

## S1000™ Thermal Cycler

Bio-Rad's S1000 thermal cycler offers premium thermal performance, whether operated as a stand-alone instrument or connected to a C1000™ thermal cycler as part of a larger multi-bay configuration.

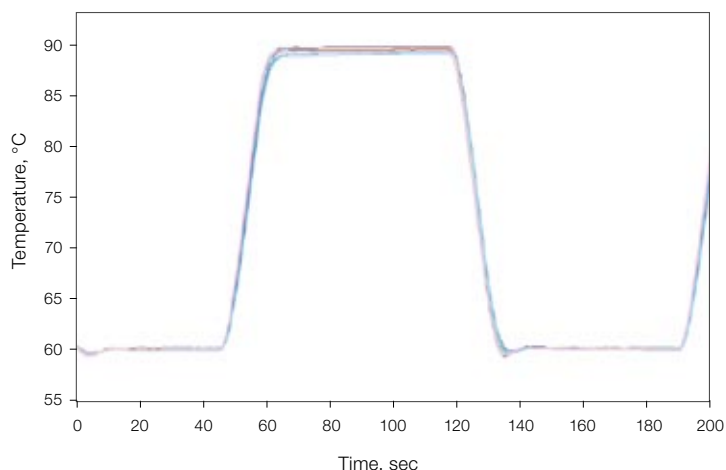
- Innovative engineering delivers exceptional performance and flexibility
- Choice of interchangeable reaction modules includes a gradient-enabled dual 48-well fast module, a gradient-enabled 96-well fast module, and a 384-well module
- Patent-pending thermal electrics and reduced-mass sample block design provide quick time to target temperature for fast protocol run times
- Fully adjustable heated lid accommodates a broad range of vessels
- Optional PC control and networking capability for up to 32 systems enables the ultimate in high throughput



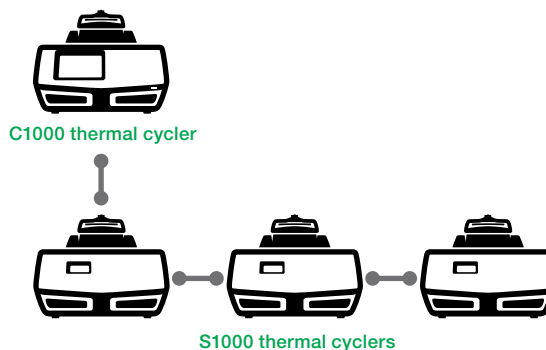
### Specifications

<b>Thermal Cycler</b>		<b>(cont.)</b>	
Input power	400 W, maximum	Temperature control modes	Calculated and block
Frequency	50–60 Hz, single phase	PCR license	Yes
Display	LCD	Programming options	Semigraphical, text-based
Ports	4 USB A, 1 USB B	Security features	Password-protected folders
Fuses	2 6.3 A, 250 V, 5 x 20 mm	PC compatibility	With C1000 cycler; Windows XP or higher
Memory	>1,000 typical programs	Instant incubation	Yes
Dimensions (W x D x H)	13 x 18 x 8"		
Weight	24 lb		
<b>Reaction Modules</b>		<b>Dual 48/48 Fast</b>	<b>384-Well</b>
Sample capacity	<b>96-Well Fast</b> 96 x 0.2 ml tubes	2 x 48 x 0.2 ml tubes	1 x 384-well PCR plate
Maximum ramp rate	5°C/sec	4°C/sec	2.5°C/sec
Average ramp rate	3.3°C/sec	3°C/sec	2°C/sec
Temperature range	0–100°C	0–100°C	0–100°C
Temperature accuracy	±0.2°C of programmed target at 90°C	±0.2°C of programmed target at 90°C	±0.2°C of programmed target at 90°C
Temperature uniformity	±0.4°C well-to-well within 10 sec of arrival at 90°C	±0.4°C well-to-well within 10 sec of arrival at 90°C	±0.4°C well-to-well within 10 sec of arrival at 90°C
Gradient capability	Yes	Yes	No
<b>Gradient*</b>			
Gradient accuracy	±0.2°C of programmed temperature at end rows		
Row uniformity	±0.4°C well-to-well (within row) within 10 sec of arrival at target temperature		
Gradient range	30–100°C		
Temperature differential range	1–24°C		

\* Gradient specifications apply to 96-well and dual 48-well reaction modules.



**Rapid arrival at target temperature and superior uniformity.** Graph shows temperature measured by probes in 15 wells across sample block of a 1000-series thermal cycler. Traces are nearly indistinguishable due to high uniformity. Note the consistent ramp rate throughout heating and cooling. 1000-series thermal cyclers exhibit 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.



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

## Ordering Information

Catalog #	Description
185-2096	<b>S1000 Thermal Cycler With 96-Well Fast Reaction Module</b> , includes thermal cycler chassis, 96-well fast reaction module, power cord, reagent and consumable samples, instructions
185-2048	<b>S1000 Thermal Cycler With Dual 48/48 Fast Reaction Module</b> , includes thermal cycler chassis, dual 48/48 fast reaction module, power cord, reagent and consumable samples, instructions
185-2384	<b>S1000 Thermal Cycler With 384-Well Reaction Module</b> , includes thermal cycler chassis, 384-well reaction module, power cord, reagent and consumable samples, instructions
184-2000	<b>S1000 Thermal Cycler Chassis</b> , includes power cord, instructions; does not include reaction module
184-1000	<b>C1000 Thermal Cycler Chassis</b> , includes USB flash drive, power cord, instructions; does not include reaction module
184-0048	<b>Dual 48/48 Fast Reaction Module</b> , 2 independent 48-well blocks, fits C1000 and S1000 thermal cyclers
184-0096	<b>96-Well Fast Reaction Module</b> , fits C1000 and S1000 thermal cyclers
184-0384	<b>384-Well Reaction Module</b> , fits C1000 and S1000 thermal cyclers
170-8870	<b>iTaq™ DNA Polymerase</b> , 5 U/μl, includes 250 U polymerase, 1.25 ml of 10x PCR buffer (200 mM Tris-HCl, pH 8.4, 500 mM KCl), 1.25 ml of 50 mM MgCl <sub>2</sub> solution
170-8891	<b>iScript™ cDNA Synthesis Kit</b> , 100 x 20 μl reactions, includes 5x iScript reaction mix, iScript reverse transcriptase, nuclease-free water
172-5301	<b>iProof™ High-Fidelity DNA Polymerase</b> , 2 U/μl, 100 U, includes 5x reaction buffers, MgCl <sub>2</sub> solution, DMSO
172-5310	<b>iProof HF Master Mix</b> , 100 x 50 μl reactions, includes 2x master mix (0.04 U/μl), DMSO (for highest fidelity with most templates)
TWI-0201	<b>0.2 ml Tubes With Domed Caps</b> , natural, 1,000
TBC-0802	<b>0.2 ml 8-Tube Strips and Domed Cap Strips</b> , natural, 20 bags of 12 8-tube strips and 12 8-cap strips (1,920 tubes and 1,920 caps)
MSB-1001	<b>Microseal® 'B' Adhesive Seals</b> , 100
TCS-0801	<b>Domed 8-Cap Strips</b> , for 0.2 ml tubes and plates, natural, 120
MLP-9601	<b>Multiplate™ 96-Well 0.2 ml Unskirted PCR Plates</b> , natural, 25 plates

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.

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.

Windows XP is a trademark of Microsoft Corporation.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://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