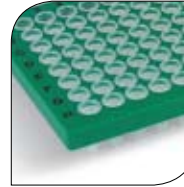


Amplification: PCR Plastic Consumables



PCR Plastic Consumables

Reaction Tubes, Plates, Sealers, and Accessories





PCR Plastic Consumables

One just right for you ...

Bio-Rad thin-wall PCR tubes, PCR plates, sealers, and accessories are precisely manufactured for optimal fit and cycling performance in Bio-Rad thermal cyclers and real-time systems. These high-quality consumables are suitable for a wide variety of applications. Bio-Rad tubes, tube caps, and PCR plates are molded, inspected, and packaged in a Class 100,000 or 10,000 cleanroom environment to prevent possible nucleic acid or nuclease contamination, then process sampled and tested to be negative for DNase, RNase, and human DNA.

When you open each box of Bio-Rad reaction vessels, you can be confident that the contents are in pristine condition. Contact your local sales representatives for sample packs.



The reaction vessels and sealing systems featured in this catalog have been designed to provide the best possible fit and performance with Bio-Rad and other brands of thermal cyclers and real-time PCR systems.



Thin-Wall PCR Tubes

Thin-wall polypropylene PCR tubes fit precisely in Bio-Rad sample blocks, ensuring accurate heat transfer to the sample. Extremely tight-sealing caps allow cycling of samples as small as 5 μ l in cyclers with a heated lid.

PCR Tube, Tube Strip, and Cap Selection Guide

Feature	0.5 ml Tubes	0.2 ml Tubes	High-Profile 0.2 ml Tube Strips	Low-Profile 0.2 ml Tube Strips	Domed Cap Strips	Optical Flat Cap Strips
Thermal cycling volume	10–200 μ l	5–125 μ l	5–125 μ l	5–125 μ l	\geq 5 μ l	\geq 5 μ l
Maximum volume	600 μ l	300 μ l	250 μ l	200 μ l		
Availability of white tubes				•		
Well profile	High	High	High	Low		

Individual PCR Tubes, 0.2 ml and 0.5 ml

These full-height (high-profile) PCR tubes have double-locking caps that won't pop open during cycling. PCR volume ranges are 5–125 μ l for 0.2 ml tubes and 10–200 μ l for 0.5 ml tubes. Tubes with flat, frosted caps for easy labeling are available in both 0.2 ml and 0.5 ml sizes (not suitable for real-time PCR).

To help prevent accidental contamination by multiple users, the 0.5 ml individual tubes with attached caps are available in resealable plastic bags of 100 tubes. Use of a capping tool is recommended for proper seating of caps on tubes.



Ordering Information

Catalog #	Description
Individual PCR Tubes With Attached Caps (0.2 ml)	
TFI-0201	PCR Tubes With Flat Caps (0.2 ml), natural, 1,000
TWI-0201	PCR Tubes With Domed Caps (0.2 ml), natural, 1,000
Individual PCR Tubes Without Caps (0.2 ml)	
TBI-0201	PCR Tubes Without Caps (0.2 ml), natural, 1,000
Individual PCR Tubes With Attached Caps (0.5 ml)	
TBI-0501	PCR Tubes With Flat Caps (0.5 ml), natural, 1,000 (2 bags of 500)
TBI-0502	PCR Tubes With Flat Caps (0.5 ml), natural, 800 (8 bags of 100)

Individual PCR Tubes With Attached Caps (0.2 ml)

TFI-0201	PCR Tubes With Flat Caps (0.2 ml), natural, 1,000
TWI-0201	PCR Tubes With Domed Caps (0.2 ml), natural, 1,000

Individual PCR Tubes Without Caps (0.2 ml)

TBI-0201	PCR Tubes Without Caps (0.2 ml), natural, 1,000
----------	---

Individual PCR Tubes With Attached Caps (0.5 ml)

TBI-0501	PCR Tubes With Flat Caps (0.5 ml), natural, 1,000 (2 bags of 500)
TBI-0502	PCR Tubes With Flat Caps (0.5 ml), natural, 800 (8 bags of 100)

PCR Tube and Cap Strips

Both tubes and caps are available in strips of 8 or 12, for use in 48-well and 96-well sample blocks.

