

Contents of Shipping Container

- ❑ Cellaca PLX Instrument
- ❑ Instrument **Power Supply** and **Power Cord**
- ❑ **USB 3.0 Connector Cable**
- ❑ **PLX Slide Holder**
- ❑ **Revvity-provided Laptop** with **Power Supply** and **Power Cord**
- ❑ **Matrix Software** and De Novo Software **FCS Express** (both pre-installed on Laptop)
- ❑ **De Novo Software FCS Express License Packet with Dongle**
- ❑ **Cellaca PLX User Manual** (PDF file on Laptop)
- ❑ **Matrix Software User Manual** (PDF file on Laptop)
- ❑ **Cellaca PLX Quick Start Guide** (PDF file on Laptop – *this document*)
- ❑ **Cellaca MX/PLX Plate Loading Template** – Indicates mixing/loading wells
- ❑ **Cellaca MX/PLX Focus Guide** – Assists with adjusting instrument focus
- ❑ Three **Revvity Counting Plates (12x2 orientation, CHM24-A100)**
- ❑ Three **Revvity Counting Plates (3x8 orientation, CHM24-B100)**
- ❑ Ten **Cellaca PLX Low Fluorescence Slides (CHM2-ACR)**
- ❑ Two **AO/PI Viability Reagent Vials (CS2-0106-5mL)**

Unboxing the Instrument

The Cellaca MX and PLX instruments are part of the same product family and share many features, including being powered by the *Matrix Software*.

Powered By  **Matrix™**

Unpack and visually inspect the PLX to ensure no physical damage has occurred during shipping. For assistance in setting up the instrument, visit the Cellaca MX/PLX pages on our website for training videos on unboxing and getting started.



Site Preparation

Instrument must be placed on a level surface and plugged directly into a power outlet. *Use of a surge protector is recommended.* Ensure all cables are free from tangles prior to powering on the Cellaca PLX.

Follow all equipment safety protocols and keep the area around instrument clean both during and post operation. Do *not* position the device so that it is difficult to disconnect from power main.

Setting Up the System

The *Cellaca PLX Image Cytometry System* is comprised of the instrument connected via USB cable to an Operating Computer (Laptop) that is used to run the Matrix software. The Operating Computer can be linked to a network for accessing external files, printers and for storing scan results.

Connect instrument Power Supply/Power Cord to the Cellaca PLX and plug Power Cord into electrical outlet. Turn PLX Power Switch *ON* and confirm that light bar on front of instrument is lit. Set up and power on Revvity-provided Laptop, and connect to instrument using USB 3.0 Connector Cable.

Note: 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.*

From desktop of Operating Computer, double-click the **Matrix** icon to launch the software.



Calibrating Instrument Prior to First Use

After initial setup of an instrument, you must calibrate it using the Matrix software prior to first-time use. 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 plate or slide holder loaded in the instrument. Contact Support for assistance with calibration.


Matrix Screen Elements

Upon launching the Matrix Software you are presented with the **Acquire** tab > *Setup* option by default. Basic screen elements are described below.


Note: If the Matrix 21 CFR Part 11 module has been enabled for your system, users must log in *before* they can begin using the software.

The *Navigation Bar* visible across top of the screen is always displayed and contains workflow tabs. It also contains the **Eject/Load** buttons that control movement of the instrument stage (for unloading/loading consumables).

Functionality associated with each tab is described below.

 Home

Home Tab: Displays the Matrix logo distinguishing the software from prior versions. Also contains the **About Matrix** button which displays version details and Revvity contact information.

 Acquire

Acquire Tab: Contains the two sequential screens in the *Data Acquisition* workflow. Users can select a *Favorite* (with all assay parameter settings pre-defined) to streamline the workflow or enter setup details before advancing to **Preview** or **Count** (if *Skip Preview* feature is enabled) the sample.

 Data

Data Tab: Contains the three sequential screens in the *Data Analysis* workflow. Users must first *Select* a scan result before viewing its count *Results*. You can also perform a *Recount* by fine-tuning assay parameter settings to be used in the analysis to create a new scan result.


 Manage

Manage Tab: Contains the various system lists of *Favorites*, *Assays*, *Cell Types*, *ACS Templates* and *Report Templates* used during *Data Acquisition* and *Data Analysis* workflows. You can manage system lists by importing/exporting, renaming, deleting or showing/hiding individual entities. In addition, you can create new or modify existing favorites, assays, cell types, and report templates.


