



A Trusted Thermal Cycler for NGS Library Preparation

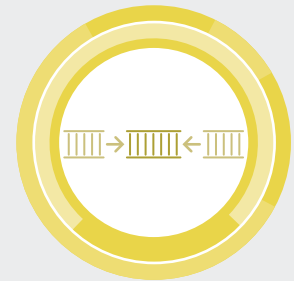
NGS library preparation workflows can include several steps, such as RNA/DNA extraction, RNA integrity analysis, fragmentation, adapter sequence addition, cDNA synthesis, target enrichment, and amplification. Robust thermal cyclers, like the C1000 Touch Thermal Cycler, are a vital part of the NGS workflow. Bio-Rad's thermal cyclers utilize Peltier technology for precise temperature control and thermal gradient technology for simplified PCR optimization.

- **Need flexibility?**
Minimize footprint with the interchangeable blocks of the C1000 Touch Thermal Cycler
- **Need higher volumes?**
Consider the C1000 Touch Thermal Cycler with 96-Deep Well Reaction Module
- **Need reliability?**
 - Rapid arrival at target temperature and superior uniformity produces reproducible results
 - Trusted thermal cyclers for over 25 years, worldwide
 - Recommended in sequencing protocols by leading sequencing companies

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



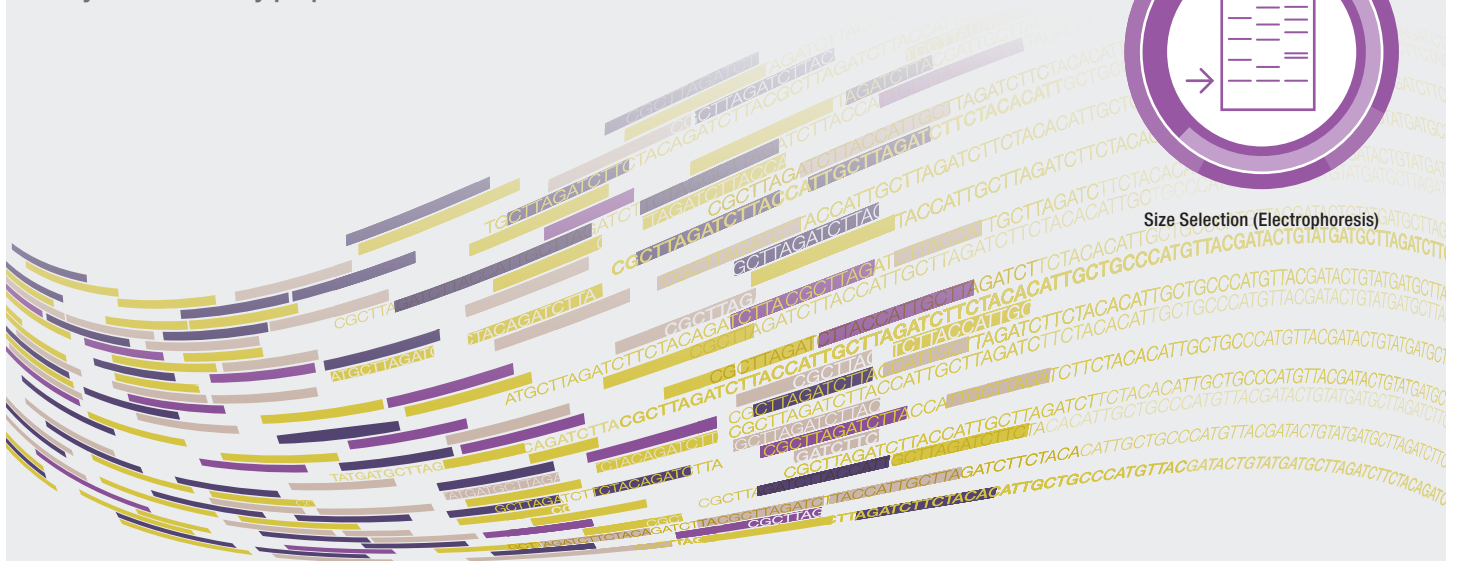
Fragmentation



Adapter Sequence Addition



Size Selection (Electrophoresis)



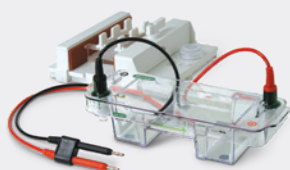
The Best Performance with the Most Flexibility



- **Easy protocol editing** — large color touch screen makes editing and running protocols easier
- **Flexible platform** — increase throughput capability by linking up to 3 additional S1000 Thermal Cyclers or add PC control for up to 32 cyclers
- **Convenient file management** — keep protocols organized using personalized folders or a USB flash drive
- **Thermal gradient** — use the programmable temperature gradient to quickly identify optimal annealing temperatures

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 multi-sample, high-throughput nucleic acid applications.

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

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



Ensure sequencing success with Hard-Shell PCR Plates and 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
1851197	C1000 Touch Thermal Cycler with 96–Deep Well Reaction Module
1851196	C1000 Touch Thermal Cycler with 96-Well Fast Reaction Module
1851148	C1000 Touch Thermal Cycler with Dual 48/48 Fast Reaction Module
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

Catalog #	Description
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

Visit bio-rad.com/C1000touch for more information.

BIO-RAD, HARD-SHELL, MICROSEAL, MINI-SUB, and SUB-CELL are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

Hard-Shell Plates are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 7,347,977; 6,340,589; and 6,528,302. All trademarks used herein are the property of their respective owner.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00 **Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30 **Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580 **Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 01923 47 1301

