

## Release Notes WAVEcontrol v4.5.17

These release notes contain a list of all new features, changes, and fixes available with the WAVEcontrol software version v4.5.17.

The release date is June 9<sup>th</sup>, 2023.

### 1. New Features

This version of WAVEcontrol includes no new features.

### 2. Fixed Issues

#### 2.1. General

2.1.1. Performance and stability improvements.

2.1.2. SFT evaluation criteria updated: Upper bound for DMSO Level at 30 ul/min increased from 160 to 200 pg/mm<sup>2</sup>. SFT template updated with respective changes.

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

2.1.3. SFT evaluation criteria updated: Upper bound for DMSO Level at 200 ul/min increased from 160 to 200 pg/mm<sup>2</sup>. SFT template updated with respective changes.

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

#### 2.2. Configuration

2.2.1. Configuration of sample rack upon *Immobilization* series cloning has been improved and now cloning maintains original rack type.

#### 2.3. Evaluation

2.3.1. *Ligand Screening* evaluation using the *Use single concentration* setting (traditional injection with single concentration) has now been fixed.

Note: The fix only works if a new evaluation is added to a *Ligand Screening* series.

2.3.2. *Ligand Screening* evaluation using the *Use single concentration* setting for traditional injection for multiple concentrations (MCK) has now been fixed.

Note: The fix only works if a new evaluation is added to a *Ligand Screening* series.

2.3.3. Manual  $R_{\max}$  adjustment has been fixed and now allows setting to lower than 1.0.

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

2.3.4. Ligand name now correctly displayed in the evaluation summary table when analysing *Ligand Screening* data and can be exported in csv format when copying from table.

Note: This only works if a new evaluation is added to a *Ligand Screening* series. After the new evaluation is added, the project must be saved otherwise no changes will be visible if navigating to another screen and back.

2.3.5. Interpolation error leading to inability to evaluate data from *Ligand Screening* series has been fixed.

### **Great news! How can I get it?**

From June 9<sup>th</sup> 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](mailto:support.creoptix@malvernpanalytical.com)

Thank you,

Creoptix Team