T100[™] Thermal Cycler

Specifications

Amplification

The T100 thermal cycler is the smart choice for your amplification needs. This modern and compact 96-well PCR instrument offers comprehensive features that make running PCR easier than ever. Start a PCR run or incubation in seconds. Manage your PCR protocols using the large, designed-for-simplicity, color touch-screen user interface. If protocol optimization is required, determine your optimal annealing temperature in a single run using the thermal gradient feature.

- Save time programming with the intuitive touch screen
- Get superior results faster by optimizing your PCR assays in a single run using a thermal gradient
- Save valuable benchspace with the compact design
- Keep your protocols organized using personalized folders or a USB flash drive
- Be confident in your results with the reliability you expect from Bio-Rad



Specifications

Thermal Cycler Input power Display Port Fuses

Memory Dimensions (W x D x H) Weight

Temperature control modes PCR license Programming options Reporting Instant incubation

Performance Sample capacity

Maximum ramp rate Average ramp rate Temperature range Temperature accuracy Temperature uniformity

Thermal Gradient

Gradient capability Gradient range Temperature differential range 100–150 VAC, 50–60 Hz; 220–240 VAC, 50–60 Hz; 700 W maximum 5.7 in. VGA color touch screen 1 USB A Two 10 A, 250 V, 5 x 20 mm 500 typical programs onboard; unlimited with USB flash drive expansion $26 \times 47 \times 23$ cm (10 x 18 x 9 in.) 9 kg (20 lb) Calculated and block Yes Step-based graphical Exportable run logs, system logs Yes

Bulletin 6060

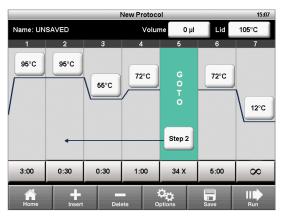
96 x 0.2 ml tubes, 0.2 ml tube strips, or 1 x 96-well plate 4°C/sec 2.5°C/sec 4−100°C ±0.5°C of programmed target ±0.5°C well-to-well within 30 sec of arrival at target temperature

Yes 30-100°C 1-25°C





T100 thermal cycler Home screen.



Intuitive graphical programming. The thermal cycler's onboard software displays an editable thermal profile of the PCR protocol, making it easy to create and run new protocols.

Saved Protocols 11:33		
Folders	Files	Preview
	IPRF15KB	Lid: 105°C Volume: 20 µl
MAIN	IPRF1KB	1. 98°C, 0:30 2. 98°C, 0:05 3. 60°C, 0:10
AVA	IPRF8KB	4. 72°C, 7:30 5. GOTO step 2, 29 X 6. 72°C, 5:00
VSB	ISCRIPT	7. 12°C,∞
	ITAQ-FST	
	~	
Home Folder Op		

Saved Protocols screen showing the preinstalled library of standard protocols in the MAIN folder.

Ordering Information

-	
Catalog #	Description
186-1096	T100 Thermal Cycler, includes 96-well thermal cycler, power
	cord, T100 tube support ring
170-8870	iTaq [™] DNA Polymerase, 5 U/µl, includes 250 U polymerase,
	1.25 ml 10x PCR buffer (200 mM Tris-HCl, pH 8.4, 500 mM
	KCl), 1.25 ml 50 mM MgCl, solution
170-8890	iScript [™] cDNA Synthesis Kit, 25 x 20 µl reactions, includes
	5x iScript reaction mix, iScript reverse transcriptase,
	nuclease-free water
172-5301	iProof [™] High-Fidelity DNA Polymerase, 2 U/µl, 100 U,
	includes 5x reaction buffers, MgCl ₂ solution, DMSO
HSS-9601	Hard-Shell [®] High-Profile 96-Well Semi-Skirted PCR
	Plates, clear shell, clear well, 25
MLP-9601	Multiplate [™] 96-Well Unskirted PCR Plates, clear, 25 plates
MSB-1001	Microseal [®] 'B' Adhesive Seals, optically clear, 100
TBS-1201	12-Tube Strips without Caps (0.2 ml), clear, 100 strips
	(1,200 PCR tubes)
TCS-1201	Domed 12-Cap Strips, for 0.2 ml PCR tubes and plates,
	clear, 20
TWI-0201	PCR Tubes with Domed Caps (0.2 ml), clear, 1,000
. = • ·	

For more information, visit

www.bio-rad.com/web/T100Specifications.



Bio-Rad's 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 iProof and iTaq DNA polymerases 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.

Hard-Shell plates are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 7,347,977; 6,340,589; and 6,528,302.



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 3065 7550 Canada 905 364 3435 China 66 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 361 459 6100 India 91 124 4029300 Israel 03 963 6050 Italy 39 02 216091 Japan 81 3 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 9 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 026 674 55 05 Taiwan 886 2 2578 7189 Thailand 1800 88 22 88 United Kingdom 020 8328 2000