



PX1™ PCR Plate Sealer and Heat Sealing Consumables



Easy and Reliable Plate Sealing



User Friendly

Quick and easy, from programming to sealed plate

- Intuitive touch-screen interface
- Quick sealing with the touch of a button
- Optimized sealing protocols for easy storage and use

Fastest Heat Up Time

Avoid delaying an experiment waiting for your sealer to heat up. The innovative heater design allows the PX1 PCR plate sealer to reach sealing temperature in less than 3 minutes.



Bio-Rad PX1



Competitor A



Competitor B

Eliminate Effects of Sample Evaporation from All Your PCR Workflows



Experimental setup



Seal your plate with the touch of a button



Real-time PCR



PCR



Droplet Digital™ PCR

Confidence in Your Results

Minimized sample evaporation improves reproducibility and productivity.



Trust your results

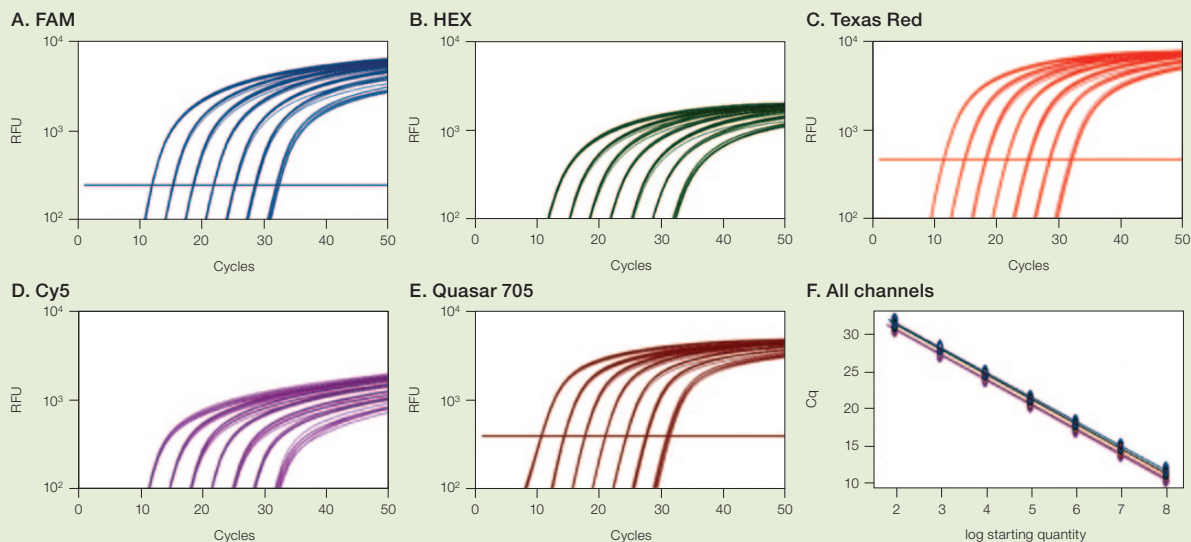
Avoid complications in data interpretation due to sample evaporation.

Avoid repeating experiments

Minimize the need to repeat experiments resulting from sample loss.

Save time

Use fully validated Bio-Rad heat seals and PCR plastics to easily integrate the PX1 PCR plate sealer into your workflow.



The optically clear heat seal prevents sample evaporation that can lead to large Cq standard deviations. Five-color multiplex linearity test demonstrates highly reliable qPCR results. **A–E**, fluorescence data from a series of tenfold dilutions of plasmid DNA (10^8 – 10^2 copies) amplified using reporter dyes to monitor five targets. The standard curves of five multiplexed genes have an $R^2 > 0.990$. **F**, standard curves generated from data in A–E. Reaction efficiencies are between 90 and 105%. Cq, quantification cycle; RFU, relative fluorescence units.

A Complete Selection of Heat Seals

Bio-Rad heat seals help deliver reliable data by minimizing sample evaporation. Choose from a selection of fully validated seals ideal for a broad range of PCR applications.



