

Real-Time PCR

# The Next Evolution in qPCR for Your Research Lab

## The Reliability You Depend on from Bio-Rad with Improved Connectivity and Performance for Your Modern 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

- Optional WiFi connectivity — connect to network drives and the BR.io cloud platform
- Works with the BR.io cloud platform for remote monitoring and analysis
- Our best thermal uniformity and accuracy yet
- Sensitive modern touch screen — responsive even with double gloves
- Updated interface is easier to use and more intuitive than ever
- Uniform optical system free of regular calibration requirements

### Improved Performance

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



### 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 — perfect for training labs and scientists transitioning from other platforms. Our new larger status LED shows you from across the lab whether your instrument is available or in use.

### Analysis How You Want, Where You Want

CFX Maestro Software Desktop Software and our BR.io cloud platform allow multiple users to design and review data anywhere, especially when you set up the CFX Opus System to save your runs to networked file storage or our BR.io cloud platform.

### Ready Out of the Box

CFX Opus Systems are factory calibrated for many common dyes and do not require consistent recalibration. WiFi connectivity allows you to place the instrument anywhere, so you can stay connected without cables.

## Ordering Information

Catalog #	Description
12011319	<b>CFX Opus 96 Real-Time PCR System</b> , includes CFX Opus 96 System and communication cables. Order CFX Maestro Software separately
12011452	<b>CFX Opus 384 Real-Time PCR System</b> , includes CFX Opus 384 System and communication cables. Order CFX Maestro Software separately
12013758	<b>CFX Maestro Software</b> , compatible with Windows 10, includes USB installation drive, video quick guides, PDF instruction manual

Search **CFX qPCR Systems** 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](https://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. © 2023 Bio-Rad Laboratories, Inc.



**Bio-Rad  
Laboratories, Inc.**

---

Life Science  
Group

**Website** [bio-rad.com](https://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

