

# Permanent Clear Heat Seal

Catalog #	Description
1814035	<b>Permanent Clear Heat Seal</b> , package of 100

---

**For research purposes only.**

---

The Permanent Clear Heat Seal provides the strongest Bio-Rad heat sealing option and is ideal for applications requiring secure seals, such as water bath cycling. This seal is recommended for real-time PCR. The seal is nonpeelable and nonpierceable.

**Features**

- Forms a permanent seal with polypropylene materials
- Has seal integrity from –80 to 110°C with pressurized heated PCR lid
- Demonstrates high solvent resistance
- Is nonhazardous
- Can be stored at room temperature; avoid direct exposure to light
- Is certified RNase- and DNase-free

**Directions for Use**

**Note:** Optimize sealing parameters in an empty plate prior to completing experimental workflows.

1. Place a microplate in the plate support block and position it correctly on the open PX1 PCR Plate Sealer drawer.
2. Carefully inspect the orientation of the seals inside the bag. Once seals are removed from the packaging, the sealing surface cannot be distinguished from the nonsealing surface.
3. Remove a seal and place the sealing side down against the plate, as described on the bag.
4. **Gently place the sealing frame on top of the seal** and align the vertical supports in the grooves on the drawer.
5. Seal the plate. For suggested sealing parameters, refer to Sealing Guidelines for Permanent Clear Heat Seal.

**Sealing Guidelines for Permanent Clear Heat Seal**

Product	Catalog Number	Number of Wells	Well Color	Temperature, °C	Time, sec
Hard-Shell™ 384-Well Standard PCR Plates	HSP3xxx	384	Clear	175	3
			White	167	3
Hard-Shell 384-Well 480 PCR Plates	HSR4xxx	384	Clear	175	3
			White	167	3
Hard-Shell Low-Profile 96-Well Skirted PCR Plates	HSP9xxx	96	Clear	180	3
			White	180	3
Hard-Shell High-Profile 96-Well Semi-Skirted PCR Plates	HSS9xxx	96	Clear	165	3
			White	180	3
Multiplate Low-Profile 96-Well Unskirted PCR Plates	MLL9xxx	96	Clear	180	3
			White	160	3
Multiplate High-Profile 96-Well Unskirted PCR Plates	MLP9xxx	96	Clear	150	3
			White	165	3
iQ 96-Well PCR Plates	2239441	96	Clear	180	3

**Heat Sealing Film Comparison Chart**

Product	Catalog Number	Key Features	Compatibility
Optically Clear Heat Seal	1814030	<ul style="list-style-type: none"> <li>Validated for real-time PCR</li> <li>Peelable</li> </ul>	PP, PE,* PS
Permanent Clear Heat Seal	1814035	<ul style="list-style-type: none"> <li>Strongest Bio-Rad heat sealing option</li> <li>Ideal for water bath cycling</li> <li>Validated for real-time PCR</li> </ul>	PP
Pierceable Foil Heat Seal	1814040	<ul style="list-style-type: none"> <li>Validated for use with the QX200™, QX200 AutoDG™, QX600™, and QX600 AutoDG Droplet Digital™ PCR Systems</li> <li>Validated for PCR</li> </ul>	PP, PS*
Peelable Foil Heat Seal	1814045	<ul style="list-style-type: none"> <li>Ideal for sample storage</li> <li>Validated for PCR</li> </ul>	PP, COC*

\* Limited and variable compatibility based on plastic composition.

COC, cyclic olefin copolymer; PE, polyethylene; PP, polypropylene; PS, polystyrene.

BIO-RAD, AUTODG, DROPLET DIGITAL, DROPLET DIGITAL PCR, HARD-SHELL, QX200, and QX600 are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2024 Bio-Rad Laboratories, Inc.

Purchase of Digital PCR and/or Single-Cell NGS Sample Preparation products (the "Products") from Bio-Rad Laboratories is subject to Bio-Rad Laboratories, Inc. Standard Terms and Conditions of Sale, which can be accessed at <https://www.bio-rad.com/en-us/terms-conditions>. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for you to distribute or resell the Products. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for the development or commercialization of diagnostic assays for use with the Products without a license from Bio-Rad. It is the user's obligation to obtain a commercial license from Bio-Rad for (i) all commercial uses (not just diagnostic uses) and (ii) sale of assays for use on Bio-Rad's dPCR and ddSEQ instruments. The Products and/or their use are covered by U.S. and foreign patents and/or pending patent applications owned by or under license to Bio-Rad Laboratories, Inc. See <https://www.bio-rad.com/en-us/trademarks>.