



The Next Evolution in qPCR for Your Biopharmaceutical Lab

The reliability you depend on from Bio-Rad with improved connectivity and performance for your modern controlled lab

The CFX Opus Real-Time PCR System is the next evolution in qPCR from Bio-Rad. With improved thermal performance and our proprietary, accurate optical shuttle system, your data will be more consistent than ever. A sleek, modern design includes a refreshed and easy-to-use interface with new connectivity options and features.

In addition to Ethernet compatibility, the CFX Opus System can connect directly to WiFi for completely wireless operation

and data retrieval. Need flexibility? The CFX Opus System is capable of stand-alone operation and direct data transfer to a USB device or it can be connected to a PC.

With internet connectivity, the CFX Opus System can be programmed to send your data and any notifications directly to your email or local networked drive, enabling easy access by collaborators or from other network-connected locations.

Benefits

- Our best thermal uniformity and accuracy yet
- Consistent optics avoid need for revalidation
- Flexible service and qualification plans to fit your needs
- Audit and e-signature compatibility with CFX Maestro Security Edition Software
- Bridging information available



Service on Your Terms

The CFX Opus System comes factory-calibrated and does not require costly regular recalibration. Bio-Rad also offers thermal validation and maintenance plans to fit your needs.

Control Your Data

CFX Maestro Security Edition Software leverages Windows user profiles (local or directory) to offer audit and signature functions to comply with your quality policies.

Improved Performance

In addition to our optical shuttle system, which reads consistent path lengths across the entire plate, our new block provides our best thermal consistency yet, ensuring that you get robust data from all of your wells.

Power in an Easy-to-Use Package

Our user interface and software are intuitive and easy to use for experienced CFX Real-Time PCR System users and novices alike — training and onboarding will be a breeze.

Data Integrity Is Paramount

Avoid losing precious run data. Data for all wells is captured and retained even when a well is omitted or removed from the plate layout. In the event of a power failure, the CFX Opus System will automatically store the active protocol and resume the run when power returns.

Ordering Information

Catalog #	Description
12011319	CFX Opus 96 Real-Time PCR System , includes CFX Opus 96 System and communication cables. Order CFX Maestro Software separately
17005940	CFX Opus 96 Real-Time PCR System with Starter Package , includes CFX Opus 96 System, CFX Maestro Software, license for qbase+ Software, communication cables, reagents, consumables
12011452	CFX Opus 384 Real-Time PCR System , includes CFX Opus 384 System and communication cables. Order CFX Maestro Software separately
17005939	CFX Opus 384 Real-Time PCR System with Starter Package , includes CFX Opus 384 System, CFX Maestro Software, license for qbase+ Software, communication cables, reagents, consumables
12013758	CFX Maestro Software Version 2.0 , compatible with Windows 7 and 10, includes USB installation drive, video quick guides, PDF instruction manual, license for qbase+ Software
12012832	CFX Maestro Software Version 2.0, Security Edition , includes 1 user license, USB installation drive, HASP HL key
12013028	CFX Maestro Software Version 2.0, Security Edition , includes 5 user licenses, 5 USB installation drives, 5 HASP HL keys

Search **CFX Opus** in the Apple App Store and Android Play Store for an augmented reality tour. Contact your local sales representative for an in-person demonstration.

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

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. 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

