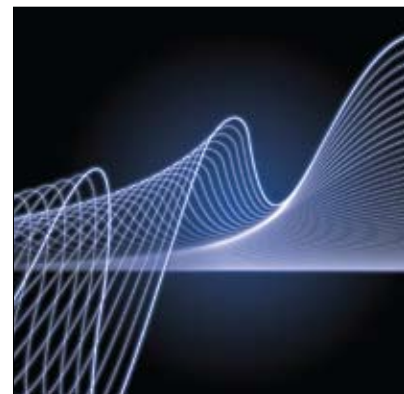


## Real-Time PCR: CFX96™ Real-Time PCR Detection System



## Designed for the Way You Work

Bio-Rad's CFX96 real-time PCR detection system meets all your real-time PCR needs — whether you are running your first experiment or analyzing complex gene expression studies. With five-target detection, unsurpassed thermal cycler performance, and easy-to-use software, the CFX96 system has been designed for the way you work.

The CFX96 real-time PCR system makes it easy for you to:

- Start doing experiments right away, with factory-calibrated optics and minimal setup
- Integrate real-time PCR into your lab with a compact footprint, quiet operation, and the ability to run a stand-alone system without a computer
- Fit experiments into your schedule by shortening run times with fast thermal cycling
- Save research time by optimizing experiments in a single run using a thermal gradient
- Conserve your samples and reagents with true 5-target multiplexing and optimal quantitative results using sample volumes as low as 10  $\mu$ l
- Streamline data analysis with built-in analysis modules and integrated software solutions for high resolution melt analysis or independent gene expression

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

**BIO-RAD**

## Simplicity Through Innovation™

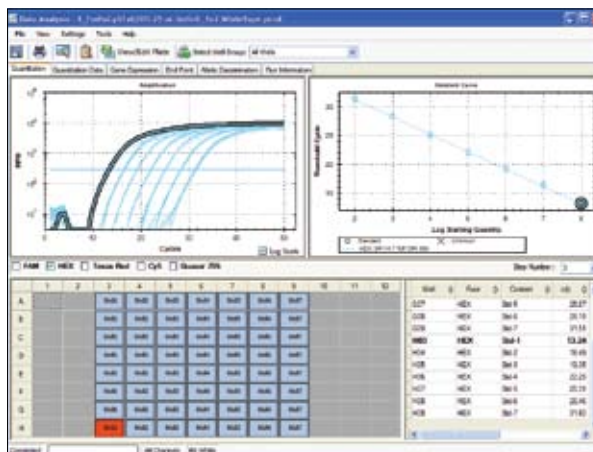
The CFX96 real-time PCR detection system is built on Bio-Rad's fast, friendly, and flexible 1000-series thermal cycling platform. The system incorporates innovative optical technologies with long-lasting LEDs and solid-state components to provide maximal reliability and flexibility.

- 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
- Several control configurations are available — run a stand-alone system or independently run up to 4 instruments from 1 computer

## Software Solutions

CFX Manager™ 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 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<sup>PLUS</sup> 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
184-1000	<b>C1000™ Thermal Cycler Chassis</b> , includes USB flash drive, power cord, instructions; does not include reaction module
184-5096	<b>CFX96 Optical Reaction Module</b> , includes CFX Manager software, license for qbase <sup>PLUS</sup> software, communication cable, reagent and consumable samples, instructions
185-5096	<b>CFX96 Real-Time PCR Detection System</b> , includes C1000 thermal cycler chassis, CFX96 optical reaction module, CFX Manager software, license for qbase <sup>PLUS</sup> software, communication cable, power cord, reagent and consumable samples, instructions
184-5001	<b>CFX Manager Software, Security Edition</b> , includes 1 user license, installation CD, HASP HL key, instructions
184-5025	<b>Precision Melt Analysis Software</b> , includes 2 user licenses, installation CD, 2 HASP HL keys, calibration kit, instructions
184-5008	<b>CFX Manager Software, Chinese Edition</b> , includes 3 user licenses, installation CD, 3 HASP HL keys, instructions
184-5028	<b>CFX Manager Software, Russian Edition</b> , includes 3 user licenses, installation CD, 3 HASP HL keys, instructions
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

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.

Purchase of SsoFast EvaGreen supermix 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. No other patent rights are conveyed expressly, by implication, or by estoppel. 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.

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.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://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