



Advancing qPCR Together

The CFX96 Touch Deep Well Real-Time PCR System offers precise quantification and target discrimination for up to five targets in large-volume reactions. The system incorporates industry-leading technology and stand-alone functionality to provide a robust and easy-to-use real-time PCR system for reliable results.

The CFX96 Touch Deep Well Real-Time PCR System provides:

- **Minimal setup** with factory-calibrated dyes
- **Easy viewing of real-time data traces** with integrated LCD touch screen
- **Fast experiment optimization** with built-in thermal gradient
- **User-friendly data analysis modules** with CFX Maestro Software
- **Effortless export of RDML files** for import into qbase+ Software

Visit bio-rad.com/CFX96DeepWell for more information.



Simplicity through Innovation

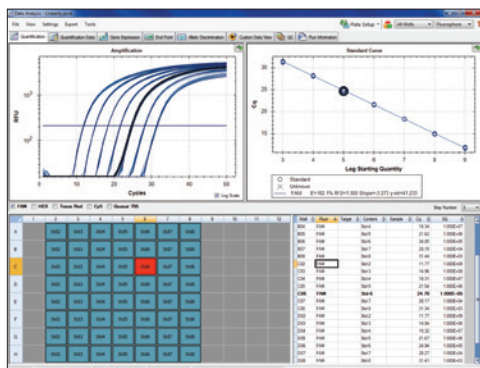
The CFX96 Touch Deep Well System is built on the Bio-Rad C1000 Touch Thermal Cycler. The system incorporates innovative optical technologies with long-lasting LEDs and solid-state components to provide maximal reliability and flexibility.

- Scanning optics read each well individually with high sensitivity and no cross talk to deliver optimal quantitative results
- Multiple data acquisition modes, including a fast scan option for SYBR® Green I or EvaGreen® users, let a run be tailored to suit the application
- Touch screen enables rapid run setup and viewing of data traces in real time
- Several control configurations are available for experiments — run a stand-alone system without a computer or up to 4 instruments from 1 computer

Software Solutions

CFX Maestro Software offers tools that simplify real-time PCR for every researcher. Start experiments in no time following the Startup Wizard and the intuitive experiment setup. Enter or edit well information before, during, or after a run. You can 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 tools to take the guesswork out of analyzing results for any real-time PCR application.

- CFX Maestro Software lets you perform gene expression analysis using multiple reference genes and individual reaction efficiencies
- Precision Melt Analysis Software lets you genotype samples based on their thermal denaturation properties
- Data can be imported into qbase+ Software for independent analysis of gene expression results with proven solutions for quality control, normalization, and inter-run calibration



CFX Maestro Software data analysis module.

Ordering Information

Catalog #	Description
Instruments	
1854096	CFX96 Touch Deep Well Real-Time PCR System with Starter Package , includes C1000 Touch Thermal Cycler Chassis, CFX96 Deep Well Optical Reaction Module, CFX Maestro Software, qbase+ Software license, cables, reagents, consumables
1844096	CFX96 Deep Well Optical Reaction Module for Real-Time PCR Systems with Starter Package , includes CFX96 Deep Well Optical Reaction Module, CFX Maestro Software, qbase+ Software license, cables, reagents, consumables
1814000	PX1 PCR Plate Sealer , includes heat sealing instrument
1845075	CFX Automation System II , includes plate handler and barcode scanner, mounting plate, automation software
Software	
	All software for Windows unless otherwise noted.
12013758	CFX Maestro Software 2.0
12004128	CFX Maestro Software for Mac
12012832	CFX Maestro Software 2.0, Security Edition , 1 license
12013028	CFX Maestro Software 2.0, Security Edition , 5 licenses
12012834	CFX Maestro Software 2.0, Chinese Edition
12012833	CFX Maestro Software 2.0, Russian Edition
1845025	Precision Melt Analysis Software , 2 user licenses, data analysis software, 2 HASP HL keys, calibration kit
Reagents	
1708841	iScript Reverse Transcription Supermix , 400 µl (4 x 100 µl vials), 100 x 20 µl reactions
1725035	iScript gDNA Clear cDNA Synthesis Kit , 100 x 20 µl reactions
1725038	iScript Advanced cDNA Synthesis Kit , 100 x 20 µl reactions
1708891	iScript cDNA Synthesis Kit , 100 x 20 µl reactions
1725271	SsoAdvanced Universal SYBR® Green Supermix , 5 ml (5 x 1 ml vials), 500 x 20 µl reactions
1725281	SsoAdvanced Universal Probes Supermix , 5 ml (5 x 1 ml vials), 500 x 20 µl reactions
1725160	SsoAdvanced PreAmp Supermix , 1.25 ml (1 x 1.25 ml vial), 50 x 50 µl reactions
12010220	Reliance One-Step Multiplex Supermix , 5 ml (5 x 1 ml vials), 1,000 x 20 µl reactions
12010221	Reliance One-Step Multiplex Supermix , 10 ml (2 x 5 ml vials), 2,000 x 20 µl reactions
1725120	iTaq Universal SYBR® Green Supermix , 2 ml (2 x 1 ml vials), 200 x 20 µl reactions
1725151	iTaq Universal SYBR® Green One-Step Kit , 500 x 20 µl reactions
1725130	iTaq Universal Probes Supermix , 2 ml (2 x 1 ml vials), 200 x 20 µl reactions
1725141	iTaq Universal Probes One-Step Kit , 500 x 20 µl reactions
Consumables	
HSP9601	Hard-Shell Low-Profile 96-Well Skirted PCR Plates , white shell/clear well, 50
MSB1001	Microseal 'B' Adhesive Seals , optically clear, 100
MSF1001	Microseal 'F' PCR Plate Seals , foil, pierceable, 100

BIO-RAD, HARD-SHELL, and MICROSEAL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. EvaGreen is a trademark of Biotium, 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. Hard-Shell Plates are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 7,347,977; 6,340,589; and 6,528,302. All trademarks used herein are the property of their respective owner.



Bio-Rad PCR reagents and analytical instruments are manufactured under an ISO 13485:2016 certified Quality Management System and are quality control tested to ensure consistent product performance you can trust.



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

