

Selected Instrument Features

- Large high-resolution 4.7" (12 cm) display
- Wide range of user-programmable preferences
- Space-saving design for easy setup and transportation
- USB communication
- Alphanumeric naming for maximum flexibility in naming and identifying user protocols
- Graphical representation of protocols for easy viewing, editing, and running
- Menu-driven software with ready-made templates for quick programming
- Heated lid with adjustable temperature control
- Self-adjusting lid accommodates standard 0.2 ml thin-wall PCR tubes, strips, and 96-well plates
- Programmable, power-saving auto standby option when instrument is idle
- Temperature monitoring and control through instrument algorithm and block measurement modes
- Printing and storage of detailed run and validation reports
- NIST-traceable temperature performance



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

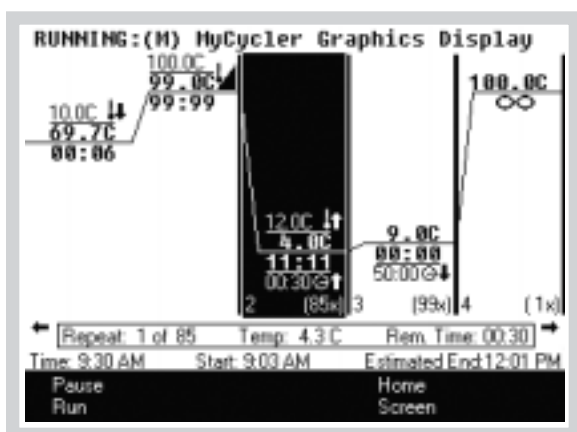
Range	4–100°C
Accuracy	±0.5°C
Uniformity	±0.5°C
Overshoot maximum	0.8°C
Heated lid	Up to 110°C

Descriptive

Footprint	23 x 26.4 x 54.6 cm (9 x 10.4 x 21.5") (H x W x D)
Weight	10 kg (22 lb)
Sample capacity/sample size	96-well x 0.2 ml reaction module
Display	4.7" (12 cm) high-resolution screen
Licensed for PCR	Yes
Operating temperature	18–28°C ambient
Communications	USB
Electrical approvals	IEC, CE

Programming

Maximum number of programs	99 onboard
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
170-9701	MyCycler Thermal Cycler, includes power cord, quick reference guide, instructions
170-9703	MyCycler Thermal Cycler System, includes MyCycler thermal cycler, PCR tubes, iTaq™ DNA polymerase, power cord, quick reference guide, instructions
223-9469	0.2 ml Thin-Wall PCR 8-Tube Strips, natural, 120
223-9472	0.2 ml Thin-Wall PCR 8-Cap Strips, natural, 120
223-9473	0.2 ml Thin-Wall Tubes, natural, 1,000
223-9476	0.2 ml Thin-Wall PCR 12-Tube Strips, natural, 80
223-9477	0.2 ml Thin-Wall PCR 12-Cap Strips, natural, 80
223-9441	96-Well 0.2 ml Thin-Wall PCR Plates, 25
223-9442	92-Well PCR Plate Sealing Mats, 5
223-9444	Optical Quality Sealing Tape, optimized for use with 223-9441, 100 sheets
170-8870	iTaq DNA Polymerase, includes 250 U polymerase, 10x PCR buffer, MgCl ₂ solution
170-8872	MgCl ₂ Solution, 50 mM, 1.25 ml
170-8874	dNTP Solution, 10 mM of each dNTP, 200 µl

Practice of the patented polymerase chain reaction (PCR) process requires a license. The MyCycler thermal cycler is licensed and may be used with PCR licenses available from Applied Biosystems. Its use with authorized reagents also provides a limited PCR license in accordance with the label rights accompanying such reagents. Some applications may require licenses from other parties.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** (800) 4BIORAD **Australia** 02 9914 2800 **Austria** (01)-877 89 01 **Belgium** 09-385 55 11 **Brazil** 55 21 507 6191
Canada (905) 712-2771 **China** (86-21) 63052255 **Czech Republic** (420) 2-4141 0532 **Denmark** 45 44 52-1000 **Finland** 358 (0)9 804 2200
France 01 47 95 69 65 **Germany** 089 318 84-177 **Hong Kong** 852-2789-3300 **India** (91-124) 6398112/113/114, 6450092/93 **Israel** 03 951 4127
Italy 39 02 216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460 **Latin America** 305-894-5950 **Mexico** 52 5 534 2552 to 54
The Netherlands 0318-540666 **New Zealand** 64-9-4152280 **Norway** 47-23-38-41-30 **Poland** (48) 22-8126 672 **Portugal** 351-21-472-7700
Russia 7 095 721 1404 **Singapore** 65-2729877 **South Africa** 00 27 11 4428508 **Spain** 590 5200 **Sweden** 46 (0)8-55 51 27 00 **Switzerland** 061 717-9555
United Kingdom 0800-181134