

## Bio-Plex Data Pro™ Software Release Notes Version 1.1

February 24, 2012

These release notes contain the following sections:

1. **What's New in Bio-Plex Data Pro 1.1**
2. Installation Issues
3. Getting Updates and Resources
4. Technical Support

### 1. What's New in Bio-Plex Data Pro 1.1

- Project Sharing — projects can be shared between users. You can export a project to share it with other Bio-Plex Data Pro users, and you can import projects exported by other users.
- Analyte Ordering and Filtering — you can change the order in which the analytes are displayed, and you can remove them from all views until you want to see them again.
- Display of Multiple Analytes in Charts and Tables — multiple analytes can be selected and displayed in all plots, supporting tables, and tabular output.

### 2. Installation Issues

Bio-Plex Data Pro software requires Microsoft .Net 4 and a related Microsoft HotFix component (HotFix-KB248432). These components are installed during installation if they are not already present on your computer. During installation, if you are prompted to restart your computer, do so. After restart, the Bio-Plex Data Pro installation will continue. If an error appears indicating that the HotFix installation failed, close any open installer windows and install Bio-Plex Data Pro again.

If you are prompted to restart your computer and you elect not to do so, the installation window will close. If this occurs, restart your computer manually. After restart, rerun the Bio-Plex Data Pro setup again. If an error appears indicating that the HotFix installation failed, close any open installer windows and install Bio-Plex Data Pro again.

During product removal, the installer may detect a process that must be closed. In rare situations, the installer may fail to close the process. If this happens, close the current installation window and rerun the Bio-Plex Data Pro setup again.

### 3. Getting Updates and Resources

Bio-Plex Data Pro updates, training resources, and tutorials are available on the Bio-Rad Laboratories website. From any view, open the Help Menu and select Go To Resources and Updates to download available resources.

#### 4. Technical Support

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

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

**Fax:** 510-741-5802

**E-mail:**

[LSG.TechServ.US@Bio-Rad.com](mailto:LSG.TechServ.US@Bio-Rad.com) (U.S.)

[LSG.TECHServ.Intl@Bio-Rad.com](mailto:LSG.TECHServ.Intl@Bio-Rad.com) (International)