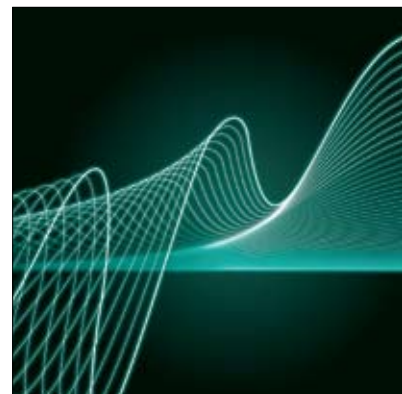


Real-Time PCR: CFX384™ Real-Time PCR Detection System



Designed for the Way You Work

The CFX384 system brings flexibility and ease of use to researchers performing high-throughput real-time PCR in a 384-well format. With the ability to run without a computer, unsurpassed performance, and powerful, yet easy-to-use software, the CFX384 system has been designed for the way you work.

The CFX384 real-time PCR detection system makes it easy for you to:

- Rely on performance — the optical system uses long-lasting solid-state technology with filtered LEDs and filtered photodiodes for precise quantitation and target discrimination
- Conserve your samples and reagents — perform 4-target multiplex reactions with optimal quantitative results, using sample volumes as low as 3 μ l
- Configure the system to fit your laboratory needs — run the CFX384 in several control configurations, including integrated with the CFX automation system
- Streamline your data analysis — use the built-in analysis module and integrated software solutions for high resolution melt analysis or independent gene expression
- Trust your results — integrate the CFX384 real-time PCR detection system with CFX Manager™ software, Security Edition to be compliant with U.S. FDA 21 CFR Part 11 regulations

For more information, visit us on the Web at www.bio-rad.com/qpcrsystems/.

BIO-RAD

Redefining Innovation

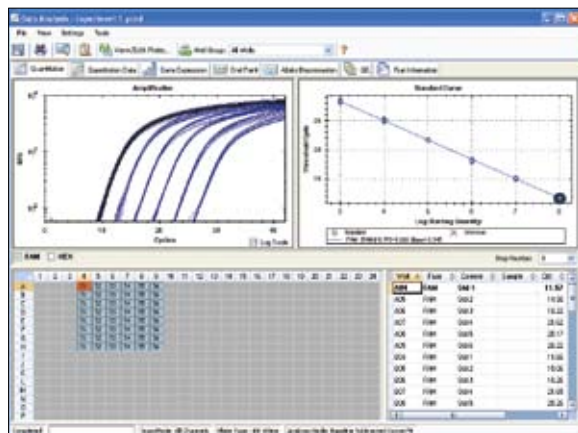
The CFX384 system incorporates innovative optical technologies with the C1000™ thermal cycler, to deliver the most sensitive, reliable detection for real-time PCR applications, including absolute quantitation, genetic variation analysis, and gene expression.

- System reads each well individually with high sensitivity and no cross talk to deliver optimal quantitative results
- Compact design fits any laboratory setting
- Several control configurations are available for experiments — run a stand-alone system with no computer, up to 4 instruments from 1 computer, or 1 instrument integrated with the CFX automation system

Software Solutions

CFX Manager software, which runs on a PC, 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. When data are in hand, use the advanced data modules to take the guesswork out of analyzing results.

- CFX Manager 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
- Import data into qbase^{PLUS} software for independent analysis of gene expression results with proven solutions for quality control, normalization, and inter-run calibration



CFX Manager software data analysis module.

Ordering Information

Catalog #	Description
184-5384	CFX384 Optical Reaction Module , includes CFX Manager software, license for qbase ^{PLUS} software, communication cable, reagent and consumable samples, instructions. Order to upgrade an existing C1000 thermal cycler.
185-5384	CFX384 Real-Time PCR Detection System , includes C1000 thermal cycler chassis, CFX384 optical reaction module, CFX Manager software, license for qbase ^{PLUS} software, communication cable, power cord, reagent and consumable samples, instructions
184-5001	CFX Manager Software, Security Edition , includes 1 user license, installation CD, HASP HL key, instructions
184-5025	Precision Melt Analysis Software , includes 2 user licenses, installation CD, 2 HASP HL keys, calibration kit, instructions
184-5008	CFX Manager Software, Chinese Edition , includes 3 user licenses, installation CD, 3 HASP HL keys, instructions
184-5028	CFX Manager Software, Russian Edition , includes 3 user licenses, installation CD, 3 HASP HL keys, instructions
184-5072	CFX Automation System , includes robotic plate handler, base tray, bar code scanner, CFX automation control software CD, instruction manual
170-8862	iQ™ Supermix , 500 x 50 µl reactions
170-8882	iQ™ SYBR® Green Supermix , 500 x 50 µl reactions
170-8891	iScript™ cDNA Synthesis Kit , 100 x 20 µl reactions, includes 5x iScript reaction mix, iScript reverse transcriptase, nuclease-free water
172-5200	SsoFast™ EvaGreen® Supermix , 200 x 20 µl reactions, 2x mix contains dNTPs, Sso7d fusion polymerase, MgCl ₂ , EvaGreen dye, stabilizers
HSP-3805	Hard-Shell® Thin-Wall 384-Well Skirted PCR Plates , clear shell, white well, 50
MSB-1001	Microseal® 'B' Adhesive Seals , 100

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 Invitrogen Corporation. Bio-Rad Laboratories, Inc. is licensed by Invitrogen Corporation to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in human in vitro diagnostics and all other applied fields under one or more of U.S. Patents 5,656,493, 5,333,675, 5,475,610 (claims 1, 44, 158, 160–163 and 167 only), and 6,703,236 (claims 1–7 only), or corresponding claims in their non-U.S. counterparts, owned by Applied Biosystems. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied's United States Patent 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

Bio-Rad's real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Nos. 6,767,512 and 7,074,367.

Practice of the patented 5' Nuclease Process requires a license from Applied Biosystems. The purchase of these products includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research when used with the separate purchase of Licensed Probe. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

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 Nos. 7,347,977, 6,340,589, and 6,528,302.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400 Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000