



## Advanced Performance You Can Depend On

### PCR Faster

- See fast run times and accurate, reliable results with superior thermal performance

### PCR Easier

- Meet your current and future needs with modular flexibility — interchangeable modules include dual 48/48-well fast, 96-well fast, 96-deep well, and 384-well reaction modules; all reaction models are gradient enabled
- Use a wide range of reaction vessels with the fully adjustable heated lid

### PCR Smarter

- Expect dependable results for years — patented\* O-ring seal protects thermal electric modules for long-lasting reliability

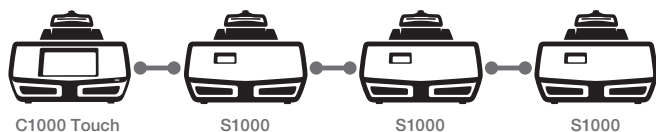
For more information, visit us on the Web at [www.bio-rad.com](http://www.bio-rad.com).

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

**BIO-RAD**

## Flexible Configuration Options

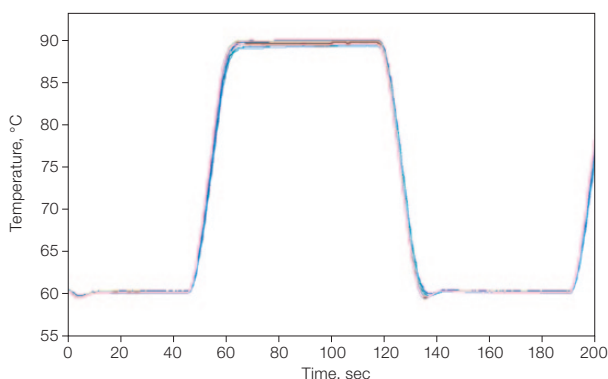
The S1000 thermal cycler can function as a stand-alone instrument or can be connected to a C1000 Touch™ thermal cycler to create a multi-bay instrument.



Expansion capabilities provide the ultimate in high throughput. Up to three S1000 cyclers can be connected to a C1000 Touch cycler to operate as a four-bay instrument. With the addition of a PC and CFX Manager™ software version 2.1 or higher, up to 32 instruments can be operated in tandem.

## Superior Uniformity and Shorter Run Times

The S1000 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 1000-series thermal cycler. Traces are nearly indistinguishable, indicating tight thermal uniformity. Note the consistent ramp rate throughout heating and cooling.

## Thermal Gradient for Optimizing Protocols

A programmable temperature gradient lets you identify optimal annealing temperature for each primer set in a single experiment. Temperature differentials of 1–24°C can be programmed across all reaction module sample blocks.



The thermal gradient feature lets you incubate each row at a different temperature to determine which temperature produces the highest yield and specificity.

## Ordering Information

Catalog #	Description
184-2000	<b>S1000 Thermal Cycler Chassis</b> , includes power cord; does not include reaction module
185-2148	<b>S1000 Thermal Cycler with Dual 48/48 Fast Reaction Module</b> , includes S1000 thermal cycler chassis, dual 48/48 fast reaction module
185-2196	<b>S1000 Thermal Cycler with 96-Well Fast Reaction Module</b> , includes S1000 thermal cycler chassis, 96-well fast reaction module
185-2197	<b>S1000 Thermal Cycler with 96-Deep Well Reaction Module</b> , includes S1000 thermal cycler chassis, 96-deep well reaction module
185-2138	<b>S1000 Thermal Cycler with 384-Well Reaction Module</b> , includes S1000 thermal cycler chassis, 384-well reaction module
184-0148	<b>Dual 48/48 Fast Reaction Module</b> , independent dual 48-well reaction module, fits C1000™, C1000 Touch, and S1000 thermal cyclers, gradient enabled
184-0196	<b>96-Well Fast Reaction Module</b> , fits C1000, C1000 Touch, and S1000 thermal cyclers, gradient enabled
184-0197	<b>96-Deep Well Reaction Module</b> , fits C1000, C1000 Touch, and S1000 thermal cyclers, gradient enabled
184-0138	<b>384-Well Reaction Module</b> , fits C1000, C1000 Touch, and S1000 thermal cyclers, gradient enabled

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 Applera 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.

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.



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://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 31 3689 6600  
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 30 210 777 4396 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 Malaysia 60 3 2117 5260 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 3170 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 66 2 6518311 United Kingdom 020 8328 2000