

## 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](http://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