

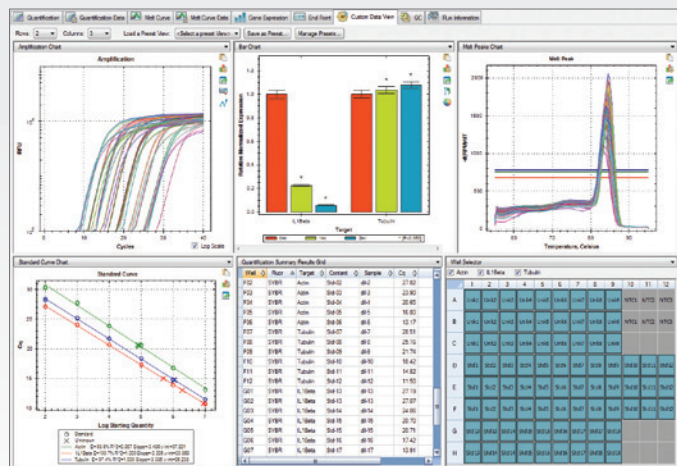


CFX Maestro Software

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

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 or a Mac
- Set up experiments and replicate easily
- Streamline selection of reference genes
- Create customized and publication-ready graphs

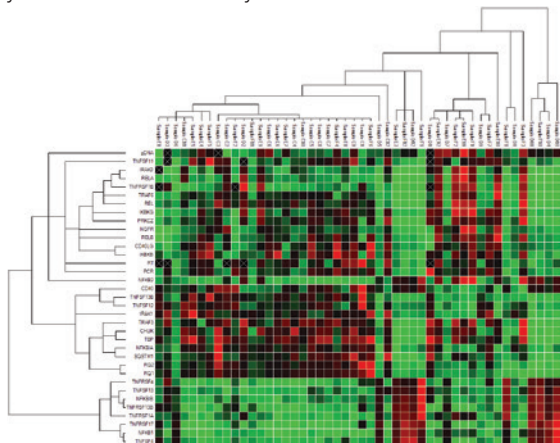


Custom data view. With custom data view, your most relevant data can be viewed and analyzed in one screen.



Perform Automatic Statistical Analysis in Seconds

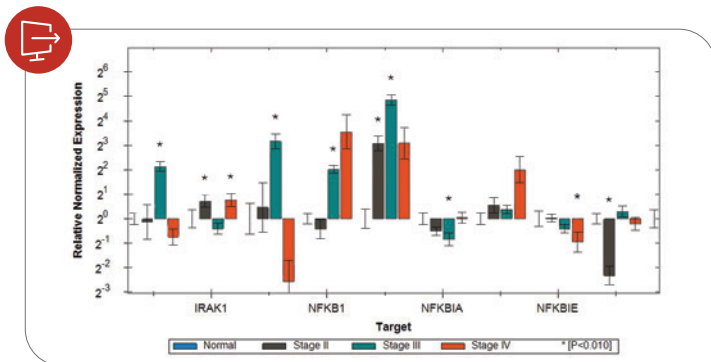
CFX Maestro Software has tools that allow you to view, evaluate, and understand your data. With just a few mouse clicks, you can view normalized expression levels, perform *t*-tests, or analyze your data with one-way ANOVA.



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 or a Mac

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

PrimePCR Assay Integration

Want to simplify your primer design, run design, and setup? With CFX Maestro and PrimePCR Real-Time PCR Assays, it's easy. Drag and drop a PrimePCR run file into CFX Maestro 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 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 2.0
12004128	CFX Maestro Software for Mac
12012832	CFX Maestro Software 2.0, Security Edition, 1 license
12013028	CFX Maestro Software 2.0, Security Edition, 5 licenses
12012834	CFX Maestro Software 2.0, Chinese Edition
12012833	CFX Maestro Software 2.0, Russian Edition

* All software for Windows unless otherwise noted.

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
1854095	CFX96 Touch Deep Well Real-Time PCR System , includes C1000 Touch Thermal Cycler Chassis, CFX96 Deep Well Optical Reaction Module, cables
1855201	CFX Connect Real-Time PCR System , includes CFX Connect Thermal Cycler Chassis, CFX Connect Optical Reaction Module, USB cable

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

BIO-RAD, HARD-SHELL, and MICROSEAL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

Hard-Shell plates are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 7,347,977; 6,340,589; and 6,528,302.

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.

All trademarks used herein are the property of their respective owner.

BIO-RAD

**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 2 3473 4460 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

