



Trusted PCR Plates and Seals for NGS Library Preparation

When nothing should be left to chance, choose Bio-Rad

Countless hours, precious samples, and an immeasurable desire for scientific advancement go into every next-generation sequencing (NGS) project. Why trust Bio-Rad PCR plates and seals?

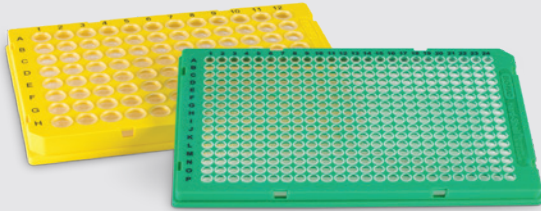
- Trusted genomics research products for more than 20 years
- Reproducibility ensured by rigorous quality control procedures
- Recommended in sequencing protocols by leading sequencing companies
- All plates and seals verified to be RNase-, DNase-, and human DNA-free

Visit [bio-rad.com/PCRplastics](https://www.bio-rad.com/PCRplastics) to learn more.

BIO-RAD

Ensure Sequencing Success with Hard-Shell PCR Plates

Hard-Shell PCR Plates are precisely manufactured for optimal fit and cycling performance, ensuring reproducibility in highly sensitive NGS library preparation protocols.



- **Maximum sample recovery** – thin-walled wells molded of virgin polypropylene facilitate low DNA binding
- **Easy well identification** – black alphanumeric labeling facilitates sample loading
- **Automation compatible** – patented rigid design provides superior stability and flatness
- **Broad cyclers compatibility** – plates specifically designed for optimal fit in thermal cyclers from all major brands

PCR Seals Ideal for NGS Library Preparation

Our foils and optically clear seals have a strong adhesive that:

- Minimizes sample evaporation
- Peels off easily for simple sample retrieval

Optically Clear Adhesive Seals: Microseal 'B' PCR Plate Seals



- **Optical seal** – same seal can also be used for qPCR, including qPCR-based library quantification assays
- **Widely adopted for research** – Bio-Rad's most popular seal

Foil Adhesive Seals: Microseal 'F' PCR Plate Seals



- **Ideal for sample storage** – highly solvent resistant and suitable for storage down to -80°C
- **Pierceable aluminum** – easily retrieve sample by piercing with pipet tip

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

Bio-Rad, Hard-Shell, and Microseal are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

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.

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

