

Trusted Thermal Cyclers for NGS Library Preparation

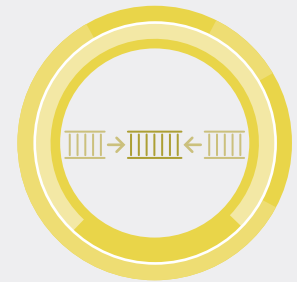
Next-generation sequencing (NGS) library preparation workflows can include several steps, such as RNA/DNA extraction, RNA integrity analysis, fragmentation, adapter sequence addition, cDNA synthesis, target enrichment, and amplification. Robust thermal cyclers, such as the PTC Tempo and PTC Harmony Thermal Cyclers, are a vital part of the NGS workflow. Bio-Rad thermal cyclers utilize Peltier technology for precise temperature control and thermal gradient technology for simplified PCR optimization.

- **Need flexibility?**
Utilize the PTC Tempo 48/48 Thermal Cycler with dual reaction blocks to simultaneously run 2 independent protocols
- **Need higher reaction volumes?**
Consider the PTC Tempo Deepwell Thermal Cycler or the PTC Harmony Deepwell Thermal Cycler to accommodate volumes up to 125 μ L, perfect for NGS library prep volume requirements
- **Need reliability?**
 - Reproducible results, due to rapid arrival at target temperature and superior uniformity
 - Trusted thermal cyclers for more than 35 years, worldwide
 - Bio-Rad thermal cyclers have been recommended in protocols by leading sequencing companies
- **Need high throughput?**
Drive sample throughput and minimize errors with the automation-compatible PTC Tempo Deepwell Thermal Cycler

Make Bio-Rad thermal cyclers a part of your NGS library prep workflow.



Fragmentation



Adapter Sequence Addition



Size Selection
(electrophoresis)



In addition to thermal cyclers, Bio-Rad offers these essential tools for the NGS library preparation workflow.

Explore the SEQuoia™ Workflow for Revolutionized RNA Sequencing



Featuring the latest advancements in enzyme technology, SEQuoia products offer superior construction of RNA sequencing libraries from low-input and low-quality samples. The SEQuoia one-tube continuous synthesis library prep reaction minimizes pipetting steps and shortens the workflow. The SEQuoia workflow can greatly reduce the total cost of library preparation without compromising data integrity or quality.

Visit bio-rad.com/NGS to learn more.

Electrophoresis and Plastic Consumables Ideal for NGS Library Preparation

Select fragment sizes and evaluate RNA integrity with horizontal electrophoresis products



Sub-Cell™ GT Cells are versatile and enable multisample, high-throughput nucleic acid applications.

Save preparation time and standardize electrophoresis runs with premixed electrophoresis running buffers that are compatible with handcast or precast gels.

Visit bio-rad.com/DNAElectro to learn more.

Ensure sequencing success with Hard-Shell™ PCR Plates and Microseal™ PCR Plate Seals



Hard-Shell PCR Plates are precisely manufactured for optimal fit and cycling performance.

Adhesive foils and optically clear seals are designed with a strong adhesive that minimizes sample evaporation.

Visit bio-rad.com/PCRPlastics to learn more.

Ordering Information

Catalog #	Description
12015392	PTC Tempo Deepwell Thermal Cycler
12015309	PTC Tempo 48/48 Thermal Cycler
12024257	PTC Harmony Deepwell Thermal Cycler
12009077	GelDoc™ Go Gel Imaging System with Image Lab Touch Software
1640301	Wide Mini-Sub™ Cell GT Horizontal Electrophoresis System
1814000	PX1 PCR Plate Sealer
17005710	SEQuoia Complete Stranded RNA Library Prep Kit, 96 reactions
12017265	SEQuoia Express Stranded RNA Library Prep Kit, 96 reactions
17006487	SEQuoia RiboDepletion Kit, 24 reactions
12011930	SEQuoia Dual Indexed Primers Plate
MSB1001B	Microseal 'B' PCR Plate Sealing Film, adhesive, optical, 100
MSF1001B	Microseal 'F' PCR Plate Seal, foil, pierceable, 100
HSP9601	Hard-Shell 96-Well PCR Plates, low profile, thin wall, skirted, white shell, clear well, 50
TFI0201	0.2 mL PCR Tubes with Flat Caps, high profile, clear, 1,000
TWI0201	0.2 mL PCR Tubes with Domed Caps, high profile, clear, 1,000

Catalog #	Description
MLP4801	Multiplate 48-Well PCR Plates, high profile, unskirted, clear, 50
MLL4801	Multiplate 48-Well PCR Plates, low profile, unskirted, clear, 50

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



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, GelDoc, Hard-Shell, Microseal, Mini-Sub, SEQuoia, and Sub-Cell are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2025 Bio-Rad Laboratories, Inc.

The Mini-Sub Cell GT, Wide Mini-Sub Cell GT, and Sub-Cell GT Cell Systems and/or their use are covered by the following U.S. patent and/or pending U.S. patent application or their foreign counterparts owned by or under license to Bio-Rad Laboratories, Inc., including, but not limited to, U.S. Patent No. 8,262,889.

Image Lab Touch Software is based in part on the work of the Cimg project (<http://cimg.eu/>). See license for details at www.cecill.info/licences/Licence_CeCILL-C_V1-en.html.

Image Lab Touch is based in part on libraries from GCC runtime and the Gnu C library.



**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 1 800 361 1808 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