

Microseal® 96-Well Skirted PCR Plates

Thin-Wall Polypropylene

Microseal 96-well thin-wall polypropylene PCR plates are ideal for high-throughput thermal cycling applications using reaction volumes of 5–125 µl. These PCR plates meet ANSI-SBS dimensional standards 1-2004 and 4-2004 for footprint and well position before and after cycling, and they can be used reliably with most robotic workstations, plate readers, and robotic arms. The automation-friendly design also features locator holes and flat vertical sidewalls for secure handling and easy bar coding (bar-coded plates are also available). The polypropylene construction confers very low protein binding, and raised rims around each well provide an excellent surface for tight sealing with a variety of sealing systems.

Microseal 96-Well PCR Plates 50 plates/pkg

MSP-9601	Natural	MSP-9611	Pink	MSP-9631	Blue
MSP-9651	White	MSP-9621	Yellow	MSP-9641	Green
MSP-9661	Black			MSP-9605	Natural, bar-coded, serially numbered in Code 39 symbology

Related Products

ADR-9611	PCR plate lifter (96-well)	Spring-arm device raises plates 5 mm above cyclers block for easier robotic or manual access
ADR-3700	PCR plate adaptor	Allows skirted 96-well plates to be used in ABI 3700 DNA analyzers

Vessel Comparison Chart

Product	PCR Volumes	Product Highlights
Microseal skirted 96- and 384-well PCR plates MSP-series	5–125 µl (96-well) 5–25 µl (384-well)	<ul style="list-style-type: none"> High-throughput, robot-friendly design in 1-component, thin-wall polypropylene 96-well format is compatible with DNA Engine Opticon® real-time PCR detection systems
Hard-Shell® skirted 96- and 384-well PCR plates HSP-series	5–125 µl (96-well) 5–30 µl (384-well)*	<ul style="list-style-type: none"> Unique 2-component design resists warping and shrinkage Rigid skirt/deck for lab automation; thin-wall polypropylene wells for optimal thermal cycling 96-well format is compatible with DNA Engine Opticon real-time PCR detection systems
Microseal semi-skirted 96-well PCR plates MSS-series	5–125 µl	<ul style="list-style-type: none"> Semi-skirt design adds stiffness and extra labeling surface Compatible with instruments limited to full-height vessels (e.g., ABI cyclers and sequencers)
Multiplate™ unskirted PCR plates, low-profile and full-height MLL-series MLP-series	5–125 µl	<ul style="list-style-type: none"> High-performance, versatile, inexpensive plates can easily be cut to desired size when running fewer than 96 samples Low-profile plates are more rigid; reduce potential for condensate formation Available in 96- and 48-well formats

* Reliable cycling of substantially smaller volumes has been demonstrated using 384-well Hard-Shell plates with Microseal 'P+' pads and high-pressure motorized lids. Please inquire for more information.



Recommended Sealing Options for Microseal 96-Well Skirted PCR Plates

Product	Product Highlights	Compatibility
Strip caps TCS-series	<ul style="list-style-type: none"> • Tight sealing for thermal cycling, storage, or transport • Choice of dome caps or ultraclear, flat optical caps • PCR volumes $\geq 5 \mu\text{l}$ 	96-well polypropylene PCR plates and 0.2 ml tubes
Microseal 'A' film MSA-5001 (50 sheets)	<ul style="list-style-type: none"> • Smooth application and release — no aerosol generation • Peel-away release liner for pristine sealing surface • PCR volumes $\geq 10 \mu\text{l}$ (96-well) or $\geq 5 \mu\text{l}$ (384-well) 	All PCR plates, tube strips
Microseal 'B' adhesive seals MSB-1001 (100 sheets)	<ul style="list-style-type: none"> • Clear, adhesive seal for thermal cycling or fluorescent assays • Plate storage or transport before or after cycling (ambient to -40°C) • PCR volumes $\geq 10 \mu\text{l}$ (96-well) or $\geq 5 \mu\text{l}$ (384-well) 	All polypropylene PCR plates
Microseal 'F' foil MSF-1001 (100 sheets)	<ul style="list-style-type: none"> • Foil can be pierced for sample recovery • Plate storage (ambient to -70°C) • PCR volumes $\geq 25 \mu\text{l}$ (96-well) or $\geq 5 \mu\text{l}$ (384-well) 	All PCR plates
96-well PCR plate sealing mats 223-9442 (5 mats)	<ul style="list-style-type: none"> • Reusable mat — ideal for sequencing • Smooth release — no aerosol generation • PCR volumes $\geq 10 \mu\text{l}$ 	96-well polypropylene PCR plates
Chill-out™ liquid wax, red or clear CHO-series	<ul style="list-style-type: none"> • Replaces mineral oil overlay — solidifies below 10°C for easy sample recovery • Protects samples from spills or aerosol formation during chilled transport • Clear Chill-out overlay is ideal for reading fluorescent assays • PCR volumes $\geq 2 \mu\text{l}$ 	All tubes, 96-well PCR plates, Hard-Shell 384-well PCR plates
Microseal 'P' and 'P+' Pads MSP-1001 MSP-1002 (10 pads)	<ul style="list-style-type: none"> • Reusable sealing pad adheres to the heated lid • Use MSP-1001 with Power Bonnet™ motorized lid (ALP-1296, ALP-1238), using PCR volumes $\geq 15 \mu\text{l}$ (96-well) or $\geq 10 \mu\text{l}$ (384-well) • Use MSP-1002 with Moto Alpha™ motorized lid (ALP-2296, ALP-2238), using PCR volumes $\geq 10 \mu\text{l}$ (96-well) or $\geq 5 \mu\text{l}$ (384-well) 	All 384-well and 96-well PCR plates with raised well rims
Auto-sealing lids MSL-2012 (flat) MSL-2022 (arched) MSL-2032 (arched, wide-tab) (4 lids)	<ul style="list-style-type: none"> • Reusable, robot-placeable sealer reduces evaporation during reaction assembly and seals tightly during thermal cycling • Flat lid can be used for postcycling storage; PCR volumes $\geq 3 \mu\text{l}$ (96-well) or $\geq 1 \mu\text{l}$ (384-well) • Arched lids can be removed by robotic arms; PCR volumes $\geq 5 \mu\text{l}$ (96-well) or $\geq 1 \mu\text{l}$ (384-well) 	Hard-Shell and Microseal skirted PCR plates

Instrument Compatibility

Microseal 96-well skirted PCR plates fit Bio-Rad 96-well Alpha™ units, including DNA Engine Opticon and Chromo4™ real-time systems. They also fit BaseStation® and AP Biotech MegaBACE DNA analyzers, and many other instruments. MSP 96-well plates fit the ABI 3700 DNA analyzer using the adaptor listed.



This product was process-sampled and has tested negative for RNase, DNase, and human DNA. Please contact Bio-Rad Laboratories, Inc. for details of the testing procedure.

Practice of the patented polymerase chain reaction (PCR) process requires a license. Bio-Rad and MJ brand thermal cyclers and systems include an Authorized Thermal Cycler and may be used with PCR licenses available from Applied Biosystems. Their use with Authorized Reagents also provides a limited PCR license in accordance with the label rights accompanying such reagents. Some applications may also require licenses from other third parties. ABI is a trademark of Applied Biosystems.

