



Thermal Cycling

Modern. Intuitive. Robust. PCR for Your Research Laboratory

The Reliability You Depend on with Streamlined User Experience and Connectivity for Your Modern 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 the thermal accuracy and uniformity you can trust to get the consistent performance you need.

Intuitive Experience

The modern user interface with enhanced touch sensitivity and the automated lid of the PTC Tempo 96, Deepwell, and 384 Thermal Cyclers are intuitive for experienced and new users alike — perfect for training labs and scientists transitioning from other platforms.

Instant Usability

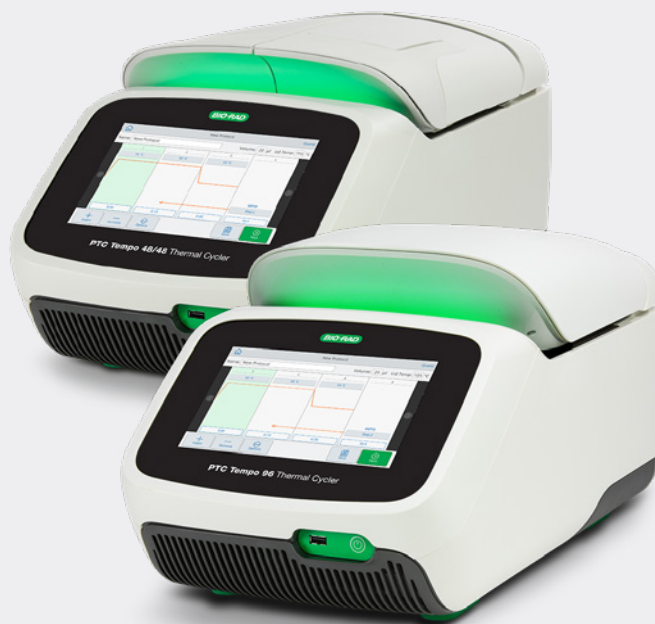
Out of the box, PTC Tempo Thermal Cyclers are factory calibrated and ready to use.

Economical Service Options

Thermal qualification and preventive maintenance services are available to fit your needs.

Mutli-Block Flexibility

The PTC Tempo 48/48 Thermal Cycler, with a non-automated lid, can simultaneously run two independent protocols.



Contemporary Connectivity

WiFi connectivity allows you to place the instrument anywhere in the lab. PCR protocols can be stored and accessed when connected to your network drive, and system logs and run reports can be downloaded via USB to store on your network. The PTC Tempo Thermal Cycler can be connected with the BR.io cloud platform to utilize cloud-enabled features such as remote protocol setup and instrument monitoring.

Visible Status

The instrument status light is visible at a glance from a distance, allowing you to walk away with confidence.

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 |  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. | |



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 Poland 00 800 00 24 67 23 Portugal 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