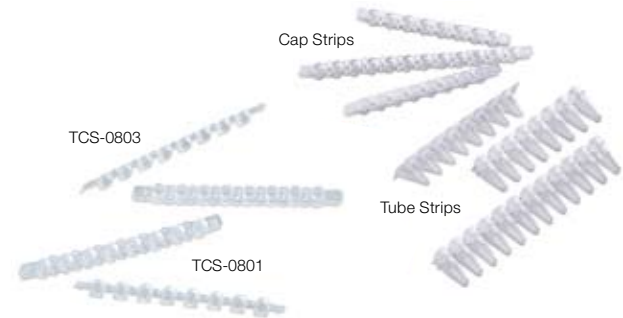
- Tight sealing and convenient handling for multiple samples
- Choice of domed or flat optical cap strips

Full-Height (High-Profile) 0.2 ml PCR Tube Strips

Recommended reaction volumes are 5–125 μ l. Tube strips and domed cap strips are also available packaged together in convenient bags sufficient for 96 samples. The resealable bags protect unused tubes and caps from accidental contamination.

Low-Profile 0.2 ml PCR Tube Strips

These tubes reduce the potential for condensate formation and allow greater light capture in fluorescence assays, such as those used in real-time PCR. Low-profile tubes are ideal for use in fast and low-volume PCR reactions. Overall height, including flat optical caps and 96-place rack, is 18.3 mm. Tube height is 15.5 mm. Low-profile tubes are available in opaque white for optical applications.



Cap Strips for 0.2 ml PCR Tubes and PCR Plates

These cap strips provide extremely tight sealing of all Bio-Rad 0.2 ml PCR tubes and plates during thermal cycling and cold storage. Flat cap strips feature ultraclear upper surfaces, ideal for fluorescence applications. Average light transmittance is 1.7-fold higher than with standard-clarity domed cap strips. Flat caps are available in strips of 8, and domed caps are available in strips of 8 or 12. Use of a capping tool is recommended for proper seating of caps on tubes or plates.

Ordering Information

Catalog #	Description
Full-Height Tube Strips Without Caps (0.2 ml)	
TBS-0201	8-Tube Strips Without Caps (0.2 ml) , 120
TBS-1201	12-Tube Strips Without Caps (0.2 ml) , natural, 100 strips (1,200 PCR tubes)
Low-Profile 8-Tube Strips Without Caps (0.2 ml)	
TLS-0801	Low-Profile 8-Tube Strips Without Caps (0.2 ml) , natural, 120 (960 PCR tubes)
TLS-0851	Low-Profile 8-Tube Strips Without Caps (0.2 ml) , white, 120 (960 PCR tubes)
Domed Cap Strips	
TCS-0801	Domed 8-Cap Strips , for 0.2 ml PCR tubes and plates, natural, 120
TCS-1201	Domed 12-Cap Strips , for 0.2 ml PCR tubes and plates, natural, 200
Optical Flat Cap Strips	
TCS-0803	Optical Flat 8-Cap Strips , for 0.2 ml PCR tubes and plates, ultraclear, 120
Full-Height Tube Strips With Domed Cap Strips (0.2 ml)	
TBC-0802	8-Tube Strips and Domed Cap Strips (0.2 ml) , natural, 20 bags of 12 x 8-tube strips and 12 x 8-cap strips (1,920 PCR tubes and 1,920 caps)
TBC-1202	12-Tube Strips and Domed Cap Strips (0.2 ml) , natural, 20 bags of 8 x 12-tube strips and 8 x 12-cap strips (1,920 PCR tubes and 1,920 caps)

Capping Tools and Racks

96-Place PCR Tube Rack and Cover

These stackable storage units for tubes and PCR plates provide a stable platform for preparing or centrifuging reactions.

Easy Cap™ Tool

The Easy Cap tool provides the pressure necessary to achieve a tight seal, one tube at a time, when capping individual or strip tubes. The narrow end securely fastens domed caps to 0.2 ml tubes or 96-well PCR plates. The wide end firmly holds thin-wall 0.5 ml tubes to prevent accidental crushing when opening or closing. The side slot allows easy opening of tight-fitting caps without generating aerosols.



Strip Cap Tool

This tool quickly and easily seats 8- and 12-cap strips on PCR plates or tubes. A grooved channel on one side is designed to seal domed caps, while the flat edge on the opposite side seats flat caps. For best results, seal tube strips while they are in the thermal cycler block or in a 96-place rack.

Ordering Information

Catalog #	Description
TRC-0501	96-Place Racks , with covers, for PCR tubes and unskirted and semi-skirted microplates, assorted colors, 5
ECT-1000	Easy Cap Tool , ensures tight seal for 0.2 ml PCR tubes or 96-well microplates
ECT-2000	Strip Cap Tool , for seating 8- and 12-cap strips on PCR plates or tubes

48-Well PCR Plates

Multiplate™ 48-Well PCR Plates

The versatile, unskirted design and 48-well format make these Multiplate unskirted PCR plates ideal for laboratories using 48-well blocks on Bio-Rad instruments. The plates are suitable for reaction volumes of 5–125 µl. Two plate styles are available:

- **Full-height (high-profile, 20.70 mm) wells, natural color** — designed to fit in most thermal cyclers
- **Low-profile (15.50 mm) wells, natural color or white** — optimized for fast PCR and low-volume reactions

Compatible Instruments

C1000™, S1000™, DNA Engine® family, MJ Mini™, MiniOpticon™ (low-profile white plates recommended)



Multiplate Full-Height 48-Well PCR Plate



Multiplate Low-Profile 48-Well PCR Plate

Ordering Information

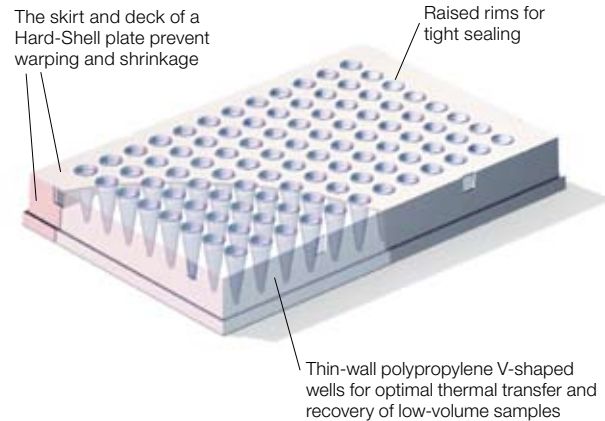
Catalog #	Description
MLP-4801	Multiplate 48-Well Unskirted PCR Plates , natural, 50
MLL-4801	Multiplate Low-Profile 48-Well Unskirted PCR Plates , natural, 50
MLL-4851	Multiplate Low-Profile 48-Well Unskirted PCR Plates , white, 50

96-Well PCR Plates

Bio-Rad offers 96-well PCR plates in several formats that are ideal for use with any PCR instrument. Unskirted, semi-skirted, and skirted plates are available for various PCR and real-time PCR needs.

Hard-Shell® 96-Well PCR Plates

Hard-Shell PCR plates are specifically designed to withstand the stresses of heat sealing, thermal cycling, and robotic handling. The patented two-component design features a skirt and deck molded from a rigid, thermostable polymer. In a separate step, the thin-wall wells are molded of virgin polypropylene selected for low DNA binding. This design prevents problems due to the warping, shrinkage, and sticking that may occur when single-component polypropylene PCR plates are exposed to the high temperatures of thermal cycling or heat sealing. Thus, performance is improved in many applications. In addition, these plates can withstand -80°C storage and high centrifugation forces, convenient for alcohol precipitations.



Hard-Shell 96-Well Skirted PCR Plates, Low-Profile

Features of the skirted plates include:

- Reaction volumes of 5–125 µl (200 µl maximum)
- Low-profile (16.05 mm) wells optimized for low-volume reactions and fast PCR
- Superior stability and flatness, allowing precise positioning for automation
- Full skirt for robotic handling and labeling surface
- Footprint and well spacing that match ANSI/SBS standard dimensions



Hard-Shell 96-Well Skirted PCR Plate

Compatible Instruments

C1000, S1000, DNA Engine family, PTC-100®, CFX96™, Chromo4™, Opticon™, Eppendorf Mastercycler series, Idaho Technology LightScanner

Ordering Information			
Description	Clear Wells	White Wells	Black Wells
Hard-Shell Thin-Wall 96-Well Skirted PCR Plates			
White Shell, 50	HSP-9601	HSP-9655	—
Red Shell, 50	HSP-9611	—	—
Yellow Shell, 50	HSP-9621	—	—
Blue Shell, 50	HSP-9631	HSP-9635	—
Green Shell, 50	HSP-9641	HSP-9645	—
Black Shell, 50	HSP-9661	HSP-9665	HSP-9666
White Shell, bar-coded, 50	HSP-9901	HSP-9955	—
White Shell, bulk pack of 400 plates	HSP-9601B	—	—

96-Well PCR Plates

www.bio-rad.com/pcrplastics/

Hard-Shell 96-Well Semi-Skirted PCR Plates, Full-Height (High-Profile)

This Hard-Shell PCR plate fits most thermal cyclers. Features of semi-skirted plates include:

- Reaction volumes