

AMPLIFICATION

CFX Manager™ Software, Security Edition

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

Secure Real-Time PCR Analysis

CFX Manager software, Security Edition gives you the power of convenience at your fingertips, making it easy to get started and get great real-time PCR results. As your research needs change and grow, the software can be customized to your applications and preferences, and it gives you the flexibility to expand because it can independently control up to four real-time PCR instruments at once.

Features of CFX Manager System Software

- User preferences give you the power to customize experiment setup and data analysis
- Express loading of protocol and plate files allows you to get started faster
- Automatic data collection allows you to enter or edit plate information before, during, or after an experimental run
- Enhanced data analysis modules perform absolute quantitation, melt-curve analysis, end-point analysis, and allelic discrimination (for multiplex data), with the automatic baseline subtraction and threshold setting
- Export options let you quickly export results and images to Excel spreadsheets and PowerPoint presentations
- Well Groups feature allows analysis of multiple experiments from a single plate
- Advanced analysis tools let you perform normalized gene expression using multiple reference genes and individual reaction efficiencies
- E-mail notification feature alerts you with an attached data file when the run is complete



Multifile gene expression analysis. Compare C_T values from different data files using the Gene Study module in CFX Manager software.



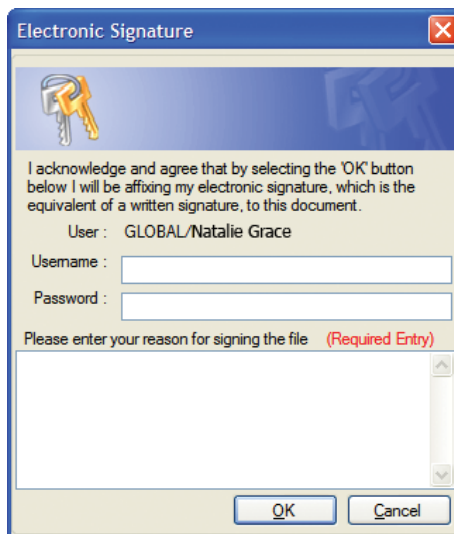
CFX Manager Software, Security Edition

Security Features of CFX Manager Software, Security Edition

CFX Manager software, Security Edition works with the security features of Windows XP and Windows Vista operating systems and offers several major functions and features that are not available in the standard version, including:

- **Embedded tools for compliance with U.S. FDA 21 CFR Part 11 regulations** — tools include electronic signature and added traceability features
- **Mandatory password-protected log-in** — valid Windows XP or Windows Vista username and password are 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 have access to data folders used by the software
- **File encryption** — files created by CFX Manager software, Security Edition cannot be opened or edited using other programs
- **Automatic file checking**— integrity and validity are checked each time a file is opened
- **Multiple electronic signatures* application** — more than one electronic signature can be applied to a software 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 #	Description
184-5000	CFX Manager Software , includes installation CD, instructions
184-5001	CFX Manager Software , Security Edition, includes 1 user license, installation CD, HASP HL key, instructions
184-5005	CFX Manager Software , Security Edition, includes 5 user licenses, 5 installation CDs, 5 HASP HL keys, instructions
184-5010	CFX Manager Software , Security Edition, includes 10 user licenses, 10 installation CDs, 10 HASP HL keys, instructions
185-5096	CFX96™ Real-Time PCR Detection System , includes C1000™ thermal cycler chassis, CFX96 optical reaction module, CFX Manager software, communication cable, power cord, reagent and consumable samples, instructions
184-5096	CFX96 Optical Reaction Module , includes CFX96 optics shuttle, CFX Manager software, communication cable, reagent and consumable samples, instructions
184-1000	C1000 Thermal Cycler Chassis , includes USB flash drive, power cord, instructions; does not include reaction module

Excel, PowerPoint, Windows XP, and Windows Vista 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

**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 905 364 3435 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 6361 7000 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