



Cellometer K2 Matrix software improvements.

Modernized software to improve your workflow with an option for 21 CFR Part 11

Upgrade your existing Cellometer™ K2 fluorescent cell counter to the Matrix™ software to make viewing, analyzing, record-keeping, and reporting on cells even more accessible. The Matrix software also operates the Cellaca™ MX high-throughput cell counter, and Cellaca PLX high-throughput image cytometer, allowing for less training time when moving between instruments and the ability to harmonize and streamline cell-based assays between groups with low and high-throughput needs.

- Guided workflow with easy selection and editing of assay settings and touchscreen operation
- Improved usability with larger on-screen images, more zooming steps, customizable brightfield, fluorescence, and counted overlays
- Improved customizable data reporting options including histograms, scatter plots, and custom calculations
- Centralized database storage of raw and quantified data, assay settings, and report templates
- Future-proof Windows 11 software overhaul for better accessibility and reliability



- New export functions including Excel and PDF files
- Name editing capacity for data sets after they have been run
- Ability to enable 21 CFR Part 11 ready module

Indicates the number of images captured

Select which channels to view

Guided workflow steps

Auto export results for additional analysis

Quickly view results on screen

The screenshot shows the Cellometer K2 Matrix software interface. On the left is a sidebar with a 'Guided workflow steps' box highlighting the 'Results' button. The main area shows a 'Wells' view with a grid of wells; one well is highlighted with a red box and an arrow pointing to the text 'Indicates the number of images captured'. Above the main view are tabs for 'Well View' and 'All Wells', and a row of buttons: 'BR', 'FL1', 'FL2', 'Zoom', and 'Counted'. An arrow points from the text 'Select which channels to view' to the 'FL1' and 'FL2' buttons. The main view displays a microscopic image of cells. At the bottom, a table shows results: 'Count', 'Concentration (cells/mL) at Dilution Factor: 2', and 'Mean Size (um)'. An arrow points from the text 'Auto export results for additional analysis' to the 'Export' button. Another arrow points from the text 'Quickly view results on screen' to the table.

Count	Concentration (cells/mL) at Dilution Factor: 2	Mean Size (um)
Live: 1028	3.570 x10 ⁶	16.7
Dead: 668	2.310 x10 ⁶	14.9
Total: 1696	5.880 x10 ⁶	16.0

21 CFR Part 11 with Matrix software

We have taken our rigorous GMP/GLP standards and extended them to encompass FDA 21 CFR Part 11 acceptance criteria as well as the recommendations for acceptance, maintenance, submission, and storage of electronic records and electronic signatures.

	Roles with permissions	User login	Audit trails	E-Signature
Cellometer K2 with Matrix and 21 CFR	Customizable	✓	✓	✓
Cellometer K2 with Cellometer Software GMP/GLP	Predefined	✓	✓	x

21 CFR Part 11 ready advantages

- Revvity offers a scalable 21 CFR Part 11 ready option that supports:
- Audit trails, time stamping, user access control, and electronic signatures
 - Assay locking and change tracking

- Creation of new customizable defined roles with multiple configurations to fit your workflow and environment compared to three limited options in GMP/GLP software
- A close-looped database system to aid with data integrity and security compared to previous folder/file structure with the potential to be modified

Understanding role privileges

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

- **Data** – Control, create, manage, export/import, and approve/reject data during the counting and analysis workflow
- **Settings** – Access the **Assays**, **Cell Types**, and **Report Templates** options under the Manage tab to create, manage, and export/import these entities
- **Users, Roles, and Audits** - 21 CFR Part 11 ready module allows users creation and management of the Matrix software **Users**, **Roles**, and **Audits** options under the **Manage** tab to monitor the audit trail automatically generated by the system

Privileges

Data:	☐ Create	☐ Read	☐ Update	☐ Delete	☐ Export	☐ Import	☐ Approval
Settings:	☐ Create	☐ Read	☐ Update	☐ Delete	☐ Export	☐ Import	
Users:	☐ Create	☐ Read	☐ Update	☐ Delete			
Roles:	☐ Create	☐ Read	☐ Update	☐ Delete			
Audits:	☐ Read	☐ Export					

