



T100™ Thermal Cycler

The T100 thermal cycler is the smart PCR choice. This compact, 96-well thermal cycler offers a comprehensive package of features, including an easy-to-use touch screen, thermal gradient, and reliable performance. Running PCR experiments has never been easier.

With the T100 thermal cycler, you can:

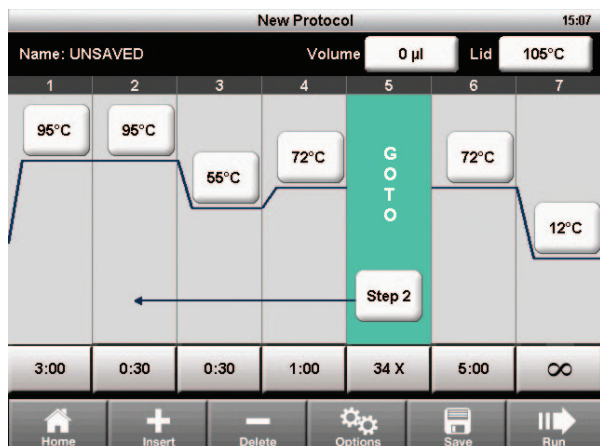
- Save time programming with the intuitive touch screen
- Get superior results faster by optimizing your PCR assays in a single run using a thermal gradient
- Save valuable benchspace with the compact design
- Keep your protocols organized using personalized folders or a USB flash drive
- Be confident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

For more information, visit us on the Web at www.bio-rad.com/T100.

BIO-RAD

Easier Programming

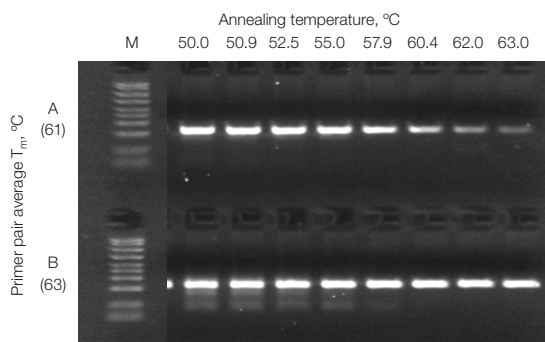
The intuitive touch screen makes it easy to program the T100 thermal cycler. A high-resolution touch screen (5.7") presents a button-driven graphical interface to help you personalize your thermal cycling conditions. It is easy to create a protocol, start a run, or manage files, so you spend less time programming.



Easy graphical programming shows the thermal profile of a new protocol.

Faster Optimization

A programmable temperature gradient lets you identify the optimal annealing temperature for a primer set in a single experiment. Temperature differentials of 1–25°C can be programmed across the sample block. Save research time by quickly optimizing reaction conditions, so you can focus on your results.



Optimization of an assay results in better yields. A superior yield is obtained at the optimal annealing temperature. In primer pair A, a reduced yield is seen above the optimal melting temperature (T_m). In primer pair B, nonspecific products are formed below the optimal T_m . M, markers.

Better Performance

For those who simply need dependable performance, the T100 cycler delivers the reliability you expect from Bio-Rad. You can be confident in the results obtained from your most crucial experiments.



Obtain consistent results with low sample volumes. A 300 bp product was amplified from genomic DNA in 5 μ l reactions. M, markers.

Ordering Information

Catalog #	Description
186-1096	T100 Thermal Cycler , includes 96-well thermal cycler, power cord, T100 tube support ring
170-8890	iScript™ cDNA Synthesis Kit , 25 x 20 μ l reactions, includes 5x iScript reaction mix, iScript reverse transcriptase, nuclease-free water
170-8870	iTaq™ DNA Polymerase , 5 U/ μ l, 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 ₂ solution
172-5301	iProof™ High-Fidelity DNA Polymerase , 2 U/ μ l, 100 U, includes 5x reaction buffers, MgCl ₂ solution, DMSO
MLP-9601	Multiplate™ 96-Well Unskirted PCR Plates , natural, 25 plates
MSB-1001	Microseal® 'B' Adhesive Seals , optically clear, 100
TBS-1201	12-Tube Strips without Caps (0.2 ml) , natural, 100 strips (1,200 PCR tubes)
TCS-1201	Domed 12-Cap Strips , for 0.2 ml PCR tubes and plates, natural, 200
TWI-0201	PCR Tubes with Domed Caps (0.2 ml) , natural, 1,000

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.

This product is 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

**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 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