

Real-Time PCR

# The Next Evolution in qPCR for Your Detection Workflow

## Consistency and Uniformity When Nucleic Acid Detection Workflows Require Large Volumes

The CFX Opus Deepwell Real-Time PCR System is the next evolution in quantitative PCR (qPCR) from Bio-Rad™ Laboratories, Inc. for the detection of DNA and RNA or other large-volume applications. It offers exceptional thermal uniformity and our proprietary, accurate optical shuttle system for reaction volumes up to 125 µl.

With a refreshed, easy-to-use interface and new connectivity options, the CFX Opus Deepwell System is designed to integrate seamlessly and flexibly into your workflows. Connecting via Ethernet or wireless network (WiFi), it can be operated remotely through our BR.io cloud platform or locally by a PC running CFX Maestro Software. With its sensitive touch screen, it can also run in stand-alone mode. The system can send your data to your connected, networked, or remote drives, or save it directly to a USB.

### Benefits

- 5-channel multiplexing plus fluorescence resonance energy transfer (FRET) for protein characterization applications
- One of the largest reaction volumes in the industry
- Optional WiFi connectivity — connect to network drives and the BR.io cloud platform
- Sensitive, modern touch screen — responsive even with double gloves
- Uniform optical system free of regular calibration requirements and passive reference



### Focus on Detection

Some labs are using larger reaction volumes of nonprecious sample matrices to push their nucleic acid detection workflows further. The CFX Opus Deepwell System is specifically designed with exceptionally large volume compatibility along with impressive thermal performance.

### Multiplexing Capability

Our optical shuttle system, which reads consistent pathlengths across all wells, allows for five-color multiplexing without the need for passive reference.

### Flexibility

While the CFX Opus Deepwell System can manage volumes up to 125 µl, most applications will see high performance with as little as 10 µl and the well shape can physically manage volumes down to 1 µl. The system supports the widest range of volumes for many applications.

### Analysis How You Want, Where You Want

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

### Ready Out of the Box

CFX Opus Deepwell 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   |
|-----------|---|
| 12016658  | <b>CFX Opus Deepwell Real-Time PCR System</b> , includes CFX Opus Deepwell 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://www.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://www.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