Preparing Plates/Slides with Samples

Although the Cellaca PLX does not require any routine testing or calibration, counting beads are available to verify instrument functionality. *Revvity counting beads CCBM-011-2ML are recommended for use with Cellaca PLX.*

To prepare plates/slides with counting beads or cell samples:


1. From desktop of Operating Computer, launch the Matrix software by double-clicking the **Matrix** icon. 
2. Invert counting bead solution or tube containing cell sample a total of 10 times (10x). *If using counting beads, vortex bead solution for 10 seconds. Do NOT vortex cell samples.*

To Load a Counting Plate

1. Place the *Cellaca MX/PLX Plate Loading Template* on the lab bench with the appropriate layout facing up and set a counting plate on top of the template graphic, aligning the notched corner of the plate. *You should be able to view and identify the mixing/loading wells through the plate.* 
2. Set pipette to 50 µL and then pipette bead solution/cell sample up and down 10x to break up potential clumps. Load 50 µL of bead solution/cell sample into each of the individual *Loading Wells* on the counting plate. *Liquid will move into imaging area for each filled well.*

If cell samples require mixing prior to loading, pipette samples into *Mixing Wells* on the counting plate (adjusting volume as necessary to account for dilution factor). Once mixing of samples is complete, pipette 50 µL from *Mixing Wells* into associated *Loading Wells*.


To Load Slides and Insert into Slide Holder

1. Place up to four PLX slides on unused/clean Kimwipes.
2. Since two samples can be loaded per slide, it is recommended that you label individual chambers in the blue area. Take care to ensure that clear optical windows of wells are *not* touched. 

- Set Pipette to 15 μL and then pipette bead solution/cell sample up and down 10x to break up potential clumps. Load 15 μL of bead solution/cell sample into induction port of a slide well. *Liquid will move into imaging area for each filled well.*
- Insert up to four prepared slides into slide holder, aligning well A of each slide with top of holder (indicated by notched corner).



Placing Plate/Slide Holder onto Instrument Stage

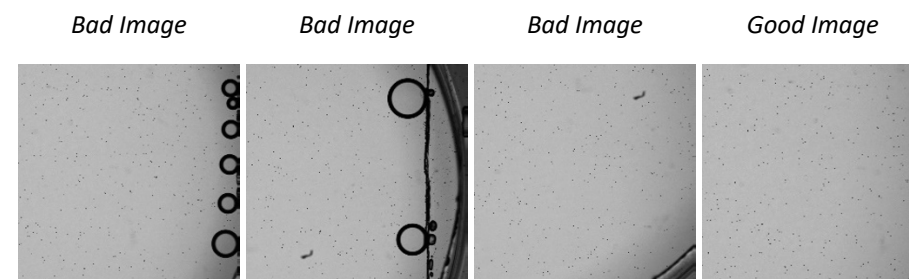
- In the Matrix software, click **Eject**  located in the Navigation Bar to eject stage from instrument.
- Place slide holder (containing up to 4 slides) or counting plate onto stage, taking care to align notched corner of holder/plate with top left corner of stage (i.e., well A1 is positioned in top left corner while facing stage).



- Click **Load**  to retract stage into instrument.

Performing Image Clipping Check

- Click the **Acquire** tab and select the *Brightfield Channel Verification* assay.
- Select a well in a corner of plate or slide holder and click **Preview**.
- Visually verify that no clipping has occurred around edges of the image. *Samples shown below display wells in a counting plate.*

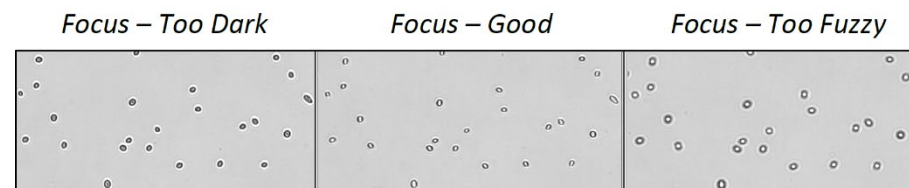


- Repeat this step for each well in outer corners of the plate/slide holder.

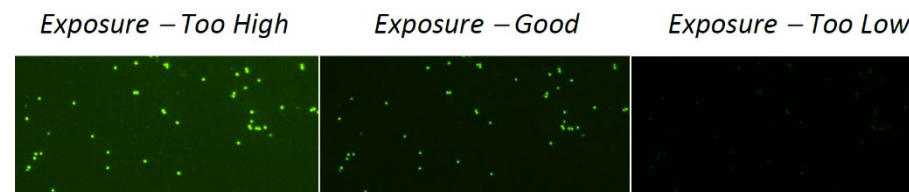
Understanding Instrument Focus

To adjust the focus of a live image being previewed, an **Auto Focus** button allows the instrument to determine the best focal position for selected well. *Focus* controls are also available if additional fine tuning is required.

