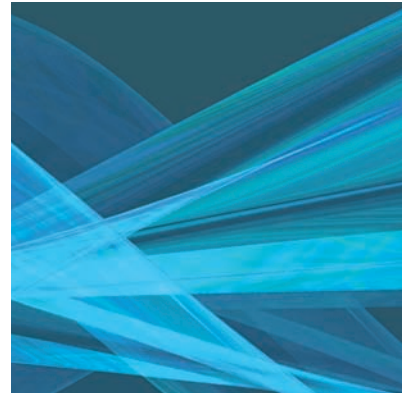


Amplification: MyCycler Thermal Cycler



A Lab's Best Friend

MyCycler™ Thermal Cycler

Exceptional Performance

- Large graphical display
- USB communication
- Self-adjusting heated lid
- Intuitive programming
- Upgradable to gradient capability

Personal Thermal Cycling

- Programmable user preferences
- Selectable power-save mode
- Sleek, compact design

Complete PCR Solutions

- iTaq™ DNA polymerase
- iProof™ high-fidelity DNA polymerase
- iScript™ cDNA synthesis kit
- PCR tubes and plates
- Unsurpassed customer service

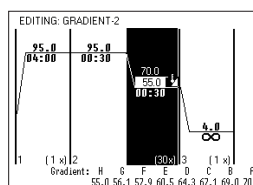
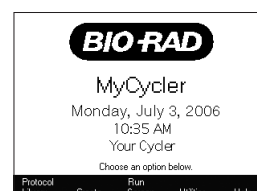
For more information, visit us on the Web at
www.bio-rad.com/products/mycycler/.



BIO-RAD

Features

- Large 12 cm (4.7") high-resolution display
- Programming using intuitive symbols and numerical entry procedures
- Graphical representation of protocols
- Menu-driven software and ready-made templates for ease of viewing, editing, and running protocols
- Innovative thermal gradient design (optional upgrade)
- Reaction module and self-adjusting lid accommodate 96 x 0.2 ml samples in tubes, strips, or plates
- Power-saving programmable automatic standby mode
- USB communication for printing of protocols and reports
- Storage of detailed run and validation reports
- Attractive, compact design



Graphical interface simplifies protocol editing.

System Specifications

Operational

Heating rate	Up to 2.5°C/sec
Cooling rate	Up to 1.5°C/sec
Method of heating/cooling	Peltier
Modes of temperature monitoring	Block and algorithm
Temperature	
Range	4–100°C
Accuracy	±0.5°C
Uniformity	±0.5°C
Overshoot maximum	0.8°C
Heated lid	Up to 110°C
Gradient	
Gradient accuracy	±0.5°C of programmed target at 90°C
Row uniformity	±0.5°C well-to-well within 30 sec of reaching target temperature
Gradient range	30–99°C
Temperature differential range	1–25°C

Descriptive

Sample capacity	96 x 0.2 ml tubes or 96-well PCR plate
Display	12 cm (4.7") high-resolution
Operating temperature	18–28°C ambient
Ports	1 USB
Electrical approvals	IEC, CE
Dimensions (W x D x H)	24 x 44 x 21 cm (9 x 17.5 x 8")
Weight	10 kg (22 lb)

Programming

Memory	99 typical programs
Cycles per protocol	Up to 9
Segments per cycle	Up to 9
Repeats per cycle	Up to 99
Segment time range	00:01–99:59 min
Pause	Yes
Battery-backed RAM	Yes
Onboard software	Menu-driven, real-time graphical display of thermal protocol, intuitive PCR language programming, automatic validation reporting, customizable auditory signals for pause and end of run
Key programming features	Onboard protocol templates for cycle sequencing, touchdown, RT-PCR, nested PCR, and long PCR, auto standby option

Ordering Information

Catalog #	Description
170-9703	MyCycler Thermal Cycler System , includes MyCycler thermal cycler, PCR tubes, iTaq DNA polymerase, 2 sample loading trays, power cord, quick reference guide, instructions
170-9705	MyCycler Firmware Upgrade Kit , includes USB cable, current operating firmware, instructions
170-9707	MyCycler Gradient Upgrade Kit , includes USB cable, current operating firmware, USB dongle activation key, instructions
170-8870	iTaq DNA Polymerase , 5 U/μl, includes 250 U polymerase, 1.25 ml 10x PCR buffer, 1.25 ml 50 mM MgCl ₂ solution
170-8891	iScript cDNA Synthesis Kit , 100 x 20 μl reactions, includes 5x iScript reaction mix, iScript reverse transcriptase, nuclease-free water
172-5301	iProof High-Fidelity DNA Polymerase , 2 U/μl, 100 U, includes 5x reaction buffers, MgCl ₂ solution, DMSO
172-5310	iProof HF Master Mix , 100 x 50 μl reactions, includes 2x master mix, DMSO (for highest fidelity with most templates)
TWI-0201	0.2 ml Tubes With Domed Caps , natural, 1,000
TBC-0802	0.2 ml 8-Tube Strips and Domed Cap Strips , natural, 20 bags of 12 8-tube strips and 12 8-cap strips (1,920 tubes and 1,920 caps)
MSB-1001	Microseal® 'B' Adhesive Seals , 100
TCS-0801	Domed 8-Cap Strips , for 0.2 ml tubes and plates, natural, 120
MLP-9601	Multiplate™ 96-Well Unskirted PCR Plates , natural, 25 plates

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.

This product is 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.

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.



**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