



Thermal Cycling

Modern. Intuitive. Robust. PCR for Your Biopharmaceutical Laboratory

The Reliability You Depend on with Streamlined User Experience and Connectivity for Your Modern Controlled Lab

Engineered to maintain thermal performance, the PTC Tempo Thermal Cycler offers the flexibility to scale with your laboratory. This conventional PCR thermal cycler features an intuitive user interface and flexible connectivity options for ease of protocol management. A motorized lid enables automation, taking your PCR to the next level with modern instrument control.

Consistent Performance

PTC Tempo Thermal Cyclers offer thermal accuracy and uniformity you can trust to get the consistent performance you need.

Intuitive Experience

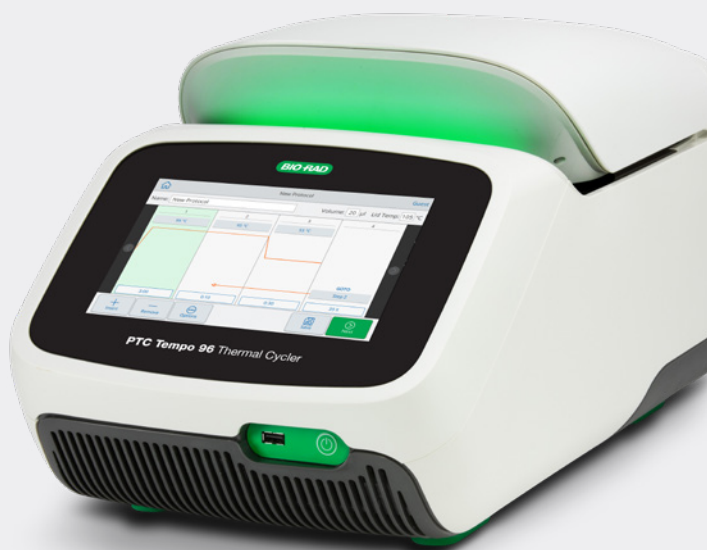
The modern user interface with enhanced touch sensitivity is intuitive for experienced and new users alike — perfect for training labs and scientists transitioning from other platforms. Additionally, PCR protocols can be stored and accessed when connected to your network drive, and the PTC Tempo Thermal Cycler connects with the BR.io cloud platform via WiFi to utilize cloud-enabled features such as remote protocol setup and instrument monitoring.

Flexible Service Options

Installation Qualification (IQ), Operational Qualification (OQ), Thermal Qualification (TQ), and preventive maintenance services are available to fit your needs.

Instant Usability

Out of the box, PTC Tempo Thermal Cyclers are factory calibrated and ready to use.



Automation Efficiency

The motorized lid of the PTC Tempo 96, PTC Tempo Deepwell, and PTC Tempo 384 Thermal Cyclers increases efficiency and ensures consistent results. This self-adjusting lid removes the need to turn knobs and levers. In addition, the PTC Tempo Thermal Cycler API Reference Guide provides information on integration with robotics systems.

Secure User Mode

User access is protected to enhance security in regulated environments.


Search your mobile app store for **Bio-Rad AR** to get an augmented reality tour. Contact your local sales representative for more information.



Visit bio-rad.com/PTCTempo for more information.

