

Release Notes WAVEcontrol v4.9.3

These release notes list all new features, changes, and fixes available with the WAVEcontrol software version v4.9.3.

The release date is March 4, 2026.

1. Fixed Issues

- 1.1. Instrument communication improvements implemented in release v4.9.1 for WAVEdelta system (4-channel instruments) are also applied to WAVEsystem (2-channel instruments).
- 1.2. Fixed minor delay at the start of dissociation as introduced in v4.9.1.
- 1.3. Improved communication with updated firmware of syringe pump.

Great news! How can I get it?

From the 4th of March 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 correcting functioning of the PC.

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

Thank you,

Creoptix Team