Live cells should have a bright center and dark, crisp clearly defined edges.

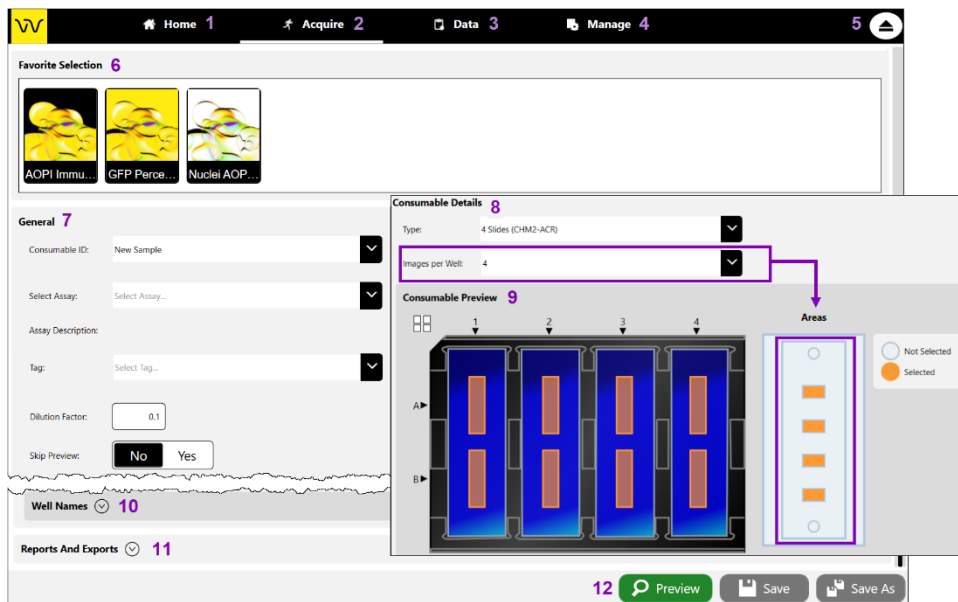


Fluorescent signals should be strong with a low, dark background.



Counting and Analysis Workflow

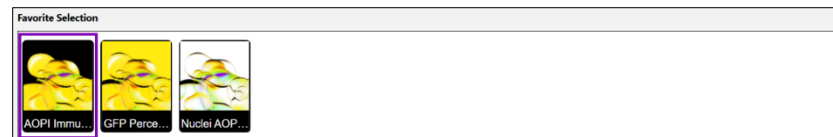
Performing a count consists of either *Selecting a Favorite* or *Entering Parameter Settings*. Upon launch of the software, the **Acquire** tab > *Setup* option is displayed representing the start of data acquisition workflow.



- 1 Home Tab** – Displays Home screen with software/instrument details.
- 2 Acquire Tab** – Use to select favorites and enter parameter settings.
- 3 Data Tab** – Use to view current scans/select saved scans for re-analysis.
- 4 Manage Tab** – Displays *Favorites, Assays, Cell Types, ACS Templates,* and *Report Templates* options for managing system lists.
- 5 Eject Button** – Controls instrument stage (toggles between **Eject/Load**).
- 6 Favorites Selection Area** – Displays favorites for quick re-use of assays.
- 7 General Area** – Use to enter consumable ID, select an assay, add a tag, specify a dilution factor or and choose whether to *Skip Preview* screen.
- 8 Consumable Details Area** – Use to enter *Type* and *Images Per Well*.
- 9 Consumable Preview Area** – Displays a Well Map for selecting wells.
- 10 Well Names Area** – Use to enter or import names for selected wells.
- 11 Reports and Exports Area** – Use to specify auto export location.
- 12 Preview Button** – If plate is loaded, *Setup* details are entered and a well is selected, displays live images to adjust focus/fluorescence exposure.


Selecting a Favorite

Scroll across the *Favorite Selection* panel to select a favorite. Simply click **Preview** (if enabled) or **Count** to proceed to the next step in the process.



Skip to *Previewing Live Images* on page 5 (if **Preview** button is enabled) or *Clicking the Count Button* on page 5.

