

## Bio-Plex Manager™ Software

### Certification of Software Validation

This is to certify that the product identified at right has been developed, tested, documented, and maintained in compliance with the guidelines and procedures of Bio-Rad's quality management system. Bio-Rad Laboratories, Inc. is certified as compliant with ISO 9001.

The software specification and validation documentation and source code are available for review under a confidential disclosure agreement by government or regulatory agencies. This is conducted through Bio-Rad Laboratories' Life Science Group, located in Hercules, CA USA. This certification has not been validated for clinical diagnostic application.

### Product Identification

#### Bio-Plex Manager Software

Software version	4.1.1
Release date	June 2006
Software version	5.0
Release date	November 2007
Software version	6.0
Release date	April 2010
Software version	6.1
Release date	November 2011
Software version	6.1.1
Release date	March 2013

Joey Lin  
Manager, Software Quality Assurance  
Protein Technologies R&D



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://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 11 5044 5699  
**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 9532 220 **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 **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 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05  
**Taiwan** 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000