

## Advancing qPCR Together

The CFX384 Touch Real-Time PCR System brings flexibility and ease of use to researchers performing high-throughput real-time PCR in a 384-well format. With its ability to run without a computer, unsurpassed performance, and powerful yet easy-to-use software, the CFX384 Touch System has been designed to advance your qPCR.

The CFX384 Touch Real-Time PCR System makes it easy for you to:

- Rely on performance with an optical system that uses long-lasting solid-state technology with filtered LEDs and filtered photodiodes for precise quantification and target discrimination
- Conserve your samples and reagents by performing 4-target multiplex reactions with optimal quantitative results using sample volumes as low as 3  $\mu$ l
- Quickly set up runs and monitor amplification traces in real time on the integrated LCD touch screen
- Configure the system to fit your laboratory needs by running the CFX384 Touch System in several control configurations, including integrated with the CFX Automation System II
- Trust your results by integrating the CFX384 Touch Real-Time PCR System with CFX Maestro Software, Security Edition to be compliant with U.S. FDA 21 CFR Part 11 regulations

Visit [bio-rad.com/qPCRSystems](http://bio-rad.com/qPCRSystems) for more information.



### Redefining Innovation

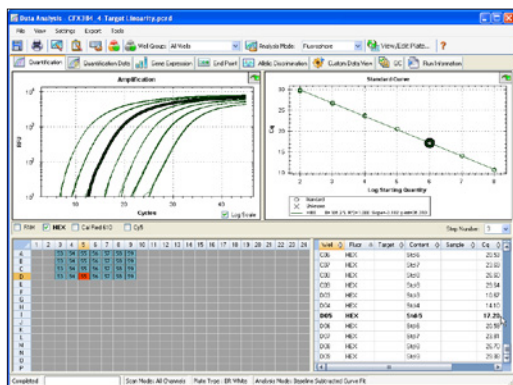
The CFX384 Touch System incorporates innovative optical technologies with the C1000 Touch Thermal Cycler to deliver the most sensitive, reliable detection for real-time PCR applications, including absolute quantification, genetic variation analysis, and gene expression.

- System reads each well individually with high sensitivity and no cross talk to deliver optimal quantitative results
- Compact design fits any laboratory setting and integrated touch screen allows viewing of data traces in real time
- Several control configurations are available for experiments — run a stand-alone system without a computer, up to 4 instruments from 1 computer, or 1 instrument integrated with the CFX Automation System II

### Software Solutions

CFX Maestro Software provides numerous features and tools to streamline data management, from experiment setup to analysis. Enter or edit well information on your own time — before, during, or after a run. Integrate with your laboratory information management system (LIMS) using built-in LIMS file management. When data are in hand, use the advanced data modules to take the guesswork out of analyzing results.

- 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
<b>Instruments</b>	
1855484	<b>CFX384 Touch Real-Time PCR System with Starter Package</b> , includes C1000 Touch Thermal Cycler Chassis, CFX384 Optical Reaction Module, CFX Maestro Software, qbase+ Software license, cables, reagents, consumables
1845384	<b>CFX384 Optical Reaction Module for Real-Time PCR Systems with Starter Package</b> , 384-well optical module for real-time PCR, includes CFX Maestro Software, qbase+ Software license, cables, reagents, consumables, for use with C1000 Touch Thermal Cycler Chassis
1814000	<b>PX1 PCR Plate Sealer</b> , heat sealing instrument
1845075	<b>CFX Automation System II</b> , includes plate handler and barcode scanner, mounting plate, automation software
<b>Software</b>	
	All software for Windows unless otherwise noted.
12013758	<b>CFX Maestro Software 2.0</b>
12004128	<b>CFX Maestro Software for Mac</b>
12012832	<b>CFX Maestro Software 2.0, Security Edition</b> , 1 license
12013028	<b>CFX Maestro Software 2.0, Security Edition</b> , 5 licenses
12012834	<b>CFX Maestro Software 2.0, Chinese Edition</b>
12012833	<b>CFX Maestro Software 2.0, Russian Edition</b>
1845025	<b>Precision Melt Analysis Software</b> , includes 2 user licenses, data analysis software for all CFX Real-Time PCR Systems, 2 HASP HL keys, calibration kit
<b>Reagents</b>	
1708841	<b>iScript Reverse Transcription Supermix</b> , 400 µl (4 x 100 µl vials), 100 x 20 µl reactions
1725035	<b>iScript gDNA Clear cDNA Synthesis Kit</b> , 100 x 20 µl reactions
1725038	<b>iScript Advanced cDNA Synthesis Kit</b> , 100 x 20 µl reactions
1708891	<b>iScript cDNA Synthesis Kit</b> , 100 x 20 µl reactions
1725271	<b>SsoAdvanced Universal SYBR® Green Supermix</b> , 5 ml (5 x 1 ml vials), 500 x 20 µl reactions
1725281	<b>SsoAdvanced Universal Probes Supermix</b> , 5 ml (5 x 1 ml vials), 500 x 20 µl reactions
1725160	<b>SsoAdvanced PreAmp Supermix</b> , 1.25 ml (1 x 1.25 ml vial), 50 x 50 µl reactions
12010220	<b>Reliance One-Step Multiplex RT-qPCR Supermix</b> , 5 ml (5 x 1 ml vials), 1,000 x 20 µl reactions
12010221	<b>Reliance One-Step Multiplex RT-qPCR Supermix</b> , 10 ml (2 x 5 ml vials), 2,000 x 20 µl reactions
1725120	<b>iTaq Universal SYBR® Green Supermix</b> , 2 ml (2 x 1 ml vials), 200 x 20 µl reactions
1725151	<b>iTaq Universal SYBR® Green One-Step Kit</b> , 500 x 20 µl reactions
1725130	<b>iTaq Universal Probes Supermix</b> , 2 ml (2 x 1 ml vials), 200 x 20 µl reactions
1725141	<b>iTaq Universal Probes One-Step Kit</b> , 500 x 20 µl reactions
<b>Consumables</b>	
HSP3805	<b>Hard-Shell 384-Well Skirted PCR Plates</b> , clear shell/white well for higher fluorescent signals, 50
MSB1001	<b>Microseal 'B' Adhesive Seals</b> , optically clear, 100
MSF1001	<b>Microseal 'F' PCR Plate Seals</b> , foil, pierceable, 100

BIO-RAD, HARD-SHELL, and MICROSEAL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. 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

