

Release Notes WAVEcontrol v4.5.14

These release notes contain a list of all new features, changes, and fixes available with the WAVEcontrol software version v4.5.14.

The release date is November 29th, 2022.

1. New Features

1.1. General

1.1.1. Revised WAVEcontrol and WAVEdelta manuals: revised layout.

1.2. Configuration

1.2.1. The *CFCA wizard* has been improved with the introduction of regeneration step between different flow-rate cycles.

1.2.2. Sample pick up improved to reduce potential air gaps during sample injection.

1.3. Evaluation

1.3.1. *Quality Panel* in the *CFCA wizard* has been adjusted to CFCA data, where absence of curvature is desired as opposed to kinetics data where curvature is essential for high quality results.

2. Fixed Issues

2.1. General

2.1.1. Performance and stability improvements.

2.2. Evaluation

2.2.1. Evaluation model display has been fixed in *CFCA wizard* to avoid user error on model selection during evaluation.

2.2.2. The display of the labels in the *parameter table* when copying into CSV in *CFCA wizard* has been fixed and aligned with label displayed in WAVEcontrol.

2.2.3. Calculation of *QC ratio* in *CFCA wizard* has been improved. Previously good fitting data would often result in a low QC ratio inconsistent with data quality due a bug in the calculations which has now be resolved resulting in increased reliability of parameter calculation.

Great news! How can I get it?

From the 29th of November 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.

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 upgrade your current version: simply contact support: support.creoptix@malvernpanalytical.com

Thank you,

Creoptix Team