

Designed for the Way You Work

Fast

- Get accurate, reliable results with shorter run times and superior thermal performance
- Save time by optimizing annealing temperature in a single run using a temperature gradient

Friendly

- Create protocols quickly with 3 easy programming options: rich graphical, text-based, and automatic with the protocol autowriter
- Get reliable results for years with a patented* O-ring seal that protects thermal electric modules

Flexible

- Choose the best setup for your needs from 4 interchangeable reaction modules, including an optical module for real-time PCR
- Expand the platform as your research grows — control of 2–32 instruments gives you the throughput you need
- Use a wide range of reaction vessels with the fully adjustable heated lid



C1000 Thermal Cycler

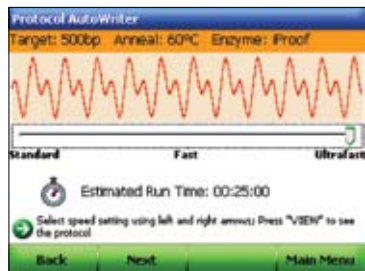
For more information, visit us on the Web at www.bio-rad.com/1000-series/

* U.S. patent 7,051,536.



Writing Protocols Is As Easy As 1-2-3

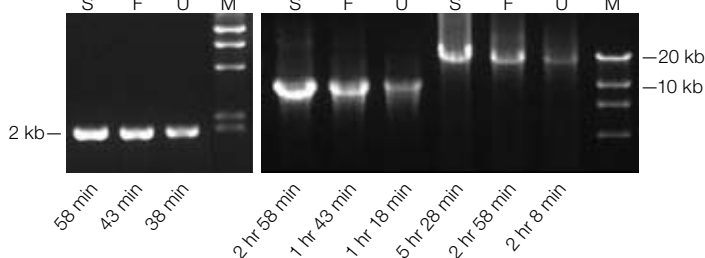
The C1000 thermal cycler's onboard software includes the protocol autowriter — a wizard-based programming option



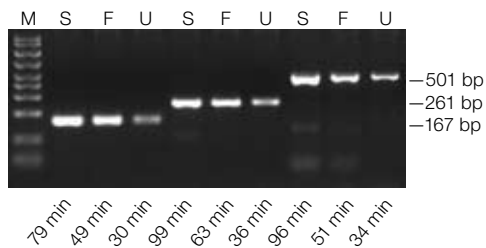
that simplifies protocol creation.

The protocol autowriter automatically suggests a cycling protocol. Enter PCR product length, primer annealing temperature, and polymerase, and select one of three speed settings.

iProof™ High-Fidelity DNA Polymerase



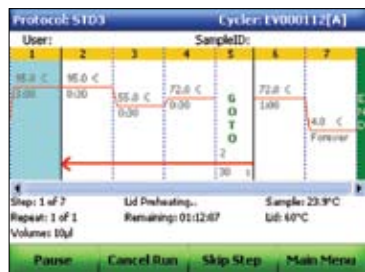
iTaq™ DNA Polymerase



Protocols generated by the protocol autowriter at standard (S), fast (F), and ultrafast (U) settings yield comparable results. To generate fast and ultrafast protocols, the protocol autowriter adjusts annealing temperatures and reduces the total number of protocol steps, so results are obtained in <30 min. M, marker.

Intuitive Graphical Interface

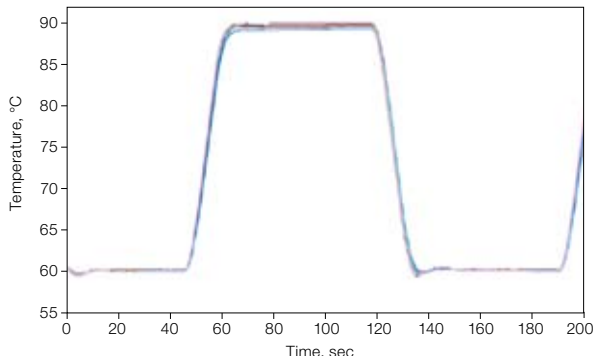
A large full-color display enhances viewing of status, programming screens, and file trees.



Run status view.

Superior Uniformity and Shorter Run Times

The C1000 thermal cycler exhibits high average ramp rates, rapid settling time, and tight thermal uniformity throughout the ramp, resulting in rapid arrival at target temperature and enabling faster protocol run times.



Rapid arrival at target temperature and superior uniformity. Graph shows temperature measured by probes in 15 wells across sample block of the C1000 thermal cycler. Traces are nearly indistinguishable, indicating high uniformity. Note consistent ramp rate throughout heating and cooling.

Ordering Information

Catalog #	Description
184-1000	C1000 Thermal Cycler Chassis , includes USB flash drive, power cord, instructions; does not include reaction module
185-1096	C1000 Thermal Cycler With 96-Well Fast Reaction Module , includes thermal cycler chassis, 96-well fast reaction module, USB flash drive, power cord, reagent and consumable samples, instructions
185-1048	C1000 Thermal Cycler With Dual 48/48 Fast Reaction Module , includes thermal cycler chassis, dual 48/48 fast reaction module, USB flash drive, power cord, reagent and consumable samples, instructions
185-1384	C1000 Thermal Cycler With 384-Well Reaction Module , includes thermal cycler chassis, 384-well reaction module, USB flash drive, power cord, reagent and consumable samples, instructions
184-0048	Dual 48/48 Fast Reaction Module , 2 independent 48-well blocks, fits C1000 and S1000™ thermal cyclers
184-0096	96-Well Fast Reaction Module , fits C1000 and S1000 thermal cyclers
184-0384	384-Well Reaction Module , fits C1000 and S1000 thermal cyclers
184-5096	CFX96™ Optical Reaction Module , includes CFX96 optics shuttle, CFX Manager™ software, communication cables, power cord, reagent and consumable samples, instructions

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's U.S. Patent No. 6,814,934 B1 for use in research and for all other fields except the fields of human diagnostics and veterinary diagnostics. 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