user regardless of any other *Roles* privileges assigned to that user.



Create Allows users to *create* a new role (i.e., **Create** button is enabled) to be displayed in the Roles List and save the role (i.e., **Save and Back** and **Save As** buttons are enabled).

Read Allows users to *view* all roles available in the Roles List. *This privilege must be granted to all users needing to view the **Manage tab > Roles** option regardless of any other Roles privileges assigned.*

Update Allows users to modify a role by editing the role name, changing privileges selected and to save their changes (i.e., **Save and Back** and **Save As** buttons are enabled). *Renaming a role automatically updates the name of the role for all users to which the role has been assigned.* Also enables the **Rename** and **View** buttons in the **Manage tab > Roles** option.

Delete Allows users to select roles available in the Roles List and delete them (i.e., **Delete** button is enabled) as necessary.

AUDITS PRIVILEGES

The following privileges control whether users can access the **Manage tab > Audit Trail** option allowing them to search for and monitor the actions of user accounts, and to confirm the *Checksum Valid* column contains a checkmark (✓) for all actions indicating records have not been copied or falsified in any manner.

Note: The *Audits: Read* privilege must be selected for the **Manage tab > Audit Trail** option to be accessible to a user regardless of any other *Audits* privileges assigned to that user.



Read Allows users to *view* the automatically generated audit log captured by the system and displayed for the *Audit Trail* option under the **Manage tab**. *This privilege must be granted to all users needing to view the **Manage tab > Audit Trail** option regardless of any other Audits privileges assigned.*

Export Allows users to highlight one or more lines in the audit log and export them (i.e., **Export** button is enabled) to a user-defined location such as the Operating Computer desktop. You must use the Validator application that comes with the Matrix software to view the captured data.

FIRST-TIME LOGIN AS DEFAULT ADMIN USER

When logging in to Matrix software for the first-time *after* the *21 CFR Part 11* module has been enabled, you must log in using Default Administrator User credentials to create an Administrator User. Follow the steps listed below.

Note: If first-time login has already occurred on your system, you will need access to Administrator User credentials to create a *new* Admin User. Skip to *Creating an Administrator User* on page 94.

1. From the desktop of the Operating Computer, double-click the **Matrix** icon to launch the software.
2. Log in as the Default Administrator User by entering the username **admin** and password **admin**, then click the **Login** button. *You will immediately be prompted to create a new Administrator User.*
3. Respond to the *Create New User* dialog by accepting the **Matrix**  user type option selected by default (i.e., creates an Administrator User to reside locally on Operating Computer where Matrix software is installed) and click the **Continue** button.

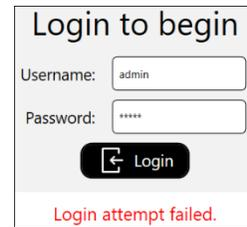


Note: If your system is connected to a network, you can create a *Windows* Administrator User by clicking the **Windows** user type and selecting a user account in your domain to be assigned the *Administrator* role.

4. In the *Account Information* area, enter a username and password for the new user. The *Administrator* role and enabled value of **Yes** will be pre-selected by default. *For Windows users, this area populates with existing username/password.*
5. In the *Personal Information* area, enter a first name, last name and title for your user. *For Windows users, this area populates with existing First Name/Last Name.*
6. Click **Save and Back** to create the user. You will automatically be logged out of the software upon creation of the new user.

- To confirm the Default Administrator User was disabled upon creation of your new Administrator User, attempt to log in again by entering username of **admin** and password of **admin**, then clicking **Login**. *An error message should be displayed.*
- Log in as your new Administrator User.

Continue with *Managing Roles* on page 95.



CREATING AN ADMINISTRATOR USER

If first-time login has already occurred on your system, you must log in using Administrator User credentials (i.e., user with *Administrator* role assigned) to create a *new* Administrator User. Follow the steps listed below.

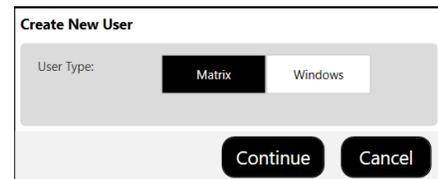
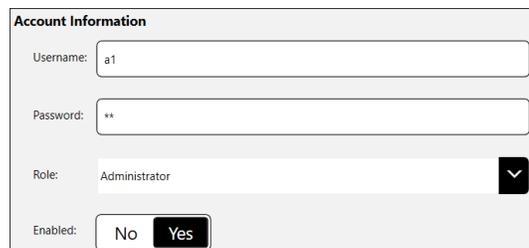
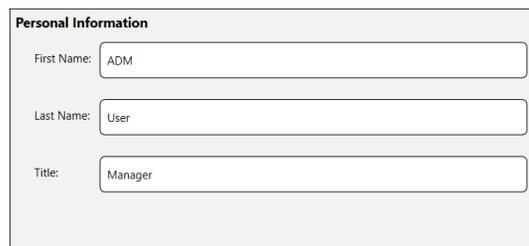
Note: If you have just created an Administrator User using the *First-Time Login as Default Admin User* section, skip to *Managing Roles* on page 95.

