

Release Notes WAVEcontrol v4.5.18

These release notes contain a list of all new features, changes, and fixes available with the WAVEcontrol software version v4.5.18.

The release date is October 3rd, 2023.

1. New Features

1.1. General

1.1.1. Additional SFT evaluation criteria “Noise (RMS) 30 ul/min @ 10Hz” has been added.

Note: The feature only works if a new SFT evaluation is added to a series.

1.2. Configuration

1.2.1. On the *Autosampler* view it is now possible to use 384 deep-well plates. Plate type “384 Wells” has been renamed to “384 Wells Low” and a new plate type “384 Wells High” has been introduced with a max volume of 240µL.

1.2.2. On the *Autosampler* view max volume of plate type “96 Wells High” has been increased from 1000µL to 2000µL to accommodate for larger volume plates.

1.2.3. In *Regeneration Free Kinetics* wizard, it is now possible to set multiple blocks of Blanks before every Sample.

1.3. Evaluation

1.3.1. The *Summary Table* now includes a new “QC” column which displays a summary of the *Quality Panel*. The same summary QC indicator is also present on the left upper corner of each sensorgram.

The color coding is as follows: Red (Fail) > Orange (Warning) > Blue (Verify) > Green (Pass).

Please note that thorough analysis of the data is still advised, the QC indicator is only a tool to aid in the analysis but does not replace the user.

1.3.2. The *Summary Table* now includes a “Rating” column which allows to rate single evaluations by clicking on the stars displayed. The default setting displays 3 empty stars (☆☆☆) which can upon selection become filled (★) allowing to rate evaluations with 1, 2 or 3 stars.

1.3.3. An additional column “Comment” has been added to the *Summary Table* which allows you to add a comment to each evaluation. Max length is 20 characters.

Fixed Issues

1.4. General

1.4.1. Performance and stability improvements.

1.5. Configuration

1.5.1. In the *Regeneration Free Kinetics* wizard, the order of blocks of Blanks and Analytes has now been fixed when multiple analytes are used.

1.5.2. The default tubing wash option for Kinetics and RAPID Kinetics series wizards is set to "Loop only" now. Therefore, by default the flow channels over the sensing area are not flushed with the washing agent.

1.6. Evaluation

1.6.1. "Export to Excel" under *Evaluations* has been fixed.

1.6.2. *Summary Table* export has been improved. When copying the *Summary Table* contents into Excel, the table view is now more transparent and organized.

- Evaluations are grouped by BioModel.
- Summary QC indicator detailed in 1.3.1 is exported as Red = 4 > Orange = 3 > Blue = 2 > Green = 1 to enable filtering on Excel file.
- Star rating detailed in 1.3.2 is exported as 0, 1, 2, 3 depending on how many starts are selected (filled) by the user.

Great news! How can I get it?

From October 3rd 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 are auto-updates to be executed.

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 allows for the correcting functioning of the PC.

If you do not get the update notification:

- For customers eligible to software updates: simply select *About* under the *Help* menu on your WAVEcontrol and click on Download and Install.
- If not sure whether you are eligible for software updates or are looking to update your current version: simply contact support: support.creoptix@malvernpanalytical.com

Thank you,

Creoptix Team