Entering Parameter Settings

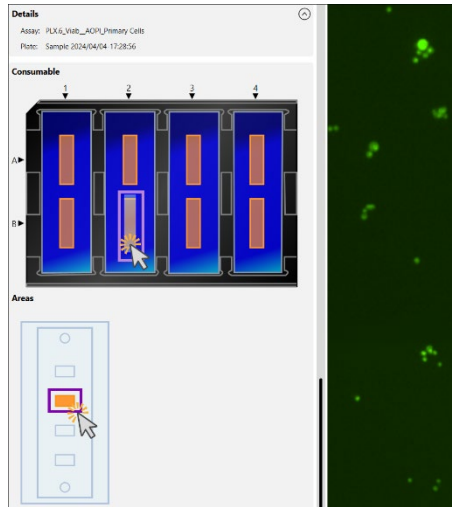
- In the **Acquire** tab, Setup screen, enter a **Consumable ID**. *If you do not enter a consumable ID, a date/time stamp will be appended to the “New Sample” default (e.g., New Sample 2024/03/25-11:58:09).*
- Select an assay from the dropdown. The **View** button is available if you find it necessary to edit assay details prior to previewing the live image.
- 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.
- Select consumable **Type** and **Images Per Well** to be taken in the scan. *Images Per Well indicator appears to the right of the Well Map.*
- In the *Consumable Preview* area, click on wells containing samples. To select individual wells, click to select or de-select accordingly. To select a block of wells, click on a well at the beginning of the block and hold button down while dragging mouse to end of the block before releasing button. To select or de-select *all* wells, click the **All Wells**  icon.
- If desired, expand the *Well Names* area to enter/import well names or the *Reports and Exports* area to modify the default auto export location.
- Click the **Preview** (if enabled) or **Count** button.

Previewing Live Images

The Preview screen displays live images for selected wells.

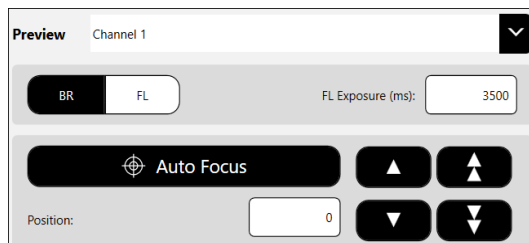
1. Expand the Details area to view selected assay and consumable ID.

2. In the *Consumable* area, click on highlighted wells to view live images of samples contained in the plate or slides. If an **Images Per Well** value other than 1 was selected, use the scroll bar to view this indicator and click on an area. As you move from well to well (and from quadrant to quadrant), the live image changes per selected well and quadrant.



To zoom in and out of an image, move the mouse to hover cursor over the viewing pane and turn the scroll wheel or, if using the touch screen, 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.) To move an image around, click on the image and drag to a new location.

3. Adjust focus of the image using focusing controls. *The Channel 1: Brightfield image is displayed by default.* Click the **Auto Focus** button or use the manual *Focus* controls if additional fine tuning is required.



Auto Focus Allows instrument to determine the best focal position for the selected well.

Focus Position Allows users to enter a numerical value for vertical (Z) position of the objective lens.

Fine Focus Manual Offset Allows users to finely adjust vertical (Z) position of the objective lens for optimal focus (in μm).



Click the up/down single arrows to adjust focus.

Coarse Focus Manual Offset Allows users to coarsely adjust vertical (Z) position of the objective lens for optimal focus (in μm).



Click the up/down double arrows to adjust focus.

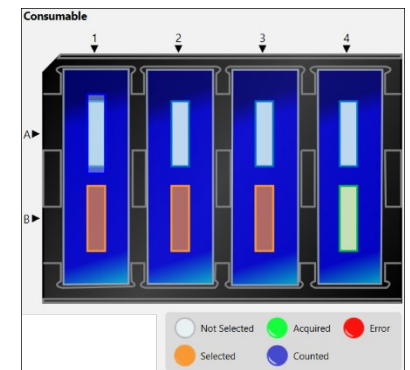
Note: For BR/FL imaging modes, each channel will be associated with two images – *Brightfield (BR)* and *Fluorescent (FL)*. When working with assays that have more than one channel, use the *Channel 1/BR* image to adjust focus and then select the *FL* image to confirm exposure. For *Channels 2-6*, you only need to select the *FL* images to confirm exposure since focus of their paired *BR* images is adjusted automatically when you performed the task for *Channel 1*.

- In the **Preview** field, click the dropdown to view available channels. *Channels displayed will depend on assay Imaging Mode.* Select a channel to change image display.
- For BR/FL imaging modes, click **BR** or **FL** buttons to change image display. *BR image will always be displayed by default.* Adjust **FL Exposure** value for FL images associated with each channel, if needed.

Clicking the Count Button

Once you have completed previewing the live image for the sample, click **Count**.