- From the desktop of the Operating Computer, double-click the **Matrix** icon to launch the software.
- Enter the username and password for an Administrator User already created by your organization, then click the **Login** button.
- Click the **Manage** tab in the Navigation Bar on left side of the screen and select the *Users* option.
- Click the **Create** button located at bottom of the screen.
- Respond to the *Create New User* dialog by accepting the **Matrix** user type option selected by default (i.e., creates an Administrator User to reside locally on the Operating Computer where Matrix software is installed) and click the **Continue** button.

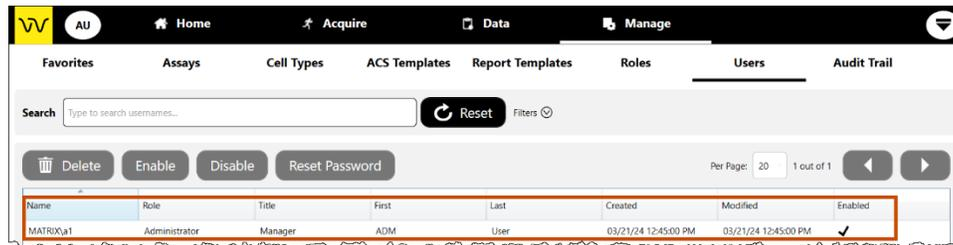



Note: If your system is connected to a network, you can create a *Windows* Administrator User by clicking the **Windows** user type and selecting a user account in your domain to be assigned the *Administrator* role.

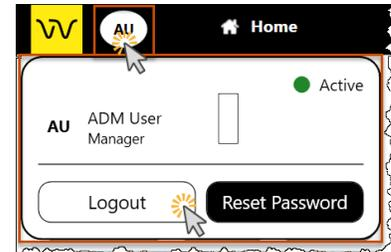
- In the *Account Information* area, enter a username and password for the user. Select the *Administrator* role from the dropdown and the enabled value of **Yes** (default value). *For Windows users, this area populates with existing Username/Password.*
- In the *Personal Information* area, enter a first name, last name and title for the user. *For Windows users, this area populates with existing First Name/Last Name.*
- Click **Save and Back** to create the Administrator User.

- Confirm the new user created appears in the *Users List*.



- Click the Administrator User ID (e.g., **AU**) in the Navigation Bar to open the User Card and select the **Logout** button.



MANAGING ROLES

Roles allow you to control varying levels of Matrix functionality to which specific groups of users will have access. As roles are created, they are added to the *Roles List* which displays a summary of privileges assigned to the role.

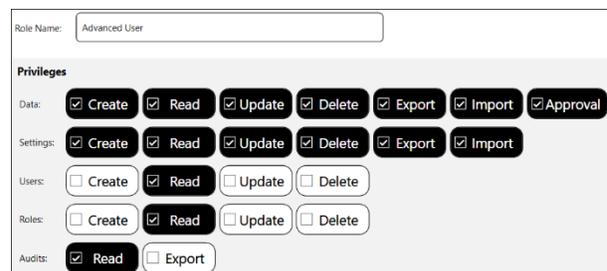
If your user account is associated with *Roles* privileges (i.e., *Roles: Create, Read, Update, and Delete*), you can create, maintain, and delete roles containing subsets of functionality to be assigned to users accordingly. *Keep in mind that the 'Administrator' role is a reserved role including full privileges and will not be displayed in the Roles List.*

To search for a role, enter a few key characters of a role name in the **Search** field and/or expand the **Filters** area to specify search criteria (e.g., *Created Between* or *Modified Between* start/end dates). As you enter criteria, the *Roles List* will automatically be updated to display matching entries.

To select a role and view its details, double-click a role entry in the *Roles List*, or click the role once to highlight it and then click the **View** button located at the bottom of the screen.

Creating a Role

- Click the **Manage** tab in the Navigation Bar on left side of the screen and select the *Roles* option.
- Click the **Create** button located at bottom of the screen.
- Enter a role name (e.g., *Advanced User*) and select privileges to be associated with the role (e.g., **Data: Create, Read; Settings: Read**).
- Click the **Save and Back** button and verify that the created role appears in the *Roles List*.



Note: When a role is selected in the *Roles List*, the **Rename**, **Delete**, and **View** buttons become enabled for that role.

Copying a Role

1. To create a role based on an existing role, double-click the role just created in the *Roles List*, or click the role once to highlight it and then click the **View** button.
2. Add and/or remove privileges, as necessary.
3. Click the **Save As** button.
4. In the dialog box displayed, you can either enter a new role name (e.g., *Technician II*) or select a name from the dropdown to override an existing role and click the **Save** button. Verify that the new role appears in the *Roles List*.

Renaming a Role

1. To rename a role, click the role once to highlight it and then click the **Rename** button.
Note: If you rename a role currently assigned to a user, the role name will be automatically updated.
2. Enter a new role name in dialog and click **Rename**. Verify that role name has been modified in the *Roles List*.

