

MyiQ™ Single-Color Real-Time PCR Detection System

The MyiQ real-time PCR detection system offers an affordable solution for the detection of common fluorescent green dyes such as FAM and SYBR Green I. This system interfaces directly with the iCycler® thermal cycler, offering superior features such as thermal gradient capabilities and Peltier-effect driven performance. The MyiQ real-time PCR detection system is a perfect solution for those just getting started with real-time PCR technology as well as those looking for additional instruments to handle increased routine assay demands. This system was developed by the same experts at Bio-Rad who have pioneered performance in real-time PCR. The MyiQ system delivers high-quality data in a single-channel instrument.



Specifications

Operational

Heating rate	Up to 3.3°C/sec
Cooling rate	Up to 2.0°C/sec
Method of heating/cooling	Peltier and Joule
Modes of temperature monitoring	Algorithm
Temperature	
Range	4–100°C
Accuracy	±0.3°C
Uniformity	±0.4°C
Overshoot maximum	<0.5°C
Heated lid	Up to 110°C

Gradient Mode

Gradient accuracy	±0.4°C
Row uniformity	±0.4°C
Gradient range	40–100°C
Temperature differential range	1–25°C

Descriptive

Sample capacity	96 wells
Sample size	15–100 µl
Licensed for PCR	Yes
Communications	USB 2.0 high-speed
Electrical approvals	IEC, CE
Dimensions (W x D x H)	29.2 x 58.4 x 38.7 cm (11.5 x 23 x 15.2")
Weight	17.6 kg (39 lb)

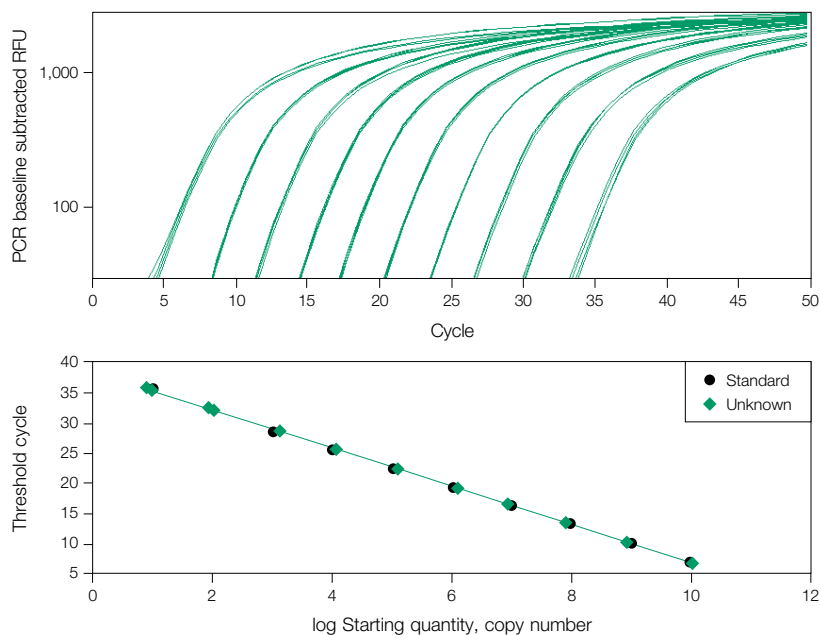
Optical System

Optical lamp	Tungsten-halogen
Fluorescence excitation range	475–495 nm
Fluorescence detection range	515–545 nm
Sensitivity	Detects one copy of IL-1β in human genomic DNA
Dynamic range	9 orders of magnitude
Excitation/emission filters	FAM/SYBR Green I filters included

MyiQ Single-Color
Real-Time PCR
Detection System

Software

Operating system	Windows XP Professional
Multissay analysis	Up to 5 standard curves per plate
Data analysis modes	PCR quantification with standard curve Melt-curve analysis End-point analysis for up to 5 assays Gene expression analysis by relative quantity (ΔC_T) or normalized expression ($\Delta\Delta C_T$) Multifile gene expression analysis for over 5,000 C_T values
Data export modes	Export data directly to Microsoft Excel file format Copy and paste graphs and charts directly into Microsoft Word or PowerPoint files Print all available data graphs and tables from software-based menu Detailed data reports export run conditions, data graphs and tables, and data analysis parameters to printer or destination file



Superb linearity over a wide dynamic range of template starting concentrations. A 10-fold serial dilution of plasmid DNA was amplified using iQ™ supermix with primers and a FAM-labeled probe specific to the IL-1β gene. Four replicates of each template concentration were monitored with the MyiQ real-time system. Standard curve had R² = 0.999, efficiency = 106.8%.

Ordering Information

Catalog #	Description
170-9770	MyiQ Single-Color Real-Time PCR Detection System , includes iCycler base, optics module, software CD-ROM, 96-well optical reaction module, optical-quality 96-well PCR plates, sealing tape, communication cables, power cords, instructions
170-9775	MyiQOptical System Upgrade Kit , includes optics module, software CD-ROM, optical lid for 96-well reaction module, optical-quality 96-well PCR plates, sealing tape, communication cables, power cords, instructions
170-9778	MyiQReal-Time PCR Detection System With Desktop PC , includes iCycler base, optics module, software CD-ROM, 96-well optical reaction module, optical-quality 96-well PCR plates, sealing tape, communication cables, power cords, instructions, desktop computer, monitor
170-9779	MyiQReal-Time PCR Detection System With Laptop PC , includes iCycler base, optics module, software CD-ROM, 96-well optical reaction module, optical-quality 96-well PCR plates, sealing tape, communication cables, power cords, instructions, laptop computer
170-9799	Real-Time PCR Applications Guide
170-8756	Halogen Lamp , replacement, for iCycler iQ®, iQ™5, and MyiQ systems
TBS-0201	0.2 ml 8-Tube Strips Without Caps , natural, 125
TCS-0803	Optical Flat 8-Cap Strips , for 0.2 ml tubes and plates, ultraclear, 120
223-9441	iQ 96-Well PCR Plates , 25
MSB-1001	Microseal 'B' Adhesive Seals , 100

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 applied fields other than Human In Vitro Diagnostics under one or more of U.S. Patents Nos. 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 No. 6,814,934 B1 for use in research and for all other fields except the fields of human diagnostics and veterinary diagnostics.

Excel, Microsoft, PowerPoint, Windows, and Windows XP are trademarks of Microsoft Corporation. SYBR is a trademark of Molecular Probes, Inc.



**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 905 712 2771 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 5811 6270 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