

Compliance of The Discovery Series™ Software From Bio-Rad Laboratories With Food and Drug Administration 21 CFR Part 11

In 1997, the United States' Food and Drug Administration (FDA) published Title 21 Code of Federal Regulations (CFR) Part 11 — Electronic Records; Electronic Signatures. The scope of the regulation is to ensure that electronic records and electronic signatures are “trustworthy, reliable, and generally equivalent to paper records and handwritten signatures executed on paper”. This regulation is now becoming mandatory for our customers, scientists in the pharmaceutical and biotechnology industries.

At Bio-Rad Laboratories, we are committed to adapt to our customer needs in a changing environment. Therefore, we are currently working diligently on implementing changes in The Discovery Series™ Quantity One® 1-D analysis software and The Discovery Series PDQuest™ 2-D analysis software to make them compliant with 21 CFR part 11. **We plan to release compliant versions of Quantity One and PDQuest during the year 2003.**

Today, there are already features in The Discovery Series software managing a number of the requirements for a software audit trail. Quantity One software version 4.4 includes:

- User login with name and password
- Ability to define time limit for automatic logout
- Ability to automatically save scan backup files in separate directory
- Complete read-only file history, logging modifications by user with a date and time stamp

For more information, contact:

Bio-Rad Laboratories, Inc.
2000 Alfred Nobel Drive
Hercules, CA 94547, USA

Phone: +1 510-741-1000

Fax: +1 510-741-5811

E-mail: lsg.techserv.us@bio-rad.com



**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 21 507 6191 **Canada** (905) 712-2771 **Czech Republic** + 420 2 41 43 05 32 **China** (86-21) 63052255 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 318 84-177 **Hong Kong** 852-2789-3300 **India** (91-124)-6398112/113/114, 6450092/93 **Israel** 03 951 4127 **Italy** 39 02 216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460 **Latin America** 305-894-5950 **Mexico** 52 5 534 2552 to 54 **The Netherlands** 0318-540666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** + 48 22 8126 672 **Portugal** 351-21-472-7700 **Russia** 7 095 721 1404 **Singapore** 65-62729877 **South Africa** 00 27 11 4428508 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717-9555 **Taiwan** (8862) 2578-7189/2578-7241 **United Kingdom** 020 8328 2000