Deleting a Role

1. To delete a role, click the role once to highlight it and then click the **Delete** button.
Note: You cannot delete a role currently assigned to a user.
2. Verify that the role name has been removed from the *Roles List*.

Guidelines for Managing Roles

The following guidelines apply when managing *21 CFR Part 11* module roles.

- The *Administrator* role (i.e., containing *all* privileges) is not displayed in the *Roles List* as it cannot be modified. In addition, you cannot create a role with the name “administrator” as it represents a locked role.
- If a user is granted access to create/update roles (i.e., *Roles: Create, Read, and/or Update*), they will have access to assigning *ALL* privileges regardless of those assigned to their own role.
- If a user is granted access to create/update users (*Users: Create, Read, and/or Update*), they will have access to assigning *ALL* roles to those users regardless of the role assigned to their own user account. *Roles will be displayed in the Roles dropdown even if the user does not have Roles: Create, Read and/or Update privileges.*
- You cannot delete a role currently assigned to a user. *You must reassign another role to the user before you can delete the role.*
- If you rename a role while it is currently assigned to a user, the role name will automatically be updated for all users to which the role is assigned.
- If a role has the *Update* privilege (i.e., *Settings: Update*) for assays, cell types, and report templates, users assigned to this role will only be able to update these entities if the selected assay, cell type, or report template is unlocked (i.e., the *Unlocked* icon is displayed for the entity).

MANAGING USERS

Users are validated upon login and their actions are tracked as they move through the system via an *Audit Trail* log. As users are created, they are added to the *Users List* which displays a summary of account details including their assigned role, personal information (*Title* and *First Name/Last Name*), date/time stamp for when the account was created and last modified, and if the account is currently enabled.

Note: Only one user may be logged into the Matrix database at a time. However, users can export scan results and upload them as files onto an alternate computer for re-analysis of captured images via *Data Analysis* mode. See *Data Analysis Mode* on page 2 for details.

If your user account is associated with *Users* privileges (i.e., *Users: Create, Read, Update, and Delete*), you can create, maintain, and delete users whose access to Matrix functionality will be controlled by their assigned roles. *Keep in mind that you must always have at least one active Administrator User for the 21 CFR Part 11 module.*

To search for a user, enter a few key characters of a user name in the **Search** field and/or expand the *Filters* area to specify search criteria (e.g., *Roles, Enabled* status, or *Created Between* and *Modified Between* start/end dates). As you enter criteria, the *Users List* is updated to display matching entries.

Name	Role	Title	First	Last	Enabled
MATRIX\A1	Administrator	Manager	ADM	User	✓
MATRIX\admin	Administrator	IT User	Admin	User	✓
MATRIX\Advanced	Advanced User	Scientist	Advanced	User	✓
MATRIX\Basic	Basic User	Technician	Basic	User	✓

To select a user and view its details, double-click a user entry in the *Users List*, or click the user once to highlight it and then click the **View** button located at the bottom of the screen.

Creating a User

1. Create a new user by selecting the **Manage** tab > **Users** option.
2. Click the **Create** button located at the bottom of the screen.
3. Respond to the *Create New User* dialog by accepting the **Matrix**  user type option selected by default (i.e., creates a user to reside locally on the Operating Computer where Matrix software is installed) and click the **Continue** button.
4. In the *Account Information* area, enter a username and password for the user. *For Windows users, this area populates with existing Username/Password.*

Note: If your system is connected to a network, you can create a *Windows* user by clicking the **Windows** user type and selecting a user account in your domain.

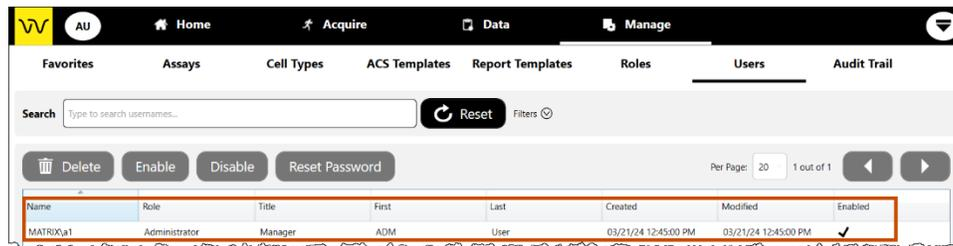
Note: Usernames must consist of at least one character and passwords are case sensitive. Once a user account is created,



The image shows a dialog box titled "Create New User". It has a "User Type:" label followed by two buttons: "Matrix" (which is highlighted with a dark background) and "Windows". At the bottom of the dialog, there are two buttons: "Continue" and "Cancel".

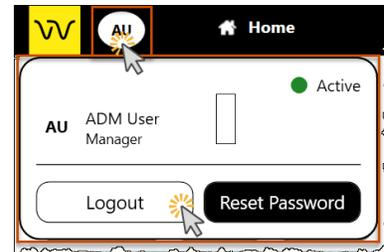
usernames and passwords cannot be changed (i.e., use the **Reset Password** button for *Matrix* users).

