

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 **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