



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 μ l
- Reaction volumes ≤ 10 μ 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 experiment 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 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 μ l (10–125 μ 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	±0.4°C (max – min ≤ 0.8°C), measured 10 sec after the block reaches target temperature

Gradient

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

CFX Maestro Data Management and Analysis Software

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

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 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. © 2023 Bio-Rad Laboratories, Inc.

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.

Ordering Information

All software for Windows unless otherwise noted.

Catalog #	Description
12016658	CFX Opus Deepwell Real-Time PCR System , includes power cord, USB cable, and Ethernet cable; does not include CFX Maestro Software or WiFi adaptor
12013758	CFX Maestro Software
12004128	CFX Maestro Software for Mac
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



**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** 82 080 007 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

