

## CFX96 Deep Well Dx System

The CFX96 Deep Well Dx System for in vitro diagnostic use offers industry-leading performance for large-volume reactions. Solid-state optical components provide sensitive detection for up to five targets. Together with unsurpassed thermal cycler performance, easy-to-use software, and the capability to support larger reaction volumes, the CFX96 Deep Well Dx is an open system offering the ultimate flexibility in commercial assay selection or rapid assay development.

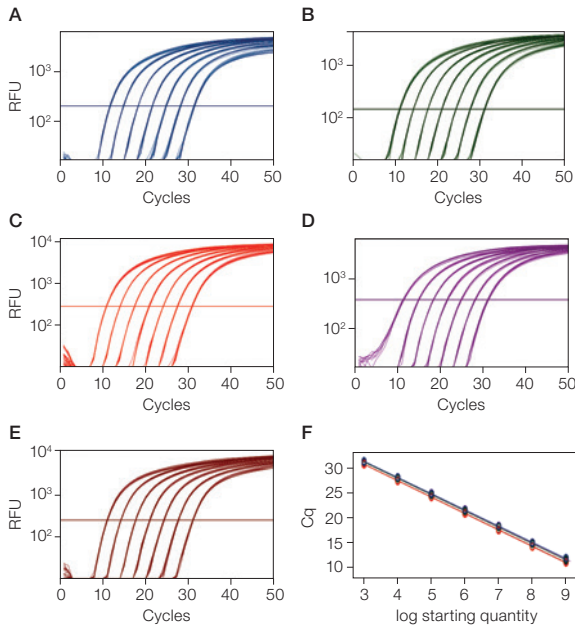
The CFX96 Deep Well Dx System makes it easy for you to:

- Create a personalized system setup with user-specific system access settings and flexible instrument configurations
- Generate robust results right away with factory-calibrated optics and fast system setup
- Streamline data analysis with built-in analysis modules and sophisticated quality control (QC) tools

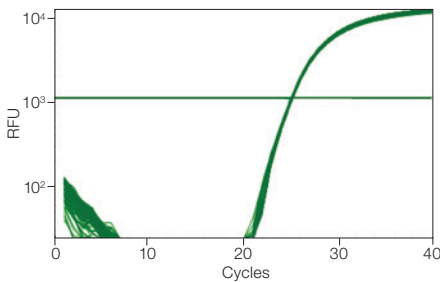


### Specifications

C1000 Dx Thermal Cycler with CFX96 Dx ORM*			
Maximum ramp rate	2.5°C/sec	Gradient	
Average ramp rate	2°C/sec	Operational range	30–100°C
Heating and cooling method	Peltier	Programmable span	1–24°C
Lid	Heats up to 105°C	Temperature range	0–100°C
		Temperature accuracy	±0.2°C of programmed target at 90°C
		Temperature uniformity	±0.4°C well-to-well within 10 sec of arrival at 90°C
Optical Detection			
Excitation	6 filtered LEDs	Dynamic range	10 orders of magnitude
Detection	6 filtered photodiodes	Scan time	
Range of excitation/ emission wavelengths	450–730 nm	All channels	12 sec
Sensitivity	Detects 1 copy of target sequence in human genomic DNA	Single-channel fast scan	3 sec
Software			
Operating systems	Windows 7 (32-bit, 64-bit), Windows 10 (64-bit)	Allelic discrimination	
Memory	Minimum of 1 GB	End-point analysis	
Multiplex analysis	Up to 5 targets per well	Data export	Save, copy, and print all graphs and spreadsheets from right-click menu
Data analysis modes	PCR quantification with standard curve		Export specified data in multiple formats
	Melt curve analysis		Copy and paste into Microsoft Excel, Word, or PowerPoint file
	Gene expression analysis by relative quantity ( $\Delta Cq$ ) or normalized expression ( $\Delta\Delta Cq$ ) with multiple reference genes and individual reaction efficiencies		Customizable reports containing run settings, data graphs, and spreadsheets can be directly printed or saved as PDFs
	Data analysis options include bar chart, clustergram, scatter plot, volcano plot, and heat map		
	Multiple file gene expression analysis for comparison of an unlimited number of quantification cycle (Cq) values		
System			
Sample capacity	96 wells	Dimensions (W x D x H)	33 x 46 x 36 cm (13 x 18 x 14 in.)
Sample size	10–125 $\mu$ l	Weight	21 kg (47 lb)
Communication	USB 2.0	Real-time PCR license	Yes
Electrical approvals	IEC, CE	In vitro diagnostic license	Yes
		CE-IVD mark	Yes



