

Upgrading Bio-Plex Manager™ 4.1, 5.0, or 6.0 Software to Bio-Plex Manager™ 6.1 Software

For technical support, call your local Bio-Rad office, or in the US, call 1-800-424-6723.

BIO-RAD

**Bio-Rad Laboratories, Inc., 2000 Alfred Nobel Drive, Hercules, CA
94547**

Copyright ©2011 Bio-Rad Laboratories, Inc. Reproduction in any form, either print or electronic, is prohibited without written permission of Bio-Rad Laboratories, Inc.

P/N 10022820 Rev B

1 Introduction

These instructions are part of an upgrade kit for computers currently running Bio-Plex Manager™ 4.1, 5.0, or 6.0, that enables them to use Bio-Plex Manager 6.1 software.

A diagram on page 5 shows the hardware connections. The connections are identical to those used for the preceding version, and are included here only for reference.

These are the main steps necessary for the upgrade, described later:

1. Ensure that your system meets the minimum requirements to run Bio-Plex Manager 6.1.
2. Uninstall the older version of Bio-Plex Manager.
3. Install Bio-Plex Manager 6.1 from the software CD.

Bio-Rad recommends uninstalling the previous version. Version 4.1, 5.0, or 6.0 will no longer work after Bio-Plex Manager 6.1 is installed.

If any problems are encountered, contact your local Bio-Rad office, or in the US, call 1-800-4BIORAD (1-800-424-6723).

The following items are provided with each upgrade kit.

Catalog #	Description
171-SUPG41	Bio-Plex Manager 6.1 software to upgrade Bio-Plex Manager 4.1 software; includes an Instrument Control license, MCV plate IV, and a Bio-Plex reservoir
171-SUPG50	Bio-Plex Manager 6.1 software to upgrade Bio-Plex Manager 5.0 software; includes an Instrument Control license
171-SUPG60	Bio-Plex Manager 6.1 software to upgrade Bio-Plex Manager 6.0 software; includes an Instrument Control license.

Use only Bio-Plex components (Validation Kit 4.0, MCV plate IV, Bio-Plex reservoir, and Calibration Kit) with your Bio-Plex system.

For information about other available upgrade kits, refer to the Bio-Plex Manager 6.1 Software Upgrade and Configuration Guide, part #10022817.

New features for Bio-Plex Manager 6.1 include multi-assay protocols, the capability to define standard lots, and an automatic standard curve optimizer.

For information about using Bio-Plex Manager 6.1, refer to the Bio-Plex Manager 6.1 Software User Manual.

2 Software Upgrade

Bio-Plex® LX200 suspension array system components (and LX100 system components manufactured after January 1, 2002) meet minimum requirements to run Bio-Plex Manager™ 6.1 software. For systems manufactured prior to that date, refer to the Bio-Plex Manager 6.1 Software Upgrade and Configuration Guide (part #10022817) for more information.

Uninstall Bio-Plex Manager 4.1, 5.0, or 6.0

1. Ensure that the array reader, microplate platform and HTF are off.
2. Select Start | Control Panel.
3. Select Add or Remove (or Add/Remove) to populate a list of currently installed programs.
4. Select Bio-Plex Manager from the list.

Upgrading Bio-Plex Manager 4.1, 5.0, or 6.0 Software

5. Select Remove and click Yes to the question “Are you sure you want to remove Bio-Plex Manager from your computer?”

If the InstallShield Wizard window is displayed, select the radio button for Remove, and click Next. It should take less than a minute to uninstall.

6. If the Shared File Detected window opens, select the box to not display the message again, and click Yes.
7. Select Finish when you see “Maintenance Complete” and exit to the desktop.
8. Remove the license (HASP) key.

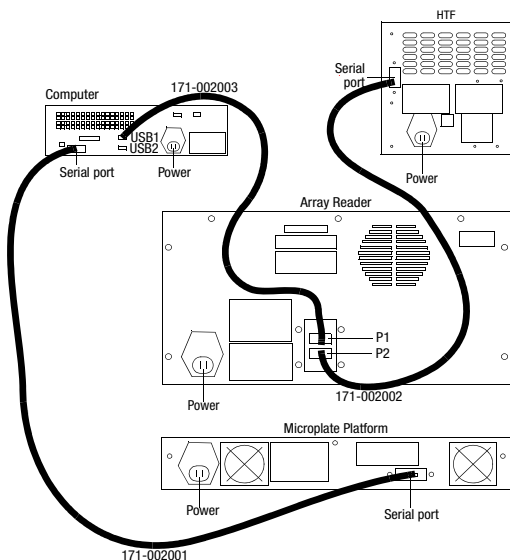
Install Bio-Plex Manager 6.1 Software

1. Insert the Bio-Plex Manager 6.1 software disk into the CD drive. The installation will start automatically.
2. Select Install Bio-Plex Manager. The Wise Installation Wizard will start and the Windows Installer will open.
3. When the Welcome to the Bio-Plex Manager 6.1 Installation Wizard window appears, click Next.
4. Enter name and organization, if desired, and select Next.
5. To accept the default Destination Folder (recommended), select Next.
6. To accept the default Database Destination Folder (recommended), select Next.
7. When queried for the installation type, select Typical, then Next.
8. When you are notified that the installer is Ready to Install the Application, select Next to begin installation.

CAUTION: Under some conditions, a Bio-Plex Manager window may open, noting that the database installation directory already contains a database file.

If you are offered a choice to replace your current database, select “No, do not replace the existing file”. Replacing your database would cause you to lose previously-saved data.

9. When prompted that Bio-Plex Manager 6.1 has been installed, select Finish. The software installation is complete.
10. Insert the new software license key (HASP key) provided with this kit into a USB port in your computer.





BIO-RAD

**Bio-Rad
Laboratories, Inc.**

*Life Science
Group*

Web site www.bio-rad.com **USA** 800 424 6723
Australia 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11
Brazil 55 31 3689 6600 **Canada** 905 364 3435 **China** 86 21 6169 8500
Czech Republic 420 241 430 532 **Denmark** 44 52 10 00
Finland 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 31 884 0
Greece 30 210 777 4396 **Hong Kong** 852 2789 3300
Hungary 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460
Malaysia 60 3 2117 5260 **Mexico** 52 555 488 7670
The Netherlands 0318 540666 **New Zealand** 64 9 415 2280
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700
Russia 7 495 721 14 04 **Singapore** 65 6415 3170
South Africa 27 861 246 723 **Spain** 34 91 590 5200
Sweden 08 555 12700 **Switzerland** 061 717 95 55
Taiwan 886 2 2578 7189 **Thailand** 66 2 6518311
United Kingdom 020 8328 2000