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

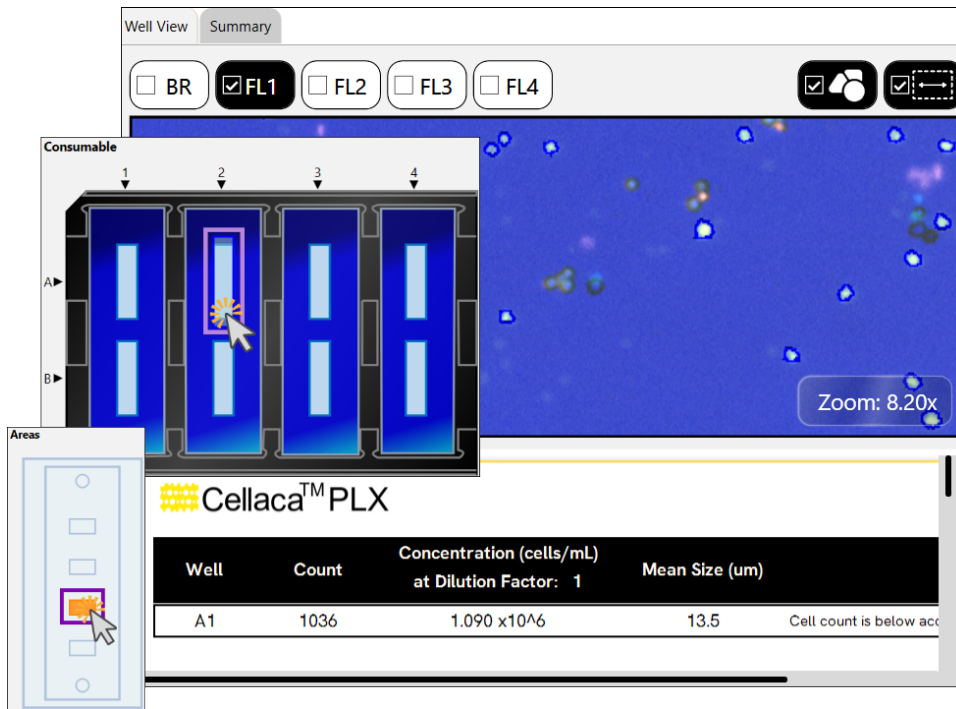


You can click on a well as soon as it is *Counted* to display its image (in the viewing pane) and its count results (below the viewing pane).

Well images are displayed in the **Well View** tab along with acquired data for the selected well. Additional tabs may be enabled for the scan result (e.g., **Summary** and **.csv**).



Analyzing Scan Results

Scan results are displayed (using assigned report tabs/templates), printed and exported based on *Reports and Exports* options defined for the assay.



When analyzing scan results, an image is displayed for the selected well. To view images from other wells, click on each well in the Well Map.

1. In *Consumable* area, click on a well to display its image in viewing pane.
2. At the top of the **Well View** tab, click the *Brightfield (BR)* or *Fluorescence* (e.g., **FL 1**, **FL2**, etc.) buttons to select/de-select channels used in display of the image. *Channel views are overlaid on top of each other.*

3. Click the **Counted Overlay** button  to show/hide a graphic overlay that highlights *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 specified cell diameter). For Expression assays, outlining is used for total cells in the masked channel but is only displayed in one color – Blue.*
4. Click the **Zoom** button  to enable/disable display of *Zoom* magnification in bottom right corner of viewing pane. *Zoom feature will still be functional even if not displayed.*
5. At the bottom of the **Well View** tab is a report containing well-level details for the sample image from the selected well.
6. Click the **Summary** tab (displayed for a Matrix v4.0+ assay) or **All Wells** tab (displayed for a v3.0 assay) to view a full page containing count results for all scanned wells.

Well	Count	Concentration (cells/mL) at Dilution Factor: 1	Mean Size (um)	
A1	1036	1.090 x10 ⁶	13.5	Cell count is below acc
A2	1036	1.090 x10 ⁶	13.5	Cell count is below acc
A3	1036	1.090 x10 ⁶	13.5	Cell count is below acc
A4	1036	1.090 x10 ⁶	13.5	Cell count is below acc
B1	1036	1.090 x10 ⁶	13.5	Cell count is below acc
B2	1036	1.090 x10 ⁶	13.5	Cell count is below acc
B3	1036	1.090 x10 ⁶	13.5	Cell count is below acc
B4	1036	1.090 x10 ⁶	13.5	Cell count is below acc

Verifying Auto Exports

Expand the *Exports* area in the Results screen to verify that automatic exports were completed and click the location link to view exported scan result files. If any output file types (e.g., *CSV*, *Excel*, *PDF* and *Word*) were selected to be generated for assay reports, verify that file type exports were also completed.

To modify default *Reports* and *Exports* options defined for an assay during a count/recount, click the **View** button associated with the selected assay and expand the *Reports and Exports* area for the assay. You can either save your changes to the assay (i.e., any changes will be applied to other scan results that use the assay) or save the changed assay as a copy with a new name and then select it to be used for data acquisition.