**Linearity of five-target multiplex detection.** A–E, fluorescence data from a series of tenfold dilutions of plasmid DNA ( $10^9$ – $10^3$  copies) amplified using reporter dyes to monitor five targets in a 75  $\mu$ l reaction volume: (■), FAM/cyclophilin; (■), HEX/GAPDH; (■), Texas Red/actin; (■), Cy5/tubulin; (■), Quasar 705/IL-1 $\beta$ ; F, standard curves generated from data in A–E, reaction efficiencies range from 100 to 102%. Cq, quantification cycle; RFU, relative fluorescence units.

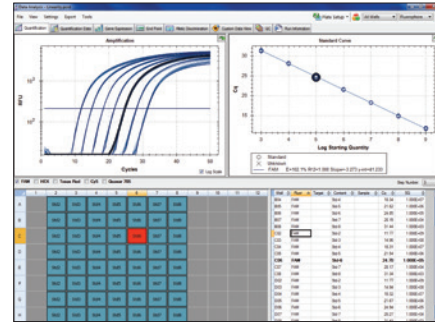


**Excellent uniformity.** IL-1 plasmid template was diluted to  $10^5$  copies/reaction and amplified in the presence of a FAM labeled detection probe with iQ Supermix. Graph shows 96 replicates of 100  $\mu$ l reactions. Average Cq =  $25.14 \pm 0.10$ . RFU, relative fluorescence units.

## Software Solutions for Accurate Results

CFX Manager Dx Software offers tools to simplify real-time PCR for every laboratory. Immediately generate results using the Startup Wizard and intuitive experiment setup. Enter or edit well information before, during, or after a run.

Analyze data when and where you want by receiving an email with an attached data file when a run is completed. When data are in hand, use the comprehensive data analysis, QC, and report tools to take the guesswork out of analyzing and reporting results for any real-time PCR application.



CFX Manager Dx Software data analysis module.

## Ordering Information

To order the CFX96 Deep Well Dx System, you must include both catalog numbers.

Catalog #	Description
1844095-IVD	<b>CFX96 Deep Well Dx ORM</b> , includes CFX Manager Dx Software, version 3.1 (catalog number 12007917)
1841000-IVD	<b>C1000 Dx Thermal Cycler</b>

USA: For in vitro diagnostic (IVD) use. The CFX96 Dx System and CFX96 Deep Well Dx System are registered with the U.S. FDA as Class II 510(k) exempt devices.

Canada: The CFX96 Dx System and CFX96 Deep Well Dx System are registered as Class I devices.

China: The CFX96 Deep Well Dx System has been certified as a Class III medical device by China's National Medical Products Administration (NMPA).

EU: For in vitro diagnostic use. The CFX96 Dx and CFX96 Deep Well Dx Systems meet the requirement of the European In Vitro Diagnostic Regulation (2017/746).

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner. © 2022 Bio-Rad Laboratories, Inc.

Notice regarding high resolution melt analysis:

No rights are granted by Bio-Rad for the use of high resolution melt analysis in the fields of human or veterinary in vitro diagnostics. In addition, it is the responsibility of the purchaser to obtain any intellectual property rights which may be required for its specific applications.

**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 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