5. Select a role from the **Role** dropdown and ensure that the **Enabled** field is set to Yes.
6. In the *Personal Information* area, enter a first name, last name and title for the user. *For Windows users, this area populates with existing First Name/Last Name.*
7. Click the **Save and Back** button to create the user.
8. Confirm the new user appears in the *Users List*.



Note: When a user is selected in the *Users List*, the **Delete**, **Enable/Disable** (based on user status as displayed in the *Enabled* column), **Reset Password** and **View** buttons become enabled for that user.

11. Click the Administrator User ID (e.g., ) in the Navigation Bar to open the User Card and select the **Logout** button.



Disabling/Enabling a User

1. To disable a user, click the user once in the *Users List* to highlight it and then click the **Disable** button.
2. Click **Yes** in response to the confirmation prompt, followed by **OK** to acknowledge the user was disabled. Verify the checkmark in the *Enabled* column is no longer displayed for the user.

Note: Once a user is disabled, they will *not* be able to log in.

3. To enable a user, click the user once in the *Users List* to highlight it and then click the **Enable** button.
4. Click **Yes** in response to the confirmation prompt, followed by **OK** to acknowledge the user was enabled. Verify the checkmark in the *Enabled* column is displayed for the user.

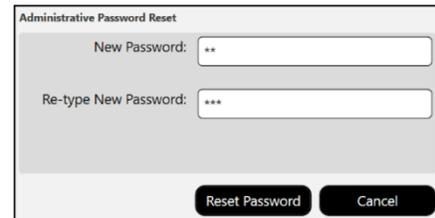
Resetting a User Password

Resetting a user password via an *Administrative Password Reset* is temporary and must be manually relayed to that user.

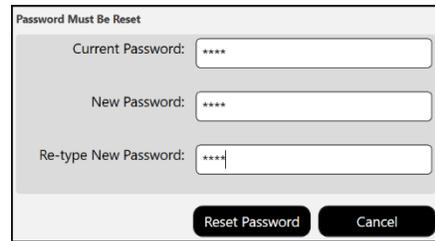
Note: You cannot reset the password for a *Windows* user as this functionality is controlled by mechanisms available in the network.

1. To reset a user password, click the user entry once in the *Users List* to highlight it and then click the **Reset Password** button.
2. Click **Yes** in response to the confirmation prompt.
3. In the *Administrative Password Reset* dialog, enter a new password for the user and confirm the password by entering it again. Click the **Reset Password** button.

Upon the user's next login with the temporary password, they will immediately be prompted to personally reset their password.



The dialog box titled "Administrative Password Reset" contains two text input fields. The first is labeled "New Password:" and has two asterisks (**) as a placeholder. The second is labeled "Re-type New Password:" and has three asterisks (***) as a placeholder. At the bottom right, there are two buttons: "Reset Password" and "Cancel".



The dialog box titled "Password Must Be Reset" contains three text input fields. The first is labeled "Current Password:" and has four asterisks (****) as a placeholder. The second is labeled "New Password:" and has four asterisks (****) as a placeholder. The third is labeled "Re-type New Password:" and has three asterisks (***) as a placeholder. At the bottom right, there are two buttons: "Reset Password" and "Cancel".

Deleting a User

1. To delete a user, click the user entry once to highlight it and then click the **Delete** button.
2. Click **OK** to acknowledge the user was deleted. Verify the user has been removed from the *Users List*.

Guidelines for Managing Users

The following guidelines apply when managing *21 CFR Part 11* module users.

- Your first-time login to Matrix software *after* this module has been enabled will be as Default Administrator using credentials provided. Upon login, you will be prompted to create a secure Administrator User and as part of its creation you will be logged out of the default account. *Default Administrator credentials will be deactivated at this time.* Log in as your newly created Administrator User to establish your user hierarchy.
- Administrator usernames cannot include "admin" or "administrator" as these represent reserved accounts.
- You must always have *at least one* active Administrator User for the *21 CFR Part 11* module.
- You cannot edit or delete the user account in which you are currently logged in to the system. If you disable an account while you are currently logged into it, you will *not* be able to log in using that account again. *An exception to this rule exists if you are logged in as the only remaining Administrator User; in this case, you will NOT be allowed to disable the account.*
- When clicking the **Reset Password** button for a user, you are performing an administrative password reset which results in the creation of a temporary password that must be manually relayed to the user. Upon the user's next login with the temporary password, they will immediately be prompted to reset the password.
- For network users, the **Reset Password** button cannot be selected as this functionality is controlled by mechanisms available in the domain.

Note: Only one user may be logged into the Matrix database at a time. However, users can export scan results and upload them as files onto an alternate computer for re-analysis of captured images via *Data Analysis* mode. See *Data Analysis Mode* on page 2 for details.

E-SIGNING COUNT RESULTS

Electronic Signatures capture the name of authorized users, the date/time when the signature was executed, and meaning of the signature (i.e., indicating it was for the *Review*, *Approval*, or *Rejection* of an action). Only users with the *Approval* privilege can electronically sign and un-sign count results.

