



## Release Notes for the ZE5™ Cell Analyzer and Everest™ Software

### Version 1.2

August 2017

## Contents

Introduction.....	1
Upgrading to Version 1.2.....	2
Documentation .....	3
Additional Information .....	4
Fixed Issues.....	4
Known Issues .....	6
Contacting Technical Support.....	7

## Introduction

The ZE5 Cell Analyzer is a compact benchtop flow cytometer that characterizes cells and their properties by streaming hydrodynamically focused cells through up to five spatially separated laser beams at varying wavelengths.

Everest Software is the comprehensive instrument control application for the ZE5 Cell Analyzer. Everest Software controls all functions of the ZE5 Cell Analyzer and provides accurate data acquisition and user-friendly data analysis. Key features of Everest Software include:

- Automated startup and shutdown
- Quality control process and onboard QC beads
- Fluorochrome selector assisting in experimental design
- Acquisition and file storage for up to 100 million events
- Multichannel triggering
- Hit detection
- Heat map

## Upgrading to Version 1.2

Upgrading Everest Software and your ZE5 Cell Analyzer requires the following general steps:

- Upgrading Everest Software on the Everest computer
- Updating the ZE5 instrument firmware

### Upgrading Everest Software

**Important:** The software upgrade process removes the currently installed version of Everest Software. Before starting the installation process, ensure that no experiments are running on the instrument and that you have saved all data and exited the software.

**Note:** You must be an administrator on the Everest Software computer in order to upgrade the software. See your system administrator for more information.

#### **To upgrade Everest software**

1. Verify that you have exited Everest Software 1.1.
2. From the ZE5 Cell Analyzer product page on the Bio-Rad website, locate the file Everest Setup v1.2.zip and download it to a folder on the Everest computer, for example the Downloads folder.
3. In Windows Explorer, navigate to and open the Downloads folder and double-click Everest Setup v1.2.zip.
4. Extract the file Everest Setup v1.2.exe.
5. Double-click Everest Setup v1.2.exe to start the Everest installer.
6. You are prompted to install the ZE5 firmware and ZE5 service tool updates. Click Install.

**Tip:** Installing the updates can take a few minutes. After the updates are installed you see the Everest Software installer wizard's Welcome screen.

7. On the Welcome screen, click Next.
8. Accept the license agreement and click Next.
9. Follow the remaining instructions in the wizard to install the software.
10. When the installation is complete, click Finish to exit the wizard.

When completed, the Everest icon appears on the desktop.

## **Updating ZE5 Cell Analyzer Firmware**

When you start Everest Software, the ZE5 Cell Analyzer detects the new firmware package and begins the update process.

### **To update the ZE5 Cell Analyzer firmware**

1. Start Everest Software on the Everest computer.

If the firmware update is required, Everest Software displays the ZE5 firmware upgrade dialog box.

2. Follow the instructions to complete the firmware update.

When the update is complete, the instrument shuts down automatically.

3. Start the ZE 5 Cell Analyzer and log in to Everest Software.

## **Documentation**

Information about the ZE5 Cell Analyzer and Everest Software is available from the following sources:

- ZE5 Cell Analyzer and Everest Software User Guide
- ZE5 Cell Analyzer and Everest Software Quick Start Guide
- Everest Software training modules

Everest Software training modules help you become familiar with basic system functionality. These modules run in your default browser, but Internet access is not required.

### **To access the ZE5 Cell Analyzer and Everest Software User Guide**

1. In Everest Software, click the Help and Information menu button in the upper right corner.
2. Select User Manual to open the PDF.

### **To access training modules**

1. In Everest Software, click the Help and Information menu button in the upper right corner.
2. Select Training Module.
3. In the Help and Training dialog box, click the link to the training module you want to view.
4. The training module opens in your default browser.

### **To access the latest product documentation**

- Visit the ZE5 Cell Analyzer product page on the Bio-Rad website ([www.bio-rad.com/ZE5](http://www.bio-rad.com/ZE5)).

## **Additional Information**

### **Monitor and Computer Operating System**

Minimum monitor resolution is 2560 x 1080.

Bio-Rad supports Everest Software only on the Microsoft Windows 10 Pro operating system.