Optically Clear Heat Seal

- Ideal for real-time PCR
- Excellent optical clarity
- Peelable for easy sample retrieval
- Compatible with PCR



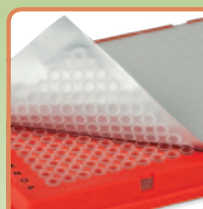
Permanent Clear Heat Seal

- Ideal for water bath cycling
- Compatible with PCR
- High solvent resistance



Pierceable Foil Heat Seal

- Fully validated for the QX100™ Droplet Digital™ PCR system workflow
- Easily pierceable with a pipet tip
- User friendly — colored stripe clearly identifies sealing surface
- Compatible with PCR



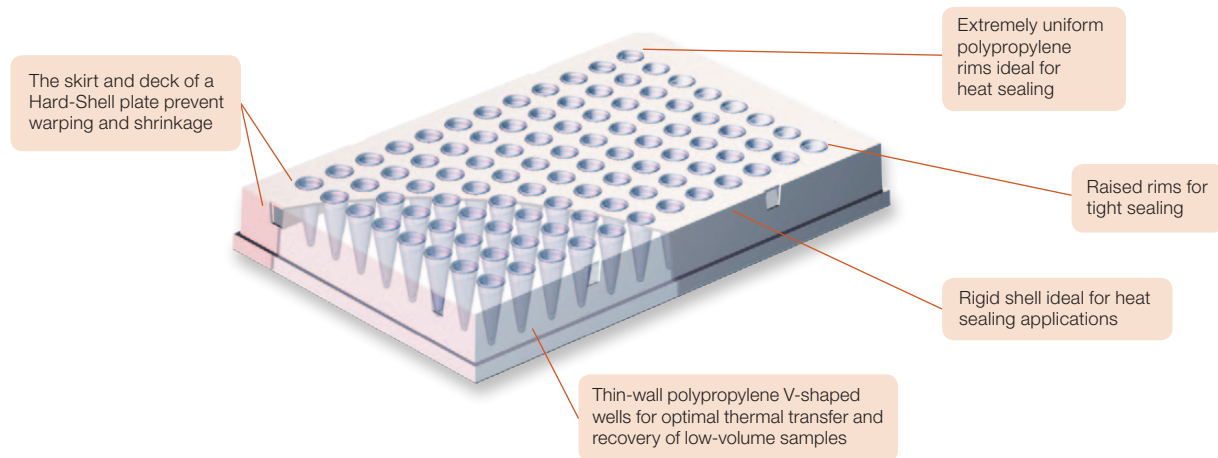
Peelable Foil Heat Seal

- Ideal for low-temperature sample storage
- Can be easily peeled from microplates stored in a -80°C freezer or in liquid nitrogen
- Compatible with PCR

Hard-Shell® Plates Are Ideal for Heat Sealing

Hard-Shell PCR plates are specifically designed to withstand the stresses of heat sealing, thermal cycling, and robotic handling. Hard-Shell PCR plate features include:

- Rigid polycarbonate shell provides superior stability and flatness that aids in the heat sealing process
- Fully validated for use with the PX1 PCR plate sealer
- Patented Hard-Shell technology prevents plate distortion during thermal cycling
- Thin-wall wells are molded from virgin polypropylene selected for low DNA binding
- Black lettering for easy well identification



Ordering Information

Catalog # Description

PX1 PCR Plate Sealer and Seals

181-4000	PX1 PCR Plate Sealer , includes heat sealing instrument, 96-well/384-well plate support block, sealing frame, power cord
181-4030	Optically Clear Heat Seal , package of 100 seals
181-4035	Permanent Clear Heat Seal , package of 100 seals
181-4040	Pierceable Foil Heat Seal , package of 100 seals
181-4045	Peelable Foil Heat Seal , package of 100 seals

Hard-Shell Plates

HSP-9601	Hard-Shell Low-Profile 96-Well Skirted PCR Plates , clear well, white shell, 50
HSS-9601	Hard-Shell High-Profile 96-Well Semi-Skirted PCR Plates , clear well, clear shell, 25
HSL-9601	Hard-Shell Low-Profile 96-Well Semi-Skirted PCR Plates , clear well, clear shell, 25

HSR-9901	Hard-Shell 96-Well 480 PCR Plates , clear well, clear shell, bar-coded, 25
HSP-3801	Hard-Shell 384-Well Standard PCR Plates , clear well, clear shell, 50
HSR-4801	Hard-Shell 384-Well 480 PCR Plates , clear well, clear shell, bar-coded, 50

For additional plate options and instrument compatibility charts, please request bulletin 5496.

Cy is a trademark of GE Healthcare group companies. FAM is a trademark of Applera Corporation. Quasar is a trademark of Biosearch Technologies, Inc. Texas Red is a trademark of Invitrogen Corporation.

Practice of the polymerase chain reaction (PCR) may require a license.

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 www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000