1. Log in as a user with a basic role (e.g., **Data: Create, Read; Settings: Read**) and use the **Acquire** tab to enter setup details and perform a count. Note that *Status* for count result is *Not Signed* and **e-Sign** button is enabled.
2. Click the **e-Sign** button. Confirm username displayed, enter a password and reason for the action, and click the **e-Sign** button again.

Note: If the user performing a count does *not* have the *Approval* privilege in their assigned role, the *User does not have Approval permissions* message is displayed.
3. To log in as a supervisor and sign the count result on behalf of another user, enter username and password for a user with a role that includes the *Approval* privilege (e.g., **Data: Create, Read, Approval; Settings: Create, Read**), add a reason and click the **e-Sign** button. Note that *Status* displayed is changed to *Signed* and **e-Unsign** button is enabled.
4. To un-sign the count result, click the **e-Unsign** button, enter username and password for a user with a role that includes the *Approval* privilege, add a reason and click the **e-Unsign** button again. Note that *Status* displayed is changed to *Unsigned* and **e-Sign** button is enabled.

Status: Not Signed e-Sign

Sign Result

Username: MATRIX\Basic

Password: *****

Reason: Valid Result

User does not have Approval permissions

e-Sign Cancel

Status: Signed - Advanced User (MATRIX\Advanced) 2024/03/2 e-Unsign

Unsign Result

Username: MATRIX\Advanced

Password: *****

Reason: Unsign Test

e-Unsign Cancel

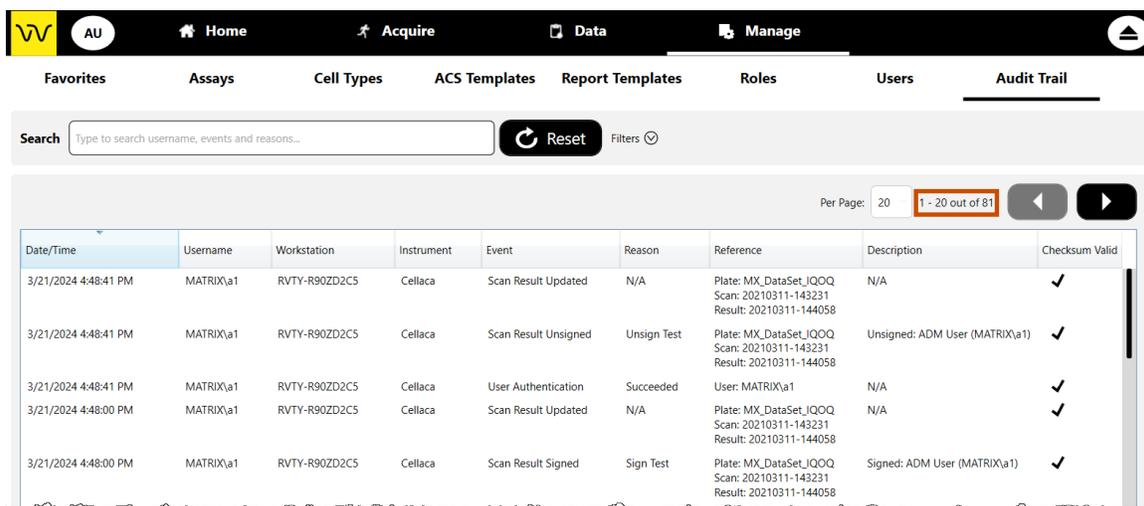
Status: Unsigned - Advanced User (MATRIX\Advanced) 2024/03 e-Sign

VIEWING AUDIT TRAIL

The *Audit Trail Log* provides assurance regarding the integrity of records, and must capture the identity of all users performing an action, the type of action, and the date/time associated with the action. Audit trails can be retained locally on the Operating Computer or on a network.

Searching the Log

1. Click the **Manage** tab > **Audit Trail** option.
2. To search for the actions of specific users, enter a few characters of a user name in the **Search** field and/or expand the *Filters* area to specify search criteria (e.g., *Usernames*, *Events*, or *Created Between* start/end dates). As you enter criteria, the *Audit Trail* is updated to display matching entries.



The screenshot shows the 'Audit Trail' section of a software interface. At the top, there is a navigation bar with tabs for 'Home', 'Acquire', 'Data', and 'Manage'. Below this, there are sub-tabs for 'Favorites', 'Assays', 'Cell Types', 'ACS Templates', 'Report Templates', 'Roles', 'Users', and 'Audit Trail'. A search bar is present with the placeholder text 'Type to search username, events and reasons...'. Below the search bar, there is a 'Reset' button and a 'Filters' dropdown. The main area displays a table of audit trail entries. The table has columns for Date/Time, Username, Workstation, Instrument, Event, Reason, Reference, Description, and Checksum Valid. The 'Checksum Valid' column contains checkmarks for each entry. The table shows five entries, all with a 'Checksum Valid' status of '✓'.