BIO-RAD

Ordering Information

Catalog #	Description	Catalog #	Description
Instruments			
12015382	PTC Tempo 96 Thermal Cycler	12001925	ddPCR™ 96-Well Plates, semi-skirted, clear/clear, 25
12015392	PTC Tempo Deepwell Thermal Cycler	MSA5001	Microseal™ 'A' PCR Plate and PCR Tube Sealing Film, nonadhesive sealing film, 50
12015309	PTC Tempo 48/48 Thermal Cycler	MSB1001	Microseal 'B' PCR Plate Sealing Film, adhesive, optical, 100
12015394	PTC Tempo 384 Thermal Cycler	MSC1001	Microseal 'C' PCR Plate Sealing Film, adhesive, optical, 100
1814000	PX1 PCR Plate Sealer	MSF1001	Microseal 'F' PCR Plate Seal, foil, pierceable, 100
Accessories			
12018650	PTC Tempo Thermal Cycler WiFi Adapter, U.S. only	1814030	PCR Plate Heat Seal, clear, optical, 100
12018509	PTC Tempo Thermal Cycler WiFi Adapter, global	1814040	PCR Plate Heat Seal, foil, pierceable, 100
1849000	Tube Frame for 0.2 ml Tubes, for use with 48/48- and 96-well reaction modules	1814045	PCR Plate Heat Seal, foil, peelable, 100
1849001	Tube Frame for 0.2 ml Tubes, for use with 96-deep well and deepwell reaction modules	Agaroses	
12020274	Tube Frame for 0.5 ml Tubes, for use with 96-deep well and deepwell reaction modules	1613102	Certified Molecular Biology Agarose, 500 g
12014294	Plate Lifter for 384-Well Reaction Module	1613107	Certified Low Range Ultra Agarose, 125 g
12019776	PTC Tempo Thermal Cycler API, 1 license	1613112	Certified Low Melt Agarose, 125 g, for DNA fragments ≥1,000 bp
12013205	Ethernet Cable	1613115	Certified PCR Low Melt Agarose, 500 g, for DNA fragments ≤1,000 bp
Reagents			
1725310	iProof HF Master Mix, 100 x 50 µl reactions, 2.5 ml	1613105	Certified PCR Agarose, 500 g, for DNA fragments ≤1,000 bp
1725320	iProof GC Master Mix, 100 x 50 µl reactions, 2.5 ml	Buffers	
1708870	iTaq DNA Polymerase, 250 U (5 U/µl), 50 µl	1610770	10x TBE (Tris/Boric Acid/EDTA) Buffer, 5 L
1708872	MgCl ₂ Solution for PCR, 50 mM, 1.25 ml	1610733	10x TBE (Tris/Boric Acid/EDTA) Buffer, 1 L
1708874	dNTP Solution Mix, 10 mM, 200 µl	1610773	50x TAE (Tris/Acetic Acid/EDTA) Buffer, 5 L
12012802	Reliance Select cDNA Synthesis Kit, 25 x 20 µl reactions	1610743	50x TAE (Tris/Acetic Acid/EDTA) Buffer, 1 L
1708840	iScript Reverse Transcription Supermix, 25 x 20 µl reactions	1610433	Ethidium Bromide Solution, 10 mg/ml
1725034	iScript gDNA Clear cDNA Synthesis Kit, 25 x 20 µl reactions	1610729	EDTA, 500 g, ≥99% pure
1725037	iScript Advanced cDNA Synthesis Kit, 25 x 20 µl reactions	1610716	Tris, 500 g, ≥99.8% pure
1708896	iScript Select cDNA Synthesis Kit, 25 x 20 µl reactions	Horizontal Electrophoresis	
1708890	iScript cDNA Synthesis Kit, 25 x 20 µl reactions	1704486	Mini-Sub™ Cell GT Horizontal Electrophoresis System
Plastics			
TLS0801	0.2 ml 8-Tube PCR Strips without Caps, low profile, clear, 120	1640300	Mini-Sub Cell GT Horizontal Electrophoresis System and PowerPac Basic Power Supply
TFI0201	0.2 ml PCR Tubes with Flat Caps, high profile, clear, 1,000	1704485	Wide Mini-Sub Cell GT Horizontal Electrophoresis System
TWI0201	0.2 ml PCR Tubes with Domed Caps, high profile, clear, 1,000	1640301	Wide Mini-Sub Cell GT Horizontal Electrophoresis System and PowerPac Basic Power Supply
TBS0201	0.2 ml 8-Tube PCR Strips without Caps, high profile, clear, 125	1704482	Sub-Cell™ GT Horizontal Electrophoresis System
TBC0802	0.2 ml 8-Tube PCR Strips and Domed Cap Strips, high profile, clear	1640302	Sub-Cell GT Horizontal Electrophoresis System and PowerPac Basic Power Supply
TCS0803	0.2 ml Flat PCR Tube 8-Cap Strips, optical, ultraclear	Rulers	
HSP9601	Hard-Shell™ 96-Well PCR Plates, low profile, thin wall, skirted, white/clear, 50	1708351	EZ Load 20 bp Molecular Ruler, 20–1,000 bp, 50 bands
MLL9601	Multiplate 96-Well PCR Plates, low profile, unskirted, clear, 25	1708352	EZ Load 100 bp Molecular Ruler, 100–1,000 bp, 10 bands
HSL9601	Hard-Shell 96-Well PCR Plates, low profile, semi-skirted, clear/clear, 25	1708353	EZ Load 100 bp PCR Molecular Ruler, 100–3,000 bp, 30 bands
HSL9605	Hard-Shell 96-Well PCR Plates, low profile, semi-skirted, clear/white, 25	1708354	EZ Load 500 bp Molecular Ruler, 500–8,000 bp, 16 bands
HSL9905	Hard-Shell 96-Well PCR Plates, low profile, semi-skirted, clear/white, barcoded, 25	1708355	EZ Load 1 kb Molecular Ruler, 1–15 kb, 15 bands
HSS9601	Hard-Shell 96-Well PCR Plates, high profile, semi-skirted, clear/clear, 25	1708200	AmpliSize™ Molecular Ruler, 50–2,000 bp, 10 bands
HSS9901	Hard-Shell 96-Well PCR Plates, high profile, semi-skirted, clear/clear, barcoded, 25	Imaging	
MLP9601	Multiplate 96-Well PCR Plates, high profile, unskirted, clear, 25	12009077	GelDoc Go Gel Imaging System with Image Lab Touch Software, includes UV/Stain-Free Imaging Tray
MLL4801	Multiplate 48-Well PCR Plates, low profile, unskirted, clear, 50	1660531	UView Mini Transilluminator
MLP4801	Multiplate 48-Well PCR Plates, high profile, unskirted, clear, 50	Visit bio-rad.com/PTCTempo for more information about the PTC Tempo Thermal Cycler.	
HSP3801	Hard-Shell 384-Well PCR Plates, thin wall, skirted, clear/clear, 50	Visit BR.io for more information about the BR.io cloud platform.	
HSP3901	Hard-Shell 384-Well PCR Plates, thin wall, skirted, clear/clear, barcoded, 50	 <p>Bio-Rad PCR reagents and analytical instruments are manufactured under an ISO 13485:2016 certified Quality Management System and are quality control tested to ensure consistent product performance you can trust.</p>	



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 080 007 7373 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

