

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