Date/Time	Username	Workstation	Instrument	Event	Reason	Reference	Description	Checksum Valid
3/21/2024 4:48:41 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Updated	N/A	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	N/A	✓
3/21/2024 4:48:41 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Unsigned	Unsign Test	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	Unsigned: ADM User (MATRIX\A1)	✓
3/21/2024 4:48:41 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	User Authentication	Succeeded	User: MATRIX\A1	N/A	✓
3/21/2024 4:48:00 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Updated	N/A	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	N/A	✓
3/21/2024 4:48:00 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Signed	Sign Test	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	Signed: ADM User (MATRIX\A1)	✓

3. Review matching entries and ensure that a checkmark appears in the *Checksum Valid* column for each entry.

Exporting Log Details

1. Click the **Export** button, navigate to an export location on the desktop and click **OK** to save the audit trail log. A *YYYYMMDD-HHMMSS-AuditTrail-MATRIX-<user>.AUDITTRAIL* file is created, where *YYYYMMDD* represents the current date, *HHMMSS* is a time stamp (hours, minutes, seconds) and *<user>* is your username.

Note: When exporting audit log details, it may be useful to record the total number of action entries in the log prior to export (e.g., 1354 as shown in the example below).



The screenshot shows the 'Audit Trail' section of the software interface. At the top, there is a navigation bar with tabs for 'Home', 'Acquire', 'Data', and 'Manage'. Below this, there are sub-tabs for 'Favorites', 'Assays', 'Cell Types', 'ACS Templates', 'Report Templates', 'Roles', 'Users', and 'Audit Trail'. A search bar is present with the placeholder text 'Type to search username, events and reasons...'. Below the search bar, there is a 'Reset' button and a 'Filters' dropdown. The main area displays a table of audit trail entries. The table has columns for Date/Time, Username, Workstation, Instrument, Event, Reason, Reference, Description, and Checksum Valid. The 'Checksum Valid' column contains checkmarks for each entry. The table shows five entries, all with a 'Checksum Valid' status of '✓'.

Date/Time	Username	Workstation	Instrument	Event	Reason	Reference	Description	Checksum Valid
3/21/2024 4:48:41 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Updated	N/A	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	N/A	✓
3/21/2024 4:48:41 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Unsigned	Unsign Test	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	Unsigned: ADM User (MATRIX\A1)	✓
3/21/2024 4:48:41 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	User Authentication	Succeeded	User: MATRIX\A1	N/A	✓
3/21/2024 4:48:00 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Updated	N/A	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	N/A	✓
3/21/2024 4:48:00 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Signed	Sign Test	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	Signed: ADM User (MATRIX\A1)	✓

2. From the Windows Applications Menu expand the **Revvity** folder and launch the *Matrix Validator* application.

- Click the **Browse** button to navigate to the exported `YYYYMMDD-HHMMSS-AuditTrail-MATRIX-<user>.AUDITTRAIL` file in the saved location.



- Double-click the file or click the file once to select it and then click **Open** to view the exported audit log.

Date/Time	Username	Workstation	Instrument	Event	Reason	Reference	Description	Checksum Valid
3/21/2024 4:48:41 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Updated	N/A	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	N/A	✓
3/21/2024 4:48:41 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Unsigned	Unsign Test	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	Unsigned: ADM User (MATRIX\A1)	✓
3/21/2024 4:48:41 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	User Authentication	Succeeded	User: MATRIX\A1	N/A	✓
3/21/2024 4:48:00 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Updated	N/A	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	N/A	✓
3/21/2024 4:48:00 PM	MATRIX\A1	RVTY-R90ZD2C5	Cellaca	Scan Result Signed	Sign Test	Plate: MX_DataSet_IQOQ Scan: 20210311-143231 Result: 20210311-144058	Signed: ADM User (MATRIX\A1)	✓

- Validate that the total number of action entries currently logged for your system (i.e., *1354* as highlighted in the sample screen shown above) matches what you recorded before the export.
- To search for the actions of specific users, enter a few characters of a user name in the **Search** field and/or expand the *Filters* area to specify search criteria (e.g., *Usernames, Events, or Created Between* start/end dates). As you enter criteria, the *Audit Trail* is updated to display matching entries.

Chapter 12. Contacting Support

This chapter presents the scope of Support services and provides contact methods. In addition, it contains instructions on how to report issues to Support and generate diagnostic reports to assist with troubleshooting.

SCOPE OF SUPPORT SERVICES

Revvity is dedicated to providing our customers with outstanding support including the following services:

- Online and in-lab customer training
- Creation of new cell types
- Optimization of counting parameters
- Creation of new report templates
- Troubleshooting via telephone
- Periodic safety checks and functional evaluations (offered as part of a separate maintenance contract)

To inquire about training, visit our website at <https://www.revvity.com/contact-us/customer-training> and choose the *Cell Counting and Image Cytometry* product. Enter your contact details and any training comments/questions.

