

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.

Bio-Rad and Bio-Plex are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

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
Software version	6.2
Release date	August 2018

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