

Certification of Software Validation

This is to certify that the products identified below have been developed, tested, documented, and maintained in compliance with the guidelines and procedures of Bio-Rad's quality system. Bio-Rad Laboratories is certified as compliant with ISO 9001.

Product Identification

Microplate Manager® Software

Software version 5.2.1
Release date February 2004

Ultramark™ OCX Development Platform

Software version 1.0
Release date February 1999

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

Ray Kraus
Software Research and Development Manager
Protein Function Division



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 712 2771 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300/5013478 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 55 5200 05 20 **The Netherlands** 0318 540666 **New Zealand** 64 9415 2280
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 095 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 0861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189/2578 7241 **United Kingdom** 020 8328 2000