



# CFX Opus Deepwell Real-Time PCR System



## Large Volumes for the Modern Detection Laboratory

The CFX Opus Deepwell Real-Time PCR System brings modern quantitative PCR (qPCR) to your nucleic acid detection lab. Whether you are pooling samples, running nonprecious matrices, or requiring large volumes for other reasons, the CFX Opus Deepwell System brings your workflow all the features and connectivity you need.

As with our other CFX Opus Systems, you can use CFX Maestro Software (for PC) locally or use your BR.io account to set up a run and access your results anytime and anywhere you have internet access.

### Benefits

- One of the industry's largest qPCR reaction volumes, up to 125  $\mu$ l
- Reaction volumes  $\leq 10$   $\mu$ l depending on the application
- Flexible connectivity and data management options for access and transfer of runs and experiment protocols
  - Operate in stand-alone mode through flexible connectivity options of wireless network (WiFi), Ethernet, and USB
  - Access the BR.io cloud platform for experimental setup and data management at any convenient location with internet access
  - Enjoy continued support for and control of up to 4 CFX instruments with CFX Maestro Software for Windows 7 (64-bit), Windows 10 (64-bit), and Security Edition

- Email your completed run
- Save protocols and completed runs in a network folder
- Set up unique user names, with optional passwords, to easily organize and find your protocols and experiments

### Specifications

#### System

Licensed for real-time PCR	Yes
Sample capacity	96 wells
Sample size	1–125 $\mu$ l (10–125 $\mu$ l recommended)
Dimensions (W x D x H)	33 x 56 x 36 cm (13 x 22 x 14 in.)
Weight	22 kg (48 lb)
Touch-screen user interface	Adjustable, with angle of rotation 12–55°
Communications	USB 2.0 or above, Ethernet, WiFi
Electrical approvals	IEC, CE
Operating system	Windows 10 IoT

#### Optical Detection System

Excitation	6 filtered LEDs
Detection	6 filtered photodiodes
Range of excitation/emission wavelengths	450–730 nm
Scan time	
All channels	12 sec
Single channel	3 sec
Dynamic range	10 orders of magnitude
Sensitivity	Detects 1 copy of target sequence in human genomic DNA
Multiplex analysis	Up to 5 targets per well

continues

## Specifications

### 96-Well Reaction Block

Sample block type	Fixed 96-well sample block
Heating and cooling method	Peltier
Lid heating	30–110°C
Temperature range	4–100°C
Maximum ramp rate	2.5°C/sec
Average ramp rate	2°C/sec
Thermal accuracy	±0.2°C
Thermal uniformity	0.8 ± 0.4°C, measured 10 sec after the block reaches the target temperature

### Gradient

Operational range	30–100°C
Programmable span	1–24°C

### CFX Maestro Data Management and Analysis Software

Operating system	Windows 7 (64-bit), Windows 10 (64-bit)
Memory	Minimum of 1 GB
Data analysis modes	PCR quantification with standard curve Melt curve analysis Gene expression analysis by relative quantity ( $\Delta Cq$ ) or normalized expression ( $\Delta\Delta Cq$ ) Multiple-file gene expression analysis for comparison of an unlimited number of quantification cycle (Cq) values Allelic discrimination Endpoint analysis Data analysis options include bar charts, box-and-whisker plots, dot plots, clustergrams, scatter plots, and volcano plots Statistics include <i>t</i> -test and analysis of variance (ANOVA)
Data export	Save, copy, and print all graphs and spreadsheets from right-click menu Export specified data in multiple formats Copy and paste into a Microsoft Excel, Word, or PowerPoint file Print directly or save as PDFs customizable reports containing run settings, data graphs, and spreadsheets Export to RDML
Image export	Export images at any pixel size and at a resolution up to 600 dpi Save images as .bmp, .jpg, or .png files

