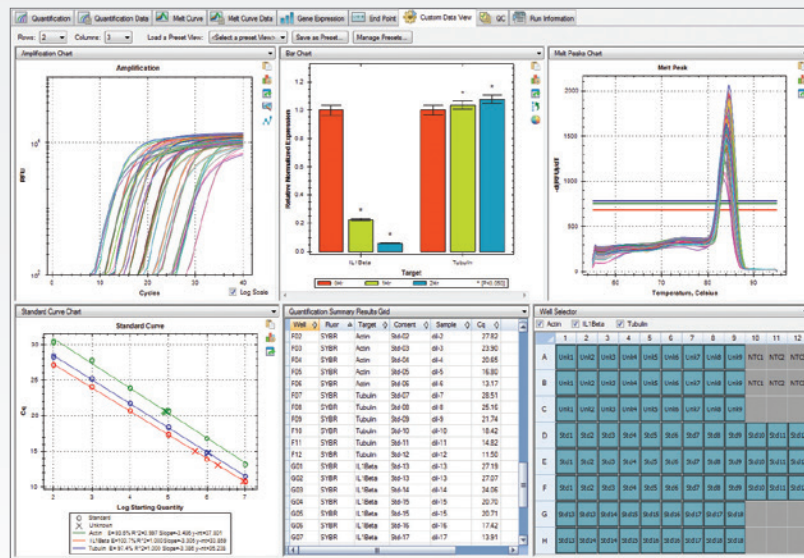


CFX Maestro Software

Real-Time PCR Data Collection, Analysis, and Statistics in One Integrated Package



Custom data view. With custom data view, your most relevant data can be viewed and analyzed in one screen.

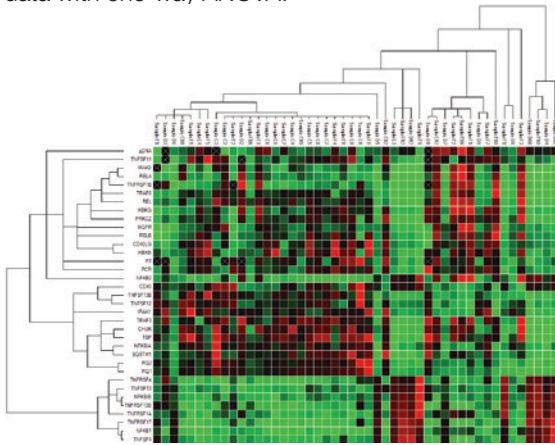
CFX Maestro Software is real-time PCR software that not only collects data, but understands your data and what you want to do with it, combining data collection, statistics, and publication tools into one comprehensive package. Whether you are a novice researcher running your first quantitative PCR (qPCR) experiment or are analyzing thousands of samples across hundreds of plates, CFX Maestro Software is built with your needs in mind and to work in whatever way is best for you.

- Perform statistical analysis of data in seconds with *t*-tests and analysis of variance (ANOVA)
- Work anywhere on a PC
- Set up experiments and replicate easily
- Streamline selection of reference genes
- Create customized and publication-ready graphs



Perform Automatic Statistical Analysis in Seconds

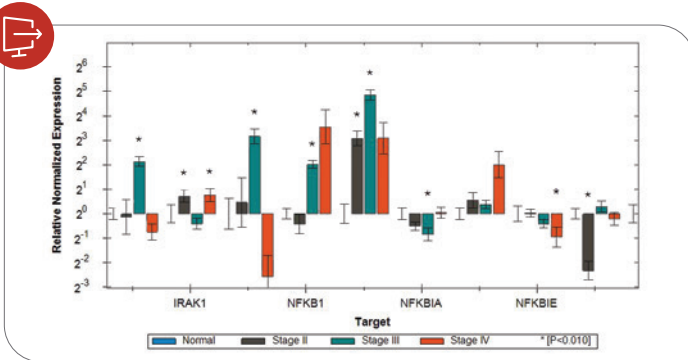
CFX Maestro Software has tools that allow you to view, evaluate, and understand your data. With just a few mouse clicks, you can view normalized expression levels, perform *t*-tests, or analyze your data with one-way ANOVA.



Automatic data analysis. Clustergram showing gene expression levels grouped by hierarchical similarity in a 24-plate study.

Create and Export Publication-Ready Graphics

Visualize your data to suit your needs using bar charts, dot plots, box and whisker plots, and more. Change colors, fonts, titles, and legends, and annotate your graphs with *P* values, text, and arrows to call out specific data. Once your data are displayed in exactly the way you want, export at any size and resolution for presentations, posters, or publication.



Easily visualize your data and export at high resolution.

Work Anywhere on a PC

CFX Maestro Software users can utilize its integrated data analysis, statistics, and publication tools from any PC. Work on your own computer, anywhere, without the need for an Internet connection.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

PrimePCR Assay Integration

Want to simplify your primer design, run design, and setup? With CFX Maestro Software and PrimePCR Real-Time PCR Assays, it's easy. Drag and drop a PrimePCR run file into CFX Maestro Software to automatically populate plate setup and run conditions. After the run, ensure run quality by viewing pass/no pass results from integrated controls (positive control, reverse transcription, genomic DNA, and RNA quality controls).

Additional Features

- Multiple user profiles — allow for easy saving of and access to custom plate setups and qPCR run data files for each user
- Reference Gene Selection Tool — CFX Maestro Software will calculate reference gene stability and help you select the ideal reference genes for your experiments
- Custom data views — view all your most important run and analysis data at a glance

Ordering Information

Catalog # Description

Software*

- 12013758 **CFX Maestro Software 2.3**
 - 12012832 **CFX Maestro Software 2.3, Security Edition**, 1 license
 - 12013028 **CFX Maestro Software 2.3, Security Edition**, 5 licenses
 - 12012834 **CFX Maestro Software 2.3, Chinese Edition**
 - 12012833 **CFX Maestro Software 2.3, Russian Edition**
- * All CFX Maestro Software is for Windows operating systems.

Instruments

- 17005940 **CFX Opus 96 Real-Time PCR System with Starter Pack**, includes integrated 96-well system; Ethernet cable; USB cable; power cord; Hard-Shell™ 96-Well PCR Plates, low profile, thin wall, skirted, white shell/clear well; SsoAdvanced Universal SYBR® Green Supermix, 2 x 1 ml; Microseal™ 'B' Adhesive Seals, 100; CFX Maestro Software
- 17005939 **CFX Opus 384 Real-Time PCR System with Starter Pack**, includes integrated 384-well system; Ethernet cable; USB cable; power cord; Hard-Shell 384-Well PCR Plates, thin wall, skirted, clear shell/white well; SsoAdvanced Universal SYBR® Green Supermix, 2 x 1 ml; Microseal 'B' Adhesive Seals, 100; CFX Maestro Software

Visit bio-rad.com/CFXMaestro for more information.

BIO-RAD, HARD-SHELL, and MICROSEAL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2024 Bio-Rad Laboratories, Inc.

SYBR is a trademark of Thermo Fisher Scientific Inc. Bio-Rad Laboratories, Inc. is licensed by Thermo Fisher Scientific Inc. to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

