

Microseal® Semi-Skirted PCR Plates

96-Well Thin-Wall Polypropylene

Microseal 96-well semi-skirted PCR plates combine the stiffness of a skirted plate with the high profile required for compatibility with some instruments. These V-bottom thermal cycling plates provide the same thin-wall polypropylene wells found in the skirted Microseal and unskirted Multiplate™ line of plates. The stiff half-height skirt can be gripped by automated plate handlers, and it is a useful surface for plate labeling. Semi-skirted plates are a convenient option for laboratories that use both Bio-Rad and ABI instruments because these 96-well plates can be used in ABI instruments without special adaptors.

Microseal semi-skirted plates are available in natural, pink, yellow, blue, and green tints. Raised rims around each well provide a superior surface for tight sealing with strip caps, sealing films and pads, and heat seals. Use a 3 mm adaptor plate (ADR-0103) to ensure proper lid contact when sealing with film or mats. For added stability while assembling reactions or centrifuging, place semi-skirted plates in a 96-place rack (TRC-0501).

Microseal Semi-Skirted 96-Well PCR Plates 8 x 12 well array, 25 plates/pkg

MSS-9601 Natural MSS-9611 Pink MSS-9631 Blue MSS-9651 White
 MSS-9621 Yellow MSS-9641 Green MSS-9661 Black

Related Products

ADR-0103 3 mm adaptor plate, required to transfer lid pressure when sealing semi-skirted plates with film or mats
 TRC-0501 96-place racks and covers, for preparing or centrifuging reactions, blue, 5 plates with covers
 TRC-0531 96-place racks and covers, for preparing or centrifuging reactions, assorted colors, 5 plates with covers

Vessel Comparison Chart

Product	PCR Volumes	Product Highlights
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
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 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



Recommended Sealing Options for Semi-Skirted 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 ultra-clear, 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
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

Note: Microseal 'B' clear adhesive seal is only recommended for use with semi-skirted plates if the end tabs are removed and a 3 mm adaptor is used (ADR-0103). Microseal 'F' foil seal is only recommended for sealing semi-skirted plates for storage, not for thermal cycling.

Instrument Compatibility

Microseal semi-skirted PCR plates fit Bio-Rad 96-well Alpha™ units that have manual or power lids, and also fit many other 96-well cyclers. These plates are compatible with instruments that require a full-height 96-well plate, including all thermal cyclers and sequencers made by Applied Biosystems. Microseal semi-skirted PCR plates do not fit in the DNA Engine Opticon and Opticon™ 2 real-time PCR detection systems. Please inquire for additional information.



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.

