



AMPLIFICATION

iQTM 5 Optical System Software, Security Edition (Version 2.0)

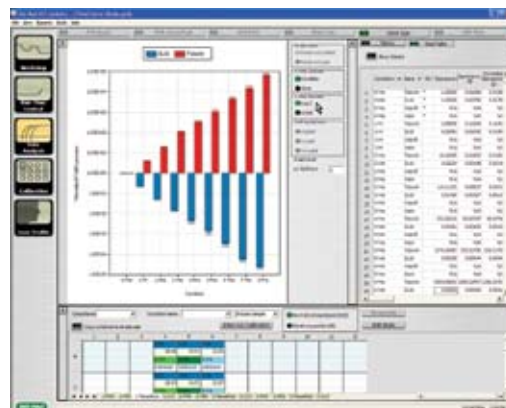
- Tools for compliance with U.S. FDA 21 CFR Part 11
- Password-protected log-in
- Audit trail
- File encryption
- Electronic signatures

Secure Real-Time Data Analysis

Making the most of a powerful real-time PCR system requires a flexible and easy-to-use software package. iQ5 optical system software meets this need for the iQ5 and MyiQTM systems, with quick-setup tools, a full suite of analysis tools, and a variety of presentation options.

Standard Features of iQ5 Optical System Software

- User profile management for creating unique user log-in names with optional password protection
- Built-in analysis modules for absolute quantitation, melt-curve analysis, endpoint analysis, and allelic discrimination (for multiplex data)
- Advanced gene expression analysis options, including comparison to multiple reference genes and correction for reaction efficiency
- Gene Study module for multiframe gene expression analysis with direct comparison of over 5,000 C_T values from iQ5 and MyiQ data files
- Ability to export data directly to a Microsoft Excel spreadsheet



Multifile gene expression analysis. Compare C_T values from different data files with the Gene Study module in iQ5 software. Over 5,000 C_T values can be analyzed in a single Gene Study file.

Additional Features of iQ5 Optical System Software, Security Edition

iQ5 Security Edition software offers all the features of the standard edition, plus important tools for compliance with U.S. FDA 21 CFR Part 11 regulations. iQ5 Security Edition software works with the security features of the Windows XP operating system to provide a secure environment for the maintenance, verification, and tracking of electronic records generated by the software.

BIO-RAD

iQ5 Optical System Software, Security Edition

iQ5 Security Edition software includes these features:

- **Mandatory password-protected log-in** — Valid Windows XP user name and password is required for all users
- **Hardware protection key (HASP HL key)** — Key must be attached to a USB port on the computer to use the software
- **Protected file directories** — Only the logged-in user and the local computer administrator can access the data folder used by the software
- **File encryption** — Files created by iQ5 Security Edition software cannot be opened or edited using other programs
- **Automatic file checking** — Integrity and validity are checked each time a file is opened
- **Electronic signatures*** — Signatures can be applied to any file that can be opened within iQ5 Security Edition software, including protocol, plate, data, Gene Study, and log files, and to calibration information
- **Multiple signature application** — More than one electronic signature can be applied to a file, for example, by both the user who ran the experiment and by another user after a review process
- **Time-and-date stamped audit trails** — Read-only information displayed in the audit trail can be viewed only while the data file of interest is open

* Electronic signatures are intended to be the "legally binding equivalent of traditional handwritten signatures" (U.S. FDA 21 CFR Part 11 Section 11.100 (c)).



Electronic signature window.

Ordering Information

Catalog #	Product Description
170-9753SE01	iQ5 Optical System Software, Security Edition , single-seat license
170-9770	MyiQ Single-Color Real-Time PCR Detection System , includes iCycler® base, optics module, software CD-ROM, 96-well optical reaction module, optical-quality 96-well PCR plates, Microseal® 'B' seals, communication cables, power cords, instructions
170-9780	iQ5 Multicolor Real-Time PCR Detection System , includes iCycler base, optics module, software CD-ROM, 5 installed filter sets, 96-well reaction module, calibration solutions, optical-quality 96-well PCR plates, Microseal 'B' seals, communication cables, power cord, quick reference cards, instructions

Excel and Windows XP are trademarks of Microsoft Corporation.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applera's United States Patent No. 6,814,934 B1 for use in research and for all other fields except the fields of human diagnostics and veterinary diagnostics.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 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 **Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000