

Microseal® 'F' Sealing Foil

for PCR Plates

This aluminized foil serves as an exceptional barrier against evaporation from -70°C to 105°C . It is thin enough to pierce with a pipet tip to recover samples from individual wells. Minimum recommended PCR sample volumes are $25\ \mu\text{l}$ in 96-well PCR plates and $5\ \mu\text{l}$ in 384-well plates.

Directions for Use

1. Remove white liner and place adhesive surface over wells.
2. Apply pressure with fingers or soft rubber roller to complete contact with each well.
3. Close the heated lid and apply pressure. When using MJ thermal cyclers with adjustable heated lids, turn the thumbwheel until the lid contacts the plate — when an increase in resistance is first felt. Continue to turn the thumbwheel about $1/2$ – $3/4$ turn past the point of first contact ($1/4$ of the thumbwheel circumference is visible, so two passes of the visible portion equals $1/2$ turn).
4. If evaporation occurs at the edges, tighten lid an additional $1/4$ turn of the thumbwheel.

Microseal 'F' Foil

MSF-1001 100 sheets per package

Related Product

MSR-0001 Sealing roller Quickly and uniformly presses seal over PCR plate wells

Sealer Comparison Chart

Product	Product Highlights	Compatibility
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
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
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
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



Microseal 'F' Foil Effectively Seals the Following Vessels

Product	PCR Volumes	Product Highlights
Multiplate™ unskirted PCR plates MLP-series	5–125 µl	<ul style="list-style-type: none"> • High-performance, versatile, and inexpensive • Full-height vessels can easily be cut to desired size when running fewer than 96 samples • Available in 96- and 48-well formats
Low-profile Multiplate unskirted PCR plates MLL-series	5–125 µl	<ul style="list-style-type: none"> • More rigid design, yet easily cut to desired size when running fewer than 96 samples • Low profile reduces potential for condensate formation • Available in 96- and 48-well formats
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)
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

* Reliable cycling of substantially smaller volumes has been demonstrated with 384-well Hard-Shell® plates using high-pressure motorized lids; please inquire for more 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

