

## Real-Time PCR: CFX Connect™ Real-Time PCR Detection System



### Advancing qPCR Together

The CFX Connect real-time PCR detection system provides excellent thermal and optical performance with powerful, easy-to-use software for all your SYBR® Green/EvaGreen and duplex experiments.

The CFX Connect real-time PCR detection system offers:

- Easy startup — obtain great results right away with factory-calibrated dyes, quick setup, and intuitive software
- Effortless optimization — save time and reduce costs by optimizing assays in a single run using the thermal gradient
- Powerful data analysis tools — quickly and accurately validate and analyze data with the advanced analysis modules of CFX Manager™ software

For more information, visit us on the Web at [www.bio-rad.com/CFXConnect](http://www.bio-rad.com/CFXConnect).



## Simplicity through Innovation™

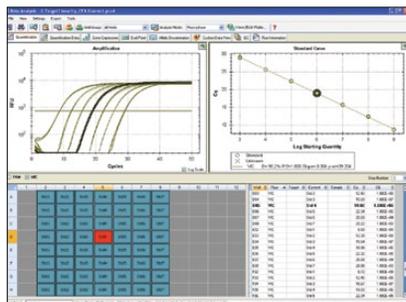
The CFX Connect real-time PCR detection system offers two-target analysis with the same reliable performance as the CFX96 Touch™ system. The system incorporates innovative optical technologies with long-lasting LEDs and solid-state components to provide maximal reliability and flexibility.

- 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 EvaGreen or SYBR® Green I users, let a run be tailored to suit the application
- Independently run up to 4 instruments from 1 computer to increase throughput and flexibility

## Software Solutions

CFX Manager software offers tools that simplify real-time PCR for every researcher. Quickly start experiments by following the intuitive experiment setup. Enter or edit well information before, during, or after a run. Perform data analysis when and where you want by receiving an email with an attached data file after a run is completed. The comprehensive data analysis tools take the guesswork out of analyzing results for any real-time PCR application.

- CFX Manager software lets you perform gene expression analysis using multiple reference genes and individual reaction efficiencies
- Precision Melt Analysis™ software accurately identifies sample genotypes based on their thermal denaturation properties
- Data can be imported into qbase<sup>PLUS</sup> software for access to wizard-guided statistical analysis, advanced copy number variation analysis, and global mean normalization methods



CFX Manager software data analysis module.

## Ordering Information

Catalog #	Description
185-5200	<b>CFX Connect Real-Time PCR Detection System</b> , includes CFX Connect thermal cycler chassis, CFX Connect optical reaction module, CFX Manager software, license for qbase <sup>PLUS</sup> software, communication cable, reagents, consumables
185-5201	<b>CFX Connect Real-Time PCR Detection System</b> , includes CFX Connect thermal cycler chassis, CFX Connect optical reaction module, CFX Manager software, license for qbase <sup>PLUS</sup> software, communication cable
184-5008	<b>CFX Manager Software, Chinese Edition</b> , includes 3 user licenses, installation CD, 3 HASP HL keys
184-5028	<b>CFX Manager Software, Russian Edition</b> , includes 3 user licenses, installation CD, 3 HASP HL keys
184-5025	<b>Precision Melt Analysis Software</b> , includes 2 user licenses, installation CD, 2 HASP HL keys, melt calibration kit
172-5200	<b>SsoFast™ EvaGreen® Supermix</b> , 200 x 20 µl reactions, 2x mix contains dNTPs, Sso7d fusion polymerase, MgCl <sub>2</sub> , EvaGreen dye, stabilizers
172-5230	<b>SsoFast Probes Supermix</b> , 200 x 20 µl reactions, 2x mix contains dNTPs, Sso7d fusion polymerase, MgCl <sub>2</sub> , stabilizers
172-5996	<b>SsoFast EvaGreen Control Assay</b> , predesigned kit for verifying performance of real-time PCR systems or training new users, includes template DNA, primers, supermix, water
HSP-9601	<b>Hard-Shell® Thin-Wall 96-Well Skirted PCR Plates</b> , white shell, clear well, 50
MSB-1001	<b>Microseal® 'B' Adhesive Seals</b> , 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.

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 U.S. 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](http://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 11 5044 5699  
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 9532 220 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 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 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 Thailand 800 88 22 88 United Kingdom 020 8328 2000