CONTACT METHODS

If there is a technical issue with your instrument or software, contact Support using the following methods:

- Visit <https://www.revvity.com/contact-us/instrument-support-and-service> and choose the *Cell Counting and Image Cytometry* product for support. Enter instrument serial number, name and model, your contact details, and a detailed description to report the issue to Support.
- Visit <https://www.revvity.com/contact-us-by-phone> to find the global phone number for your area.
- Send email to CellC-support@revvity.com

REPORTING AN ISSUE TO SUPPORT

If a technical issue encountered cannot be resolved using troubleshooting steps provided in your instrument's user manual or the issue persists after rebooting the instrument, perform the following steps *before* contacting Support to report the issue:

1. Record the error message.
2. Record the sequence of events leading up to the error, if possible.
3. If necessary, close the error message window.
4. Record the Serial Number located on the Device Label for your instrument.

Gathering these details *prior* to contacting Support will be helpful as they troubleshoot the technical issue.

GENERATING DIAGNOSTIC REPORTS

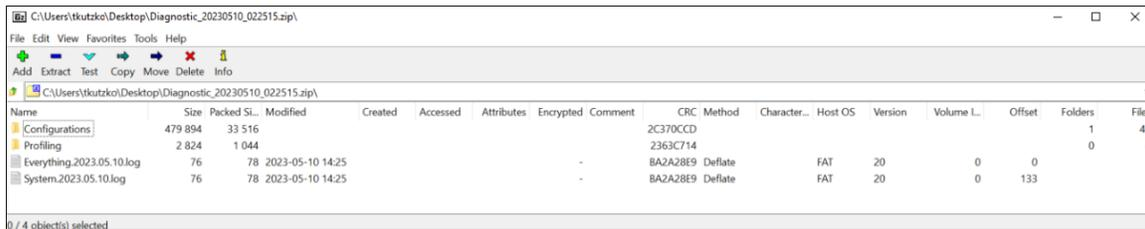
To generate a diagnostic report that can be emailed to Support when experiencing a technical issue:

1. Click the **Home** tab and then the **About Matrix** button.
2. Click the **Generate Diagnostic Report** button followed by **OK** in response to the confirmation prompt.
3. From the desktop, click the generated **Diagnostic_YYYYMMDD.zip** folder (where YYYYMMDD represents the date on which file was generated) to display files in the folder.

Generate Diagnostic Report

Successfully generated the diagnostic report on the desktop.

OK



Files in the zipped folder include logs (located in `C:\logs\Matrix\vNNN` where *NNN* represents installed version) and configuration files (located in `C:\ProgramData\Revvity\Matrix\vNNN\Configurations`).

4. Attach the zipped folder to an email, include the *Support Ticket ID* (if assigned) in the Subject line and send to: CellC-support@revvity.com

It may be helpful to clear all logs before reproducing the sequence of steps leading up to an issue and/or to reduce the size of the diagnostic report to be sent to Support. *Logs are generated automatically by your system on a daily basis. Keep in mind that clearing all logs will remove accumulated files to date for the version.*



WARNING: Logs may be required to maintain a historical archive. As using the Clear All Logs button will *permanently* remove accumulated logs for the installed Matrix software version, it is recommended that you contact IT *before* clearing logs from your system.

To clear all logs:

1. Click the **Clear All Logs** button.

Note: It is *not* recommended to clear all logs unless you are confident that they are not being archived by your organization.
2. When prompted, click **Yes** to confirm you want to clear all logs.

Clear All Logs

Are you sure you want to clear all log files?

Yes No

Click **OK** to acknowledge that all logs have been cleared.

Successfully cleared all log files.

OK

Appendix A. Report Designer for WPF Reference

Revvity uses *DevExpress* as a third-party plugin for Matrix software reporting capabilities.

Visit the following page on the DevExpress website for more information about creating Matrix software report templates using *Report Designer for WPF*.

<https://devexpress.github.io/dotnet-eud/reporting-for-desktop/articles/report-designer/report-designer-for-wpf.html>

Sample screens displaying the range of topics available for *Report Designer for WPF* functionality are shown below.

The screenshot displays the DevExpress Report Designer for WPF application. The interface includes a top navigation bar with tabs for 'Dashboard for Desktop', 'Dashboard for Web', 'Interface Elements for Desktop', and 'Interface Elements for Web'. Below this is a breadcrumb trail: 'Interface Elements for Desktop / Report Designer / Report Designer for WPF'. A search bar is located at the top left of the main content area.

On the left side, there is a vertical navigation menu with the following items:

- + Charting
- + Docking
- + Editors
- + Expression Editor
- + Filter Editor
- + Grid
- + Layout Manager
- + Map
- + Navigation Bars
- + PDF Viewer
- + Pivot Table
- + Print Preview
- Report Designer
 - + Report Designer for WinForms
 - Report Designer for WPF
 - + Report Types
 - + Creating Reports
 - + Report Elements
 - + Interface Elements
 - + Report Wizard
 - + Document Preview
- + Ribbon
- + Rich Text Editor

The main content area features a large heading 'Report Designer for WPF' and a sub-heading 'This guide contains information about the basic principles of creating reports with the Report Designer.' Below this, a paragraph states: 'The Report Designer allows you to create new reports from scratch, bind them to data and fully customize them. In addition to report editing capabilities, it allows you to display a report's Print Preview, send its outputs to a printer or export it to various formats.'

