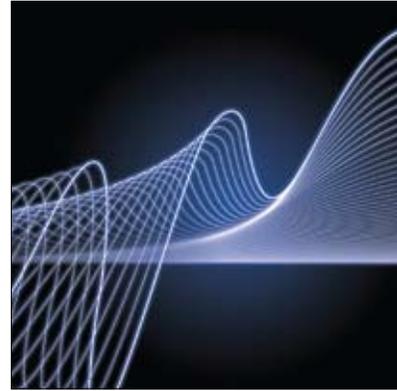


Real-Time PCR:
MiniOpticon™ Real-Time PCR Detection System



Small Size, **Big Performance**

The MiniOpticon system is the smallest real-time PCR detection system available, providing the individual researcher access to sophisticated real-time PCR techniques. When controlled by CFX Manager™ software, this compact, yet powerful package makes it easy for you to:

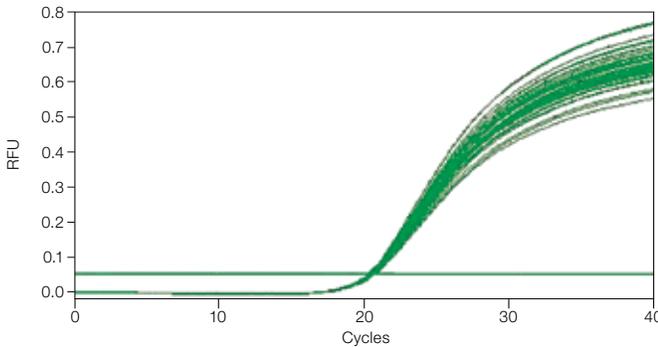
- Save research time by optimizing experiments using the thermal gradient
- Amplify up to 2 targets per well using a range of fluorophores
- Review data with built-in analysis modules, including normalized gene expression analysis using multiple reference genes and target-specific reaction efficiencies
- Analyze data where and when you want by receiving email notification with an attached data file when a run is complete

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

BIO-RAD

Precise Thermal Control

The MiniOpticon offers excellent thermal performance for fast, reproducible runs. A high degree of uniformity, made possible by combining an innovative optical system with the high performance of the MJ Mini™ thermal cycler, ensures reliable and reproducible results from sample to sample and experiment to experiment.



Uniform results. Analysis of 48 replicate 25 µL PCR samples with a $C_T \pm 0.13$ SD. RFU, relative fluorescence units.

The gradient feature allows you to simultaneously incubate samples at up to eight temperatures for any incubation step in your PCR protocol.

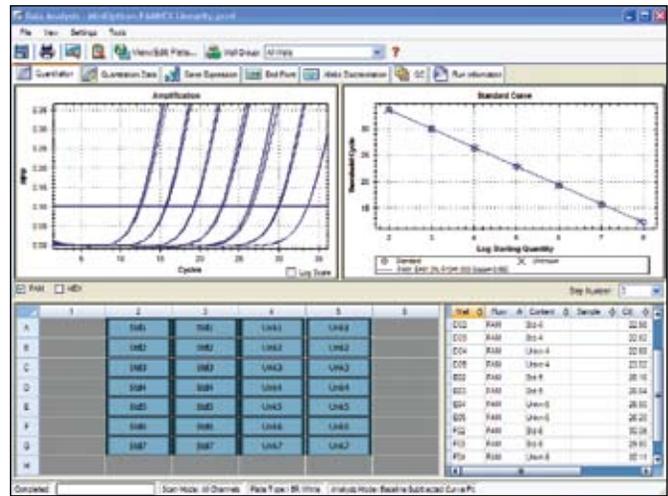


- Gradient operational range: 35–99°C
- Gradient programmable span: 1–16°C
- Gradient accuracy: $\pm 0.4^\circ\text{C}$ of programmed target at end rows within 10 sec (NIST-traceable)

CFX Manager Software

CFX Manager software version 1.5 collects, analyzes, and exports data from the MiniOpticon system. With CFX Manager software you can:

- Get to results faster using the startup wizard and intuitive experiment setup
- Perform normalized gene expression using multiple reference genes and target-specific reaction efficiencies, melt-curve analysis, end-point analysis, and allelic discrimination
- Use the Well Groups feature to analyze multiple experiments from a single plate, each with its own analysis settings
- Customize the settings to suit your needs with user preferences, including unique log-in names, gene and sample name libraries, and data analysis options
- Use the QC tools to validate the quality of your data



CFX Manager software data analysis module.

Ordering Information

Catalog #	Description
CFB-3120	MiniOpticon Real-Time PCR System , includes optical housing, MJ Mini thermal cycler, analysis software
MLL-4851	Multiplate™ Low-Profile 48-Well Unskirted PCR Plates , white, 50
TCS-0803	Optical Flat 8-Cap Strips , for 0.2 ml tubes and plates, ultraclear, 120
TLS-0851	Low-Profile 0.2 ml 8-Tube Strips Without Caps , white, 120 strips (960 tubes)
184-5003	CFX Manager Software, MiniOpticon , includes CFX Manager software version 1.5, installation CD, quick guides, instruction manual for MiniOpticon

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 Biosystems' 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 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