



Advancing PCR Together

PCR Faster

- See 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

PCR Easier

- Easily edit and run protocols using the large color touch screen
- Choose the best setup for your needs from 6 interchangeable reaction modules, including 2 optical modules for real-time PCR
- Increase throughput as your research grows — control of 2–32 instruments gives you the flexibility you need
- Use a wide range of reaction vessels with the fully adjustable heated lid

PCR Smarter

- Quickly create protocols with 2 easy programming options: graphical and automatic with the protocol autowriter
- Get reliable results for years with a patented* O-ring seal that protects thermal electric modules

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

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



Writing Protocols Is As Easy As 1-2-3

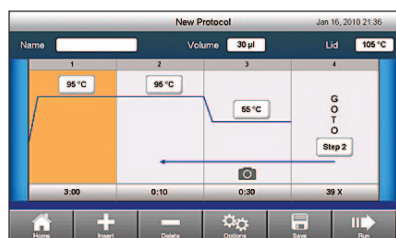
The C1000 Touch thermal cycler's onboard software includes the protocol autowriter — a wizard-based programming option that simplifies protocol creation.



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

Intuitive Graphical Interface

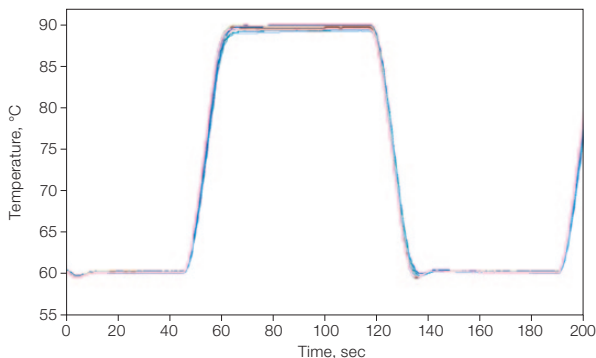
A large full-color touch screen display enhances viewing of status and programming screens.



Run status view.

Superior Uniformity and Shorter Run Times

The C1000 Touch 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 the sample block of a C1000 Touch thermal cycler. Traces are nearly indistinguishable, indicating tight thermal uniformity. Note the consistent ramp rate throughout heating and cooling.

Ordering Information

Catalog #	Description
184-1100	C1000 Touch Thermal Cycler Chassis , includes USB flash drive, power cord; does not include reaction module
185-1148	C1000 Touch Thermal Cycler with Dual 48/48 Fast Reaction Module , includes C1000 Touch thermal cycler chassis, dual 48/48 fast reaction module, USB flash drive
185-1196	C1000 Touch Thermal Cycler with 96-Well Fast Reaction Module , includes C1000 Touch thermal cycler chassis, 96-well fast reaction module, USB flash drive
185-1197	C1000 Touch Thermal Cycler with 96-Deep Well Reaction Module , includes C1000 Touch thermal cycler chassis, 96-deep well reaction module, USB flash drive
185-1138	C1000 Touch Thermal Cycler with 384-Well Reaction Module , includes C1000 Touch thermal cycler chassis, 384-well reaction module, USB flash drive
184-0148	Dual 48/48 Fast Reaction Module , independent dual 48-well reaction module, fits C1000™, C1000 Touch, and S1000™ thermal cyclers, gradient enabled
184-0196	96-Well Fast Reaction Module , fits C1000, C1000 Touch, and S1000 thermal cyclers, gradient enabled
184-0197	96-Deep Well Reaction Module , fits C1000, C1000 Touch, and S1000 thermal cyclers, gradient enabled
184-0138	384-Well Reaction Module , fits C1000, C1000 Touch, and S1000 thermal cyclers, gradient enabled
184-5096	CFX96™ Optical Reaction Module , for use with C1000 Touch thermal cycler chassis, includes CFX Manager™ software, license for qbase ^{PLUS} software, communication cable, reagents, consumables
184-5384	CFX384™ Optical Reaction Module , for use with C1000 Touch thermal cycler chassis, includes CFX Manager software, license for qbase ^{PLUS} software, communication cable, reagents, consumables

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 human in vitro diagnostics and all other applied fields under U.S. Patent Number 5,475,610 (Claims 1, 44, 158, 160-163, and 167 only), or corresponding claims in its non-U.S. counterpart, 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 Number 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

Bio-Rad's thermal cyclers and real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 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 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 11 5044 5699 Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 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 64 9 415 2280 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 Thailand 800 88 22 88 United Kingdom 020 8328 2000