

Image Analysis Software

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 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

Quantity One® 1-D Analysis Software

Software version 4.6.1
Release date October 29, 2005

PDQuest™ 2-D Analysis Software

Software version 8.0.0
Release date October 29, 2005

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