

Release Notes WAVEcontrol v4.10.0

These release notes list all new features, changes, and fixes available with the WAVEcontrol software version v4.10.0.

The release date is July 15, 2026.

1. New Features

- 1.1. A waveRAPID simulator is now available in WAVEcontrol (*Evaluate & Simulate* -> *Simulate*).
- 1.2. Copy-and-paste functionality is now available in the tables of the Wizard Rack View and Autosampler View. When placing cursor in a cell without selecting any content, paste inserts at the cursor. If selecting some content, paste replaces only the selected text.
- 1.3. The "Apply to All" and "Apply to Selected" functions in Evaluations have been expanded to support additional settings across multiple evaluations.
 - The additional parameters available for multiple evaluation selection are listed below:
 - Settings -> Kinetics
 - Kinetic parameters (table) -> Show
 - Settings -> Equilibrium
 - Kinetic parameters (table) -> Show
 - Use Log axis
 - Equilibrium graph
 - Show Legend
 - Show Kinetic Parameters
 - Display
 - Graph Content -> All
 - Colors
 - In Evaluations, manually defined graph scales can now be applied to multiple selected graphs using the "Use Custom Bounds" option. The feature is available for Kinetics (*Display* -> *Custom Bounds*) and Equilibrium graphs (on graph -> *Use Custom Bounds*).
- 1.4. The graph batch export functionality in *Evaluations* -> *Toolbox* has been expanded to support both adding all graphs to reports and exporting them as image files.
 - The "Export All" functionality under "Toolbox" was modified to include "Add to Report" and "Export to PNG"
 - The "Scope" dropdown selection list was added to allow selection of the graph types to be exported (for multicycle kinetics: Kinetics and/or Equilibrium)
 - PNG files are exported using the following naming convention:
Kinetics/Equilibrium_Cycle_Channel_Analyte

- 1.5. To ensure software performance, the .NET framework is updated to .NET 10.
- 1.6. A "Check Release Notes" link has been added to *Help -> About*, allowing users to review software changes before deciding whether to upgrade.
- 1.7. A selected colour in the tables "Channel color" and "Cycle Type colors" under "Preferences" gets saved immediately after a user's selection.

2. Fixed Issues

- 2.1. Improved software stability during "Export All" operations involving large numbers of evaluations and during long-running kinetic experiments.
- 2.2. Resolved issue that prevented simulations from being cloned.
- 2.3. In Evaluations, the manually positioned kinetic parameter box now retains its position when graphs are added to a report or exported as an image.
- 2.4. Valve errors are now clearly indicated through a red status indicator and message. New cycles are blocked until the error is resolved, and leakage-risk warnings are displayed immediately when an error is detected.
- 2.5. Corrected occasional incorrect pump flowrate when pump refill was necessary during dissociation (in pulsed injections only).
- 2.6. Resolved issue related to background-thread view changes that could cause the application to close unexpectedly.
- 2.7. Optimized the priming process to improve fluidic preparation before measurements.

Great news! How can I get it?

From the 15th of July 2026 a window will prompt once the WAVEcontrol is launched, notifying that a new update is available. Simply select Download and Install.

Please note that internet access and access to <https://github.com> is required for auto-updates to be executed.

The release can also be manually downloaded [here](#) from the Malvern Panalytical website.

To ensure optimal performance of WAVEcontrol please ensure that Windows and Dell driver updates are completed periodically. Automatic updates are recommended to be disabled to prevent automatic restarts but should be manually completed at a convenient time to allow for the correct functioning of the PC.

For any queries, please contact Creoptix support:
support.creoptix@malvernpanalytical.com.

Thank you,

Creoptix Team