

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
1640301	Wide Mini-Sub™ Cell GT Horizontal Electrophoresis System , includes 15- and 20-well combs, gel caster, 15 x 10 cm UV-transparent tray, PowerPac Basic Power Supply
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
MSB1001	Microseal 'B' PCR Plate Sealing Film , adhesive, optically clear, 100
MSF1001	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
HSP9655	Hard-Shell 96-Well PCR Plates , low profile, thin wall, skirted, white shell, white well, 50

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 for more information.

BIO-RAD, 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. © 2023 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.



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