Exports				
Exports	Status			
Raw Images	✓			
Colorized Images	✓			
Well Level CSV	✓			
Object Level ACS	✓			
Object Level CSV	✓			
DataSet	✓			
Reports	CSV	Excel	PDF	Word
PLX_Surface Marker(s)	✓	✓	✓	✓

An alternative to modifying default *Reports and Exports* options for an assay is to perform a manual export *after* image acquisition by clicking the **Export** button at the bottom of the Results screen and selecting additional options.

Export

Location: Browse Single Folder

Images Raw Images Colorized Images

Data Well Level CSV Object Level CSV Object Level ACS

Archive Data Set

Report	File Type	Auto Open	Print
PLX_Surface Marker(s)	<input type="checkbox"/> CSV	<input type="button" value="No"/> <input type="button" value="Yes"/>	<input type="button" value="No"/> <input type="button" value="Yes"/>
	<input type="checkbox"/> Excel	<input type="button" value="No"/> <input type="button" value="Yes"/>	<input type="button" value="No"/> <input type="button" value="Yes"/>
	<input type="checkbox"/> PDF	<input type="button" value="No"/> <input type="button" value="Yes"/>	<input type="button" value="No"/> <input type="button" value="Yes"/>
	<input type="checkbox"/> Word	<input type="button" value="No"/> <input type="button" value="Yes"/>	<input type="button" value="No"/> <input type="button" value="Yes"/>

Performing a Recount

If you find it necessary to fine-tune assay parameters after analyzing scan results, click the **Recount** button located at bottom of the Results screen.

Scan Details

Created On: March 30, 2024

Imaging Mode: Brightfield and Fluorescent

Recount Details

Last Used Assay: PLX.6_3SM+Total_CD3-KB + CD4-PE + CD8-APC + Hoechst View

Assay for Recount: PLX.6_3SM+Total_CD3-KB + CD4-PE + CD8-APC + Hoechst View

Assay Description: CSK-A0026

Tag: Select Tag...

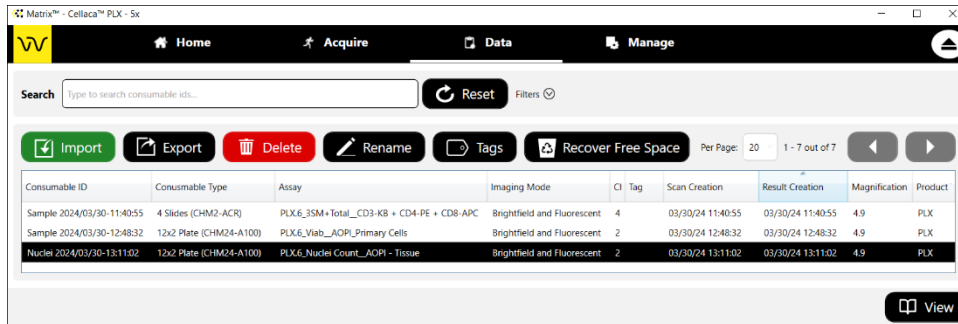
View parameter settings for the *Last Used Assay* or select a new *Assay for Recount* and view/edit its parameters.

You can de-select wells to exclude them from the recount, but you cannot add wells for the recount that were not included in the initial count.

To manage channel mappings in a recount, expand the *Channel Mappings* area. Click on individual mapping indicators and drag them to a new channel. The mapping indicator in that channel will swap positions with the one you are dragging automatically.