The central part of the screenshot shows a preview of a report titled 'Suppliers'. The report has a header section with the title 'Suppliers' and a table with the following columns: 'Company', 'Contact Name', 'Contact Title', 'Phone', 'Fax', 'Home Page', 'Address', 'Country', 'Region', 'City', and 'Postal Code'. The table is currently empty. Below the table, there is a 'Group And Sort' section with options to 'Add a Group', 'Add a Sort', 'Remove', 'Move Up', and 'Move Down'. The 'Field' column is empty, and the 'Sort Order' and 'Show Header' columns are also empty.

On the right side, there is a 'Report Explorer' pane showing a tree view of the report structure, including 'Suppliers', 'topMarginBand1', 'ReportHeader1', 'Detail', 'DetailReportBand', and 'BottomMargin'. Below the tree view is a 'Properties' pane for the selected 'Suppliers Report' object, showing various settings such as 'Data Source', 'Designer Options', 'Display Name', and 'Draw the Grid'.

Dashboard for Desktop Dashboard for Web **Interface Elements for Desktop** Interface Elements for Web

Interface Elements for Desktop / Report Designer / Report Designer for WPF

Enter here to filter...

- + Charting
- + Docking
- + Editors
- + Expression Editor
- + Filter Editor
- + Grid
- + Layout Manager
- + Map
- + Navigation Bars
- + PDF Viewer
- + Pivot Table
- + Print Preview
- **Report Designer**
 - + Report Designer for WinForms
 - Report Designer for WPF

Different aspects of using the Report Designer are covered in the following documentation sections.

- [Creating Reports](#)

The tutorials in this section provide step-by-step instructions on both basic and advanced report customization.
- [Report Types](#)

The documents in this section describe how to create reports of different types with the Report Designer.
- [Report Elements](#)

The topics in this section provide information about report controls and bands used in the Report Designer.
- [Interface Elements](#)

The documents in this section are dedicated to the elements of the Report Designer user interface.
- [Report Wizard](#)

This documentation section describes the Report Wizard, which allows you to create reports based on built-in templates.
- [Document Preview](#)

The topics in this section describe the capabilities provided by the Print Preview.

Appendix B. Software License Details

This appendix presents Revvity's *Terms and Conditions* related to the use of the Matrix software. In addition, it includes a definition of *Revvity Proprietary Information*.

TERMS AND CONDITIONS

The *Revvity, Inc. – Terms and Conditions of Sale* license agreement states the terms and conditions upon which Revvity offers to license to you the software together with all related documentation. The Matrix software is licensed to you for use only in conjunction with Revvity's family of products.

In addition, the original Matrix software and any subsequent software upgrades installed for your Revvity instrument system is protected. You may not tamper with this software (including unauthorized upgrades), disclose it to third parties or use it for any purpose other than running your Revvity instrument system.

Revvity, Inc. does not grant you any other rights to use or disclose the original Matrix software or its upgrades, and any further uses will be prosecuted by Revvity to the maximum extent possible by law. Any other use of Matrix software or its upgrades is explicitly prohibited. In addition, you may not disclose Matrix software, upgrades, or any of its features and benefits to a third party.

Revvity Proprietary Information

Cellometer and Cellaca MX products have been developed by Revvity, Inc. and include certain intellectual property of Revvity, including without limitation, software, samples, schematics, specifications, manuals, designs, and other technical, business, trade secret, proprietary and confidential information provided to Buyer by Revvity ("Revvity Proprietary Information").

Buyer is granted a non-exclusive right and license to use the Revvity Proprietary Information solely: (a) as incorporated into, and in conjunction with, the products, (b) in conformance with the specifications, and (c) for Buyer's internal use.

Buyer may not: (i) assign, sublicense, transfer, lease, rent or distribute any of its rights in the Revvity Proprietary Information; (ii) port, translate, localize or create derivative works based upon the Revvity Proprietary Information in any manner; (iii) reverse assemble, decompile, reverse engineer, translate or otherwise attempt to derive or obtain the source code, the underlying ideas, algorithms, structure or organization of the Revvity Proprietary Information; (iv) use the Revvity Proprietary Information for the benefit of any third party including as part of any service bureau, time sharing or third party training arrangement; or (v) publish any benchmark testing results on any product or the Revvity Proprietary Information without Revvity's written consent.

Revvity, Inc. retains all ownership rights in the Revvity Proprietary Information and, other than limited license set forth in this section, Buyer shall have no right in or to the Revvity Proprietary Information.

Buyer will not disclose the Revvity Proprietary Information to any third party or use it in any manner outside the scope of the license including: (1) developing, designing, manufacturing, engineering, reverse engineering, refurbishing, selling or offering for sale items, parts or components of items, derivatives of or equivalents, or (2) assisting any third party in any manner to perform such activity.

Buyer shall use reasonable care to protect the Revvity Proprietary Information, and in no event less than the care Buyer uses to protect its own like information.



www.revivity.com

revvity



Revvity, Inc.
360 Merrimack Street, Suite 200
Lawrence, MA 01843 USA