Catalog #	Description
12013758	<b>CFX Maestro Software 2</b>
12004128	<b>CFX Maestro Software for Mac</b>
12012832	<b>CFX Maestro Software 2, Security Edition</b> , 1 license
12013028	<b>CFX Maestro Software 2, Security Edition</b> , 5 licenses
12012833	<b>CFX Maestro Software 2, Russian Edition</b>
12012834	<b>CFX Maestro Software 2, Chinese Edition</b>
1845025	<b>Precision Melt Analysis Software</b> , includes 2 user licenses, installation CD, 2 HASP HL keys, melt calibration kit
1845098	<b>CFX Qualification Plate</b> , 96-well
1814000	<b>PX1 PCR Plate Sealer</b> , includes heat sealing instrument
1814030	<b>PCR Plate Heat Seal</b> , pkg of 100, optically clear seals for use with the PX1 PCR Plate Sealer
MSB1001	<b>Microseal 'B' PCR Plate Sealing Film</b> , pkg of 100, optically clear seals for PCR plates
HSP9655	<b>Hard-Shell 96-Well PCR Plates</b> , pkg of 50, low profile, thin wall, skirted, white shell/white wells
HSP9955	<b>Hard-Shell 96-Well PCR Plates</b> , pkg of 50, low profile, thin wall, skirted, white shell/white wells, barcoded
1708840	<b>iScript Reverse Transcription Supermix for RT-qPCR</b> , 25 x 20 $\mu$ l reactions
1725037	<b>iScript Advanced cDNA Synthesis Kit for RT-qPCR</b> , 25 x 20 $\mu$ l reactions
1725270	<b>SsoAdvanced Universal SYBR<sup>®</sup> Green Supermix</b> , 2 ml (2 x 1 ml), 200 x 20 $\mu$ l reactions
1725280	<b>SsoAdvanced Universal Probes Supermix</b> , 2 ml (2 x 1 ml), 200 x 20 $\mu$ l reactions
1725120	<b>iTaq Universal SYBR<sup>®</sup> Green Supermix</b> , 2 ml (2 x 1 ml), 200 x 20 $\mu$ l reactions
1725130	<b>iTaq Universal Probes Supermix</b> , 2 ml (2 x 1 ml), 200 x 20 $\mu$ l reactions
1725200	<b>SsoFast EvaGreen<sup>®</sup> Supermix</b> , 2 ml (2 x 1 ml), 200 x 20 $\mu$ l reactions
1725210	<b>SsoFast EvaGreen<sup>®</sup> Supermix with Low ROX</b> , 2 ml (2 x 1 ml), 200 x 20 $\mu$ l reactions
12010176	<b>Reliance One-Step Multiplex RT-qPCR Supermix</b> , 1 ml (1 x 1 ml), 200 x 20 $\mu$ l reactions
1725150	<b>iTaq Universal SYBR<sup>®</sup> Green One-Step Kit</b> , 100 x 20 $\mu$ l reactions
1725140	<b>iTaq Universal Probes One-Step Kit</b> , 100 x 20 $\mu$ l reactions
1725848	<b>iQ Multiplex Powermix</b> , 50 x 50 $\mu$ l reactions
1725095	<b>SingleShot SYBR<sup>®</sup> Green One-Step Kit for Cell Lysis and RT-qPCR</b> , 100 x 50 $\mu$ l reactions
1725160	<b>SsoAdvanced PreAmp Supermix</b> , 1.25 ml (1 x 1.25 ml), 50 x 50 $\mu$ l reactions
17005726	<b>SEQuoia Complete Stranded RNA Library Prep Kit</b> , 24 reactions
17005710	<b>SEQuoia Complete Stranded RNA Library Prep Kit</b> , 96 reactions
12011928	<b>SEQuoia Dual Indexed Primers Set</b> , 12 vials of unique dual indexed primers, 96 reactions
12011930	<b>SEQuoia Dual Indexed Primers Plate</b> , 96-well plate of unique dual indexed primers, 96 reactions

## Ordering Information

All software for Windows unless otherwise noted.

Catalog #	Description
12016658	<b>CFX Opus Deepwell Real-Time PCR System</b> , includes power cord, USB cable, and Ethernet cable; does not include CFX Maestro Software 2 or WiFi adaptor
17008093	<b>CFX Opus Deepwell Real-Time PCR System with Starter Pack</b> , includes CFX Opus Deepwell System, CFX Maestro Software 2, license for qbase+ Software, power cable, USB cable, Ethernet cable, reagents, consumables

Visit [bio-rad.com/CFXOpus](http://bio-rad.com/CFXOpus) 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.

EvaGreen is a trademark of Biotium, Inc. Bio-Rad Laboratories, Inc. is licensed by Biotium, Inc. to sell reagents containing EvaGreen Dye for use in real-time PCR, for research purposes only.

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  
Laboratories, Inc.**

Life Science  
Group

Website [bio-rad.com](http://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

