

Maintaining the ZE5 Cell Analyzer System When Not in Use

Flow cytometry instruments that are run and maintained on a regular basis, with no long periods between uses, have fewer performance issues.

If your instrument will be unattended for up to a month, Bio-Rad recommends enabling Vacation Mode in Everest Software. Using Vacation Mode, you can schedule automatic startups and shutdowns on a defined schedule without an operator present in the lab. Upon startup, the instrument idles for the specified period of time, running sheath liquid as a maintenance process, before shutting down.

If your instrument will be unattended for longer than a month, use the process defined for long-term storage on page 2.

This document supplements the information in the ZE5 Cell Analyzer and Everest Software User Guide, and contains updated information on how to set up and use Vacation Mode, as well as preparing your ZE5 Cell Analyzer for long term storage.

Setting up Vacation Mode

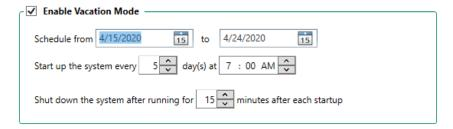
To maintain optimal system performance, Bio-Rad recommends that you run the ZE5 Cell Analyzer regularly, without long periods of rest. If your instrument will not be in use for an extended period (up to 30 days) users with administrative access can enable Vacation Mode to schedule the automatic startup and automatic shutdown processes. During the vacation period, the startup, idling, and shutdown processes occur at the specified times and intervals.

To set up Vacation Mode in Everest Software

- 1. Log into Everest Software as a user with administrative access.
- 2. Click the icon in the upper-right corner to open the main menu, and then click Preferences.
- 3. Select Enable Vacation Mode.
- 4. To set the vacation date range, click each calendar icon to set the start and end dates.

Note: At the end of the vacation period, Vacation Mode is automatically disabled.

5. Specify how often and at what time of day the system startup should occur.



6. Specify how many minutes after startup should the system shut down.

Important: Bio-Rad recommends a frequency of at least once per week, for 15 minutes of idling time. Ensure that bulk fluidics bottles contain enough fluids and space to support the number of runs occurring during the vacation period. If you adhere to the recommended idling length of 15 minutes, with fully filled DI water tanks and empty waste tanks, you can expect approximately 22 repeats of startup, idling, and shutdown.

- 7. Click OK to save the changes and close the Preferences dialog.
- 8. Log out of, but do not shut down, Everest Software.

Before and during the vacation period, a notification appears in the Login window.



Preparing the ZE5 Cell Analyzer and Everest Software for Long-Term Storage

If your ZE5 Cell Analyzer will be unused for a month or longer, Bio-Rad recommends that you follow this procedure to protect the operation of your instrument.

To set up your ZE5 Cell Analyzer for long-term storage

- 1. Follow the decontamination procedure in the Everest Software User Guide, Maintenance chapter.

 The procedure provides instructions for decontaminating both the fluidics line and sample path.
- 2. After system decontamination is finished, close Everest Software and power off the computer.
- 3. Power off the ZE5 instrument using the power button located on the back of the instrument.
- 4. If the ZE5 Cell Analyzer is not connected to an uninterrupted power supply (UPS), remove the plug from the wall outlet (otherwise, leave it plugged in).

Contacting Technical Support

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.

Legal Notices

No part of this publication may be reproduced or transmitted in any form or by any means, electronic or mechanical, including photocopy, recording, or any information storage or retrieval system, without permission in writing from Bio-Rad Laboratories, Inc.

Bio-Rad reserves the right to modify its products and services at any time. This guide is subject to change without notice. Although prepared to ensure accuracy, Bio-Rad assumes no liability for errors or omissions, or for any damage resulting from the application or use of this information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.

All trademarks used herein are the property of their respective owner.

Copyright ©2020 by Bio-Rad Laboratories, Inc. All rights reserved.