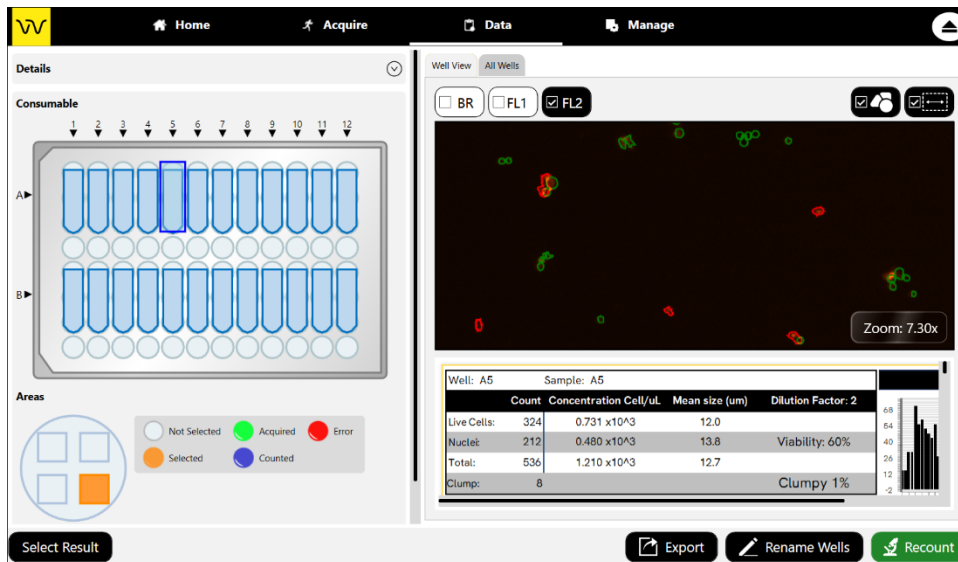
Viewing the Data Tab

The **Data** workflow tab launches the *Select*, *Results* and *Recount* screens which must be completed in sequence when analyzing scans.



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.

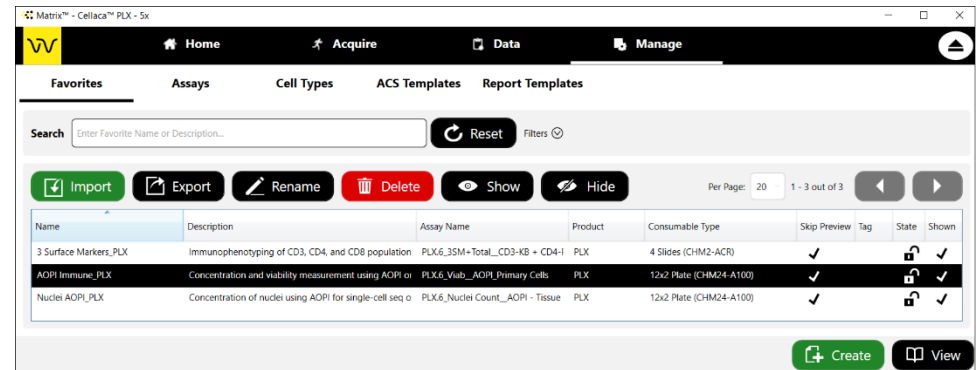
Once a scan result is displayed vary images in the **Well View** tab by changing wells and selecting channels or click the **Summary** tab to view count results.



Viewing the Manage Tab

The **Manage** tab lists *Favorites*, *Assays*, *Cell Types*, *ACS Templates* and *Report Templates* loaded on your instrument system. From this screen you can import/export, rename, delete or show/hide any of these entities in their respective lists.

The *State* column displays icons indicating whether an entity is *Locked* or *Unlocked*. Although a locked assay, cell type or report template cannot be edited, you can select it to use as a source for creating a new entity and click the **Save As** button to save it using a new name.



ACS Templates can be imported on to your system via the Matrix software. You can assign an ACS template to an assay if the *Object Level ACS* export and *Use Template* options are selected. On export, data will auto populate into the specified template and be saved as a file. *If the Auto Open option is also selected and the De Novo Software FCS Express application is installed on your system, FCS Express will be launched for viewing export data.*

Revvity provides extensive assay, cell type and report template libraries as defaults in the Matrix software. Contact Support for current listings of these libraries or for help with creating new assays, customizing cell type parameter settings and defining new cell types. For assistance with report templates, share with us the goal you are trying to accomplish, and we will create a new template to accommodate your needs.

Evaluating Viability Methods

When evaluating viability methods, it is critically important to use a single aliquot from the stock cell culture to perform *all* testing. The cell sample should be evaluated for concentration prior to staining.

Note: Cell concentrations of 1.0×10^5 – 1.0×10^7 cells/mL can be analyzed on the Cellaca PLX, with a concentration of 1.0×10^6 cells/mL being optimal.

If comparing the *Trypan Blue* and *AO/PI* methods, a portion of the sample should be stained with trypan blue and another portion stained with *AO/PI*.

Dilution or concentration of a cell sample may be required based on initial concentration. *It is recommended to use cell culture media for dilution.*

Using Trypan Blue Viability Method

Brightfield imaging and the *Trypan Blue Viability Method* can be used to determine the number, concentration and percentage of live cells for cell lines and cultured primary cells. Brightfield imaging with trypan blue staining is *not* recommended for samples containing debris, platelets or red blood cells. For accurately differentiating nucleated cells, fluorescence is required.

Preparing a Cell Sample for Trypan Blue Viability Determination

Invert the tube containing cells ten times (10x) and pipette up and down 10x to generate a homogeneous cell sample and reduce cell clumps. Do *not* shake or vortex the sample as this may damage cell membranes.

For viability measurement, stain cells by combining 50 μ L of cell sample with 50 μ L of a 0.2% trypan blue staining solution (for a final concentration of 0.1% trypan blue). Gently mix by pipetting up and down 10x.

