

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



Advancing qPCR Together

The CFX384 Touch real-time PCR detection 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 Touch system has been designed to advance your qPCR.

The CFX384 Touch real-time PCR detection 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 μ 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
- Trust your results by integrating the CFX384 Touch 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

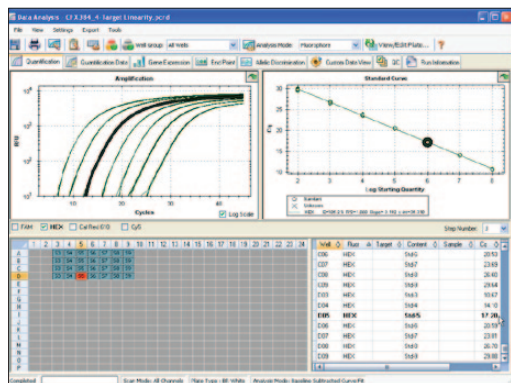
The CFX384 Touch system incorporates innovative optical technologies with the C1000 Touch™ thermal cycler to deliver the most sensitive, reliable detection for real-time PCR applications, including absolute quantification, genetic variation analysis, and gene expression.

- 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

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. 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 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
- Data can be imported 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 , for use with C1000 Touch thermal cycler chassis, includes CFX Manager software, license for qbase ^{PLUS} software, communication cable, reagents, consumables |
| 185-5484 | CFX384 Touch Real-Time PCR Detection System , includes C1000 Touch thermal cycler chassis, CFX384 optical reaction module, CFX Manager software, license for qbase ^{PLUS} software, communication cable, reagents, consumables |
| 184-5001 | CFX Manager Software, Security Edition , includes 1 user license, installation CD, HASP HL key |
| 184-5025 | Precision Melt Analysis Software , includes 2 user licenses, installation CD, 2 HASP HL keys, melt calibration kit |
| 184-5008 | CFX Manager Software, Chinese Edition , includes 3 user licenses, installation CD, 3 HASP HL keys |
| 184-5028 | CFX Manager Software, Russian Edition , includes 3 user licenses, installation CD, 3 HASP HL keys |
| 184-5072 | CFX Automation System , includes robotic plate handler, base tray, bar code scanner, CFX automation control software CD |
| 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 |
| 172-5200 | SsoFast™ EvaGreen® Supermix , 200 x 20 µl reactions, 2x mix contains dNTPs, Sso7d fusion polymerase, MgCl ₂ , EvaGreen dye, stabilizers |
| 172-5230 | SsoFast Probes Supermix , 200 x 20 µl reactions, 2x mix contains dNTPs, Sso7d fusion polymerase, MgCl ₂ , stabilizers |
| HSP-3805 | Hard-Shell® Thin-Wall 384-Well Skirted PCR Plates , clear shell, white well, 50 |
| MSB-1001 | Microseal® 'B' Adhesive Seals , optically clear, 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. HASP is a trademark of Aladdin Knowledge Systems, Ltd. SYBR is a trademark of Molecular Probes, Inc. Bio-Rad Laboratories, Inc. is licensed by Molecular Probes, Inc. to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

Notice regarding Bio-Rad thermal cyclers and real-time systems:

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 U.S. Patent Number 5,475,610 (Claims 1, 44, 158, 160-163, and 167 only), or corresponding claims in its non-U.S. counterpart, owned by Applera Corporation. 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 Applera's United States Patent Number 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 Numbers 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 Numbers 7,347,977; 6,340,589; and 6,528,302.



BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00
France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Malaysia 60 3 2117 5260 Mexico 52 555 488 7670
The Netherlands 0318 540666 New Zealand 64 9 415 2280 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 3170 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 Thailand 66 2 6518311 United Kingdom 020 8328 2000