



The CFX Opus Real-Time PCR System

Delivering Value Beyond the Initial Investment

When selecting laboratory equipment, the upfront cost is only a fraction of the total investment, with much of the expense driven by indirect factors such as maintenance, downtime, reagent waste, repeat testing, training, and workflow inefficiencies. A total cost of ownership approach looks beyond the initial price to provide a complete view of long-term value. Instruments like the CFX Opus Real-Time PCR System, designed for performance and reliability, help reduce these ongoing costs to support more efficient laboratory operations. The CFX Opus System comes in several configurations, including the CFX Opus DX Real-Time PCR System,* the CFX Opus 96 Real-Time PCR System,† and the CFX Opus 384 Real-Time PCR System.†

No Mandatory Annual Calibration and Minimal Service Needs

When you buy a CFX Opus System, you'll likely meet our service engineer at installation, and rarely after that. Built for long-term reliability with factory calibrated optics and no routine recalibration required, the system minimizes downtime, reduces service interruptions, and maintains predictable operating costs over time.



Reduced Reagent Waste and Repeat Experiments

The CFX Opus System delivers excellent thermal uniformity to produce reproducible results. This reduces failed experiments and repeat testing, conserving reagents and samples. The built-in thermal gradient enables assay optimization in a single run, saving time and resources.



Lower Cost per Data Point

The ability to analyze multiple targets in a single reaction saves reagents and consumables. The CFX Opus System discriminates up to five targets in a single well, generating more data per run while reducing the need for separate experiments.



Simplified Workflow and Data Analysis

CFX Maestro Software simplifies the entire qPCR workflow, reducing training time, manual analysis, and the need for additional tools. Key capabilities that lower cost of ownership:

- Automated setup and statistical analysis to reduce hands-on time
- All-in-one analysis and visualization to eliminate separate software
- Integrated controls and reporting to minimize repeat work and errors



*The CFX Opus Dx System is intended for in vitro diagnostic use by trained laboratory technicians.

†For Research Use Only. Not for use in diagnostic procedures.



Ordering Information

Catalog # Description

Instruments

12011319	CFX Opus 96 Real-Time PCR System, includes CFX Opus 96 System and communication cables. Order CFX Maestro Software separately
12016658	CFX Opus Deepwell Real-Time PCR System, includes CFX Opus Deepwell System and communication cables. Order CFX Maestro Software separately
12011452	CFX Opus 384 Real-Time PCR System, includes CFX Opus 384 System and communication cables. Order CFX Maestro Software separately
12016265	CFX Duet Real-Time PCR System, includes power cord, USB cable. Order CFX Maestro Software separately

Software

All software for Windows unless otherwise noted.

12013758	CFX Maestro Software, includes USB installation drive, video quick guides, PDF instruction manual
12012832	CFX Maestro Software, Security Edition, 1 license
12013028	CFX Maestro Software, Security Edition, 5 licenses
12012833	CFX Maestro Software, Russian Edition, 1 license
12012834	CFX Maestro Software, Chinese Edition, 1 license
1845025	Precision Melt Analysis Software, includes 2 user licenses, installation CD, 2 HASP HL keys, melt calibration kit



PrimePCR PCR Primers, Assays & Arrays:
view experimentally validated PCR primer and probe assays for gene expression analysis

Catalog # Description

Consumables

HSP9655	Hard-Shell™ 96-Well PCR Plates, pkg of 50, low profile, thin wall, skirted, white shell/white wells
HSP9601	Hard-Shell 96-Well PCR Plates, pkg of 50, low profile, thin wall, skirted, white shell/clear wells
MSB1001B	Microseal™ "B" PCR Plate Sealing Film, pkg of 100, optically clear seals for PCR plates
TLS0851	0.2 mL 8-Tube PCR Strips without Caps, low profile, white
TLS0801	0.2 mL 8-Tube PCR Strips without Caps, low profile, clear
TCS0803	0.2 mL Flat PCR Tube 8-Cap Strips, optical, ultraclear

Reagents

1845098	CFX Qualification Plate, 96-well
12011076	Reliance One-Step Multiplex RT-qPCR Supermix, 1 mL (1 × 1 mL), 200 × 20 µL reactions
1725270	SsoAdvanced Universal SYBR® Green Supermix, 2 mL (2 × 1 mL), 200 × 20 µL reactions
1725016	SsoAdvanced Universal Inhibitor-Tolerant SYBR® Green Supermix, 2 mL (2 × 1 mL), 200 × 20 µL reactions
1725095	SingleShot SYBR® Green One-Step Kit for Cell Lysis and RT-qPCR, 100 × 50 µL reactions
1725280	SsoAdvanced Universal Probes Supermix, 2 mL (2 × 1 mL), 200 × 20 µL reactions
1725120	iTaq Universal SYBR® Green Supermix, 2 mL (2 × 1 mL), 200 × 20 µL reactions
1725130	iTaq Universal Probes Supermix, 2 mL (2 × 1 mL), 200 × 20 µL reactions
1708890	iScript cDNA Synthesis Kit, 25 × 20 µL reactions
1708840	iScript Reverse Transcription Supermix, 100 µL, 25 × 20 µL reactions

Visit bio-rad.com/real-time-pcr-systems 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. © 2026 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.

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 55 11 3065 7550 Canada 1 800 361 1808 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 Greece 30 210 7774396 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 000 800 00 24 67 23 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 55 5488 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 7 495 721 14 04 Singapore 65 6415-3170 South Africa 27 21 531 7504 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 662 651 8311 United Arab Emirates 971 4 818 7300 United Kingdom 00 800 00 24 67 23