Using AO/PI Viability Method

Dual-fluorescence methods have been developed to accurately determine nucleated cell concentration and viability in primary cell samples containing debris and non-nucleated cells, including platelets and red blood cells.

In the *AO/PI Viability Method*, acridine orange (AO) enters all cells and stains their DNA causing nucleated cells to fluoresce *Green* (in the *470/534 Channel*), while propidium iodide (PI) only enters dead cells with compromised membranes and stains their DNA causing them to fluoresce *Red* (in the *531/655 Channel*).

- Cells stained with *both* AO and PI fluoresce *Red* due to quenching.
- Live nucleated cells are easily identified in the *Green* FL channel.
- Dead nucleated cells are easily identified in the *Red* FL channel.

As a result, debris and non-nucleated cells do not interfere with nucleated cell counts when using the *AO/PI* viability method.

Preparing a Cell Sample for AO/PI Viability Determination

Invert the tube containing cells ten times (10x) and pipette up and down 10x to generate a homogeneous cell sample and reduce cell clumps. Do *not* shake or vortex the sample as this may damage cell membranes.

For viability measurement, stain cells by combining 50 μ L of cell sample with 50 μ L of *AO/PI* staining solution. For whole blood and other viscous samples, draw sample in and out of the pipette tip at least once prior to transferring for staining. Gently mix stained solution by pipetting up and down 10x before adding sample to counting plate loading wells.

The table below shows recommended dilutions when prepping cell samples for *AO/PI* viability and final *Dilution Factor* to enter for a few sample types.

<i>Sample Type</i>	<i>Preliminary Dilution</i>	<i>Volume of Sample</i>	<i>Volume of AO/PI</i>	<i>Final Dilution Factor</i>
Whole peripheral blood or cord blood	1:10	50 μ L	50 μ L	20
PBMCs following Ficoll separation	<i>Not Required</i>	50 μ L	50 μ L	2
Mononuclear cells from processed bone marrow	<i>Not Required</i>	50 μ L	50 μ L	2
Tumor digest/ Tissue digest	<i>Not Required</i>	50 μ L	50 μ L	2
Stem cells from CD34+ separation	<i>Not Required</i>	50 μ L	50 μ L	2

Viability Staining Solution Guidelines

Use the following staining solution guidelines to prepare cell samples.

<i>Stain Type</i>	<i>Use with Cell Sample</i>	<i>Dilution Factor</i>
Trypan Blue (0.2%)	1:1	2
AO (CS1-0108-5mL)	1:1	2
PI (CS1-0109-5mL)	1:1	2
AO/PI (CS2-0106-5mL)	1:1	2
AO/PI (CS2-0106-25mL)	1:1	2

Available PLX Reagents and Kits

Revvity offers a variety of ready-to-use assay reagents and reagent kits to accurately perform fluorescence-based assays for surface marker detection, apoptosis, cell counting and viability.

Available Cellaca PLX assay reagents and kits include:

- Single Surface Marker
- Single Isotype Control
- Single Surface Marker Viability
- Two Surface Marker Total Cell
- Two Surface Marker Dead Cell
- Two Surface Marker Viability
- Three Surface Marker Total Cell
- Three Surface Marker Dead Cell
- Apoptosis
- Viability for Fluorescent Protein-expressing Cells

Visit the *Reagents and Kits for Cell Counting and Cell-Based Assays* page on our website for a current listing or contact your Revvity Sales representative to purchase Cellaca PLX assay reagents and kits directly from Revvity.

Available Product Documentation

See the following documentation for additional instrument information:

- **8003789 Cellaca PLX User Manual** for instrument operation, care and maintenance details (available as a PDF on desktop of the Revvity-provided Laptop shipped with instrument).
- **8003394 Matrix Software User Manual** for complete details on using software functionality (available as a PDF on desktop of Revvity-provided Laptop shipped with instrument).
- **8002871 Cellaca MX/PLX Plate Loading Template** for use when loading CHM24-A100 (12x2) and CHM24-B100 (8x3) Revvity counting plates.
- **8002845 Cellaca MX/PLX Focus Guide** for assistance with optimal focus.

Contacting Support

If there is a technical issue with your instrument, contact Support by visiting <https://www.revvity.com/contact-us/instrument-support-and-service> or by sending email to: CellC-support@revvity.com

Trained specialists are available to assist your team with sample analysis and optimization of assay/cell type imaging parameters.

When reporting a technical issue, it is recommended that you record any error messages generated, the sequence of steps leading up to the error, and the Serial Number of the instrument *prior* to contacting Support.