



## Release Notes for Bio-Plex Manager™ MP

Version 1.0

September 2013

The following sections are included in the Release Notes:

1. [xPONENT License Requirements](#)
2. [xPONENT 6-Month and 12-Month Maintenance Alerts](#)
3. [Known Issues and Workarounds](#)
4. [Installation Issues](#)
5. [Obtaining Updates and Resources](#)
6. [Technical Support](#)

### 1. xPONENT License Requirements

Bio-Plex Manager MP requires a specific xPONENT license. The xPONENT license must be a permanent license, it cannot be a temporary license, and it must include these modules:

- Base Software
- Automation
- Instrument Control
- Remote Web Monitoring

The following modules must *not* be licensed:

- 21 CFR Part 11
- Security

#### To view your xPONENT license

1. Start the xPONENT software and log on.
2. Click Admin.
3. Click Licensing.

The window displays information about your xPONENT license, including the modules for which you have licenses.

If you need a new license, contact Bio-Rad Technical Support.

## **2. xPONENT 6-Month and 12-Month Maintenance Alerts**

Bio-Plex Manager MP displays alerts when the 6-month and 12-month maintenance of the MAGPIX instrument is due. You need the following kits to perform the maintenance:

- Bio-Plex® MAGPIX™ 6-Month Preventive Maintenance Kit (cat #171-012010)
- Bio-Plex MAGPIX 12-Month Preventive Maintenance Kit (cat #171-012012)

Bio-Plex Manager MP periodically displays alerts until the instrument maintenance has been performed. It may take some time to order and receive the kits and complete the maintenance tasks. You can suppress the alerts.

### **To suppress the alerts**

1. Start xPONENT and log on.
2. Click Admin.
3. Click Schedule.
4. Find the 6 Month and 12 Month tasks in the Subject column.
5. Clear the Enabled checkbox for these scheduled tasks.

## **3. Known Issues and Workarounds**

### **Probe Height Procedure Fails if Heater is On**

When you run the Adjust Probe Height procedure, the heater must be turned off. In most instances, Bio-Plex Manager MP displays a warning if the heater is on, but there are occasional circumstances where you will not see a warning.

If you run the Adjust Probe Height procedure while the heater is on, the procedure may fail and the software will not inform you of the failure.

## **Instrument Status Does Not Update**

When the MAGPIX instrument is turned off, Bio-Plex Manager MP does not update the instrument status in the status bar. The status bar in Bio-Plex Manager MP continues to show that the instrument status is Connected.

If you turn the instrument on while Bio-Plex Manager MP is already running, you must reconnect to the instrument by doing one of the following:

- Click the Instrument menu, then click Reconnect.
- Restart the Bio-Plex Manager MP software.

## **Waste Container Full Error Message Displays Incorrectly**

When the waste container is empty and you continue to see the error message “The waste container is full,” replace the waste container with a different, dry waste container.

When emptying the waste container, replace it with a *different* waste container that is dry. The sensors on the emptied waste container may not be completely dry, causing the instrument to incorrectly detect the waste container as full.

## **CSV File Is Not Created**

The CSV file may not be created after a protocol reading in xPONENT. This can occur in the following scenario:

1. Bio-Plex Manager MP is running during the current session.
2. You exit Bio-Plex Manager MP.
3. You start xPONENT and run a protocol, but the CSV file is not created.

Follow the instructions below to generate the CSV file for the data result.

### **To create the CSV file**

1. Restart your computer.
2. Open the result in xPONENT.
3. Manually export the data to a CSV file.

To avoid this problem, restart your computer after you close Bio-Plex Manager MP and before you start xPONENT.

### **Standard Lots Must Match Assay Panels**

The standard lot definition for an assay panel must contain only those regions that are included in the panel. Therefore, if you modify an assay panel by adding or deleting regions, you must modify the standard lot associated with the panel or create a new standard lot so that only the regions in your panel are included.

### **Routine Log Incorrectly Indicates Prime Routine Status**

The Routine Log incorrectly shows the Prime routine completed successfully. When the Prime routine fails, the software displays a dialog box informing you that the routine did not succeed. However, the log file shows the routine succeeded.

**Hint:** Click the Routine Log link in the Maintenance view to open the Routine Commands window.

### **Selected Wells Do Not Display Correctly**

After a protocol run is interrupted, you can run the protocol again. In the Run Settings dialog, you specify which wells you want read. If you select a subset of the wells, the template in the Create/Run Protocols view usually will visually distinguish the wells that are being read from the wells that are omitted from the reading. However, sometimes the wells all appear to be included.

Note that despite the absence of visual distinction, Bio-Plex Manager MP correctly reads only the selected wells.

### **Copy Command**

In most tables in Bio-Plex Manager MP, you can right-click to open a context menu and click the Copy command. There are instances in which you will get an error message informing you that the Copy command failed. This message is incorrect and you can proceed with pasting your copied rows into Excel or some other software application.

**Note:** This only applies to the Copy command. The Copy Table command works as expected.

## **4. Installation Issues**

### **Installing Bio-Plex Manager MP**

This issue only applies only if you are an existing xPONENT user and you are installing Bio-Plex Manager MP. This issue does *not* apply to new customers of xPONENT and Bio-Plex Manager MP.

Bio-Plex Manager MP software requires Microsoft .NET 4 and a related Microsoft HotFix component (HotFix-KB248432). If these components have not previously been installed on your computer, they are automatically installed during the installation of Bio-Plex Manager MP.

During the installation of these components, you may be prompted to restart your computer. After you restart your computer, the installation of Bio-Plex Manager MP resumes.

If you do not restart your computer when prompted, the installation will not continue. The install window closes and the installation is aborted. If this occurs, restart your computer manually, then rerun the Bio-Plex Manager MP setup file.

If the HotFix installation fails, you will see an error message. If this occurs, close any open installer windows and start the Bio-Plex Manager MP installation again.

### **Uninstalling Bio-Plex Manager MP**

Before you can uninstall Bio-Plex Manager MP, you must first shut down the Bio-Plex Manager MP and xPONENT software applications. Note, shutting down Bio-Plex Manager MP automatically shuts down xPONENT.

If you attempt to uninstall the software without first shutting down Bio-Plex Manager MP, a message displays incorrectly prompting you to shut down xPONENT. You should shut down Bio-Plex Manager MP instead.

### **Other Installation Options**

There are times when you may need to use the installation options in the "Uninstall or change a program" window of the Control Panel. Neither the Change nor the Repair options in the context menu work for Bio-Plex Manager MP. The workaround is described below.

### **To modify or repair your installation**

Run the setup.exe file on the Bio-Plex Manager MP CD, follow the prompts, and choose the Modify option or the Repair option.

## **5. Obtaining Updates and Resources**

From Bio-Plex Manager MP you can view or download updates, training resources, and tutorials that are available online.

**To access updates and resources**

Click the Help menu, then click Go to Resources and Updates.

A browser opens to the Bio-Plex page, where you will find additional resources and updates.

## **6. Technical Support**

Bio-Rad Technical Support in the United States is open Monday–Friday, 5:00 AM to 5:00 PM, Pacific Time. Worldwide technical support is available on the Web at <http://www.consult.bio-rad.com>.

Phone 1-800-424-6723, option 2

Fax 1-510-741-5802

E-mail LSG.TechServ.US@Bio-Rad.com (U.S.)

LSG.TECHServ.Intl@Bio-Rad.com (International)

MAGPIX and xPONENT are trademarks of Luminex Corporation. Excel and Microsoft are trademarks of Microsoft Corporation.

10032256 Rev A