



Peltier Thermal Cycler

PTC-100®

Features

- Compact footprint
- Precise Peltier-effect heat pumps
- Time-tested reliability
- Adjustable Hot Bonnet® heated lid
- Power-failure protection
- Instant incubation



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

BIO-RAD



First released in 1988, the PTC-100 thermal cycler is a renowned workhorse used by laboratories all over the world. The PTC-100 cycler is available in two sample block formats — 96 x 0.2 ml or 60 x 0.5 ml. It has a fully adjustable heated lid that accommodates the full range of reaction vessels, from low profile to full height, including fully skirted PCR plates.

Programming options include variable ramp rates and incremental changes in incubation time and/or temperature. Advanced file management options include the ability to store programs in password-protected folders. A universal power supply accommodates voltages of 100–240 V, and the power-failure protection feature allows program execution to resume after any interruptions to the power supply.

Specifications

Speed of ramping	60-well: up to 1°C/sec 96-well: up to 1.4°C/sec
Temperature range	0–100°C
Temperature accuracy	±0.5°C of NIST-traceable standard within 30 sec of arrival at 60°C
Temperature uniformity	±0.5°C within 30 sec of arrival at 60°C
Display	20 x 4 alphanumeric LCD
Input power	100–240 VAC, 50–60 Hz, 350 W maximum
Memory	360 typical programs
Operating environment	Ambient temperature: 4–32°C Relative humidity: 10–90%, noncondensing
Heated lid	Hot Bonnet lid included
Dimensions (W x D x H)	24 x 28 x 27 cm
Weight	7.5 kg

Ordering Information

Catalog #	Description
PTC-1196G	PTC-100 Thermal Cycler With Hot Bonnet Heated Lid , holds one 96-well microplate or 96 x 0.2 ml tubes
PTC-1160G	PTC-100 Thermal Cycler With Hot Bonnet Heated Lid , holds 60 x 0.5 ml tubes

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 Corporation. 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.

Appearances and specifications are subject to change without notice.

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 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 4029300/5013478 **Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Mexico** 55 5200 05 20 **The Netherlands** 0318 540666 **New Zealand** 64 9415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 095 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 0861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189/2578 7241 **United Kingdom** 020 8328 2000