### **System Updates for Security**

Bio-Rad recommends keeping Windows software current by scheduling automatic operating system updates. Instructions can be found here: <https://www.lifewire.com/how-to-change-windows-update-settings-2625778>. Best practices:

- Allow Windows to automatically check for, download, and install updates.
- Schedule operating system updates during hours when the computer will not be needed for use with the ZE5 Cell Analyzer.
- Do not require notification to restart. Allow restarts to occur without user intervention so that the computer will be ready for use after automatic updates.

### **Gate Limit Added to New Samples**

After adding a gate limit to a sample using Assign Gate Limit Regions in the pop-up toolbar of a plot, if you edit the experiment and add more samples, the gate limit is applied to the new samples.

This is as designed. If you do not want to apply the gate limit to any new samples you might add to an experiment, apply the gate limit using the batch tools in the workspace Settings screen instead of using Assign Gate Limit Regions in the pop-up toolbar of a plot.

## **Fixed Issues**

- Shutting the system down with Autostart up enabled functions properly.
- Users can cancel the Add Polygon action correctly.
- Users can now enter the QC bead lot number when replacing bead bottles.
- While editing a multicolor experiment, plots are now added for all parameters for the negative compensation control.
- The system no longer displays a warning when it starts up successfully.
- Time plots function properly on the Analysis screen.
- The Shutdown When Complete feature now works as intended. The instrument shuts down after completion when a user records events for a stat tube add-on.
- The system no longer displays a warning if the sample has not run dry.
- The Filter Configuration screen is editable and displays the filter layout properly.

- The software works as intended when time plots are deleted on the Analysis tab.
- The system selects the proper tanks during startup.
- The Loader door is operational during shutdown after the probe is in the Home position.
- The system displays a warning that the Waste bottle is full during start up and shut down.

## Known Issues

- **Color Gating**

Color gating is not available in Everest Software version 1.2.

- **Analysis of Multipanel Experiments**

All experiments are analyzed in global mode in Everest Software version 1.2.

- **The software crashes when Apply to All is selected in the Analysis tab**

The software can crash when filters are applied to all parameters in the Analysis tab.

- **In a multipanel experiment, data and settings for the second panel do not appear during acquisition**

The software displays the data and settings for the second panel in the Analysis tab after the experiment completes.

- **The software turns on sheath while instrument is in the Pause state**

If the instrument is left in the Pause state for a period of time, the software might start sheath flow.

- **The software does not display the final rows of the QC report**

The scroll bar does not scroll down to the end of the screen, therefore the final rows of the QC report are not visible.

**Workaround:** Move to the Analysis tab and then return to the QC report. The screen refreshes and the final rows are visible.

- **Incorrect Information in the ZE5 Cell Analyzer and Everest Software User Guide**

- Information in the User Guide pertaining to color gating is incorrect. Color gating is not available in Everest Software version 1.2.
- Information in the User Guide pertaining to the analysis of multipanel experiments is incorrect. All experiments are analyzed in global mode in Everest Software version 1.2.
- In the Analyzing, Saving, and Printing chapter of the User Guide, the Publish Toolbar and Preparing Reports topics are incorrect:
  - The Reset toolbar button is not included in Everest Software version 1.2.
  - The Print Size and Print Layout toolbar buttons are not included in Everest Software version 1.2; default settings are used instead. The default page size is 8.5x11 inches and the default layout is Landscape.
  - The Annotation toolbar button is not included in Everest Software version 1.2. Users can add annotations in the Analysis tab.

## **Contacting Technical Support**

The Bio-Rad Technical Support department in the U.S. is open Monday through Friday, 5:00 AM to 5:00 PM, Pacific time.

**Phone:** 1-800-424-6723, option 2

**Email:** Support@bio-rad.com (U.S./Canada Only)

For technical assistance outside the U.S. and Canada, contact your local technical support office or click the Contact Us link at [www.bio-rad.com](http://www.bio-rad.com).

Copyright © 2017, Bio-Rad Laboratories, Inc. All Rights Reserved. Microtiter is a trademark of Thermo Fisher Scientific Inc. Windows is a trademark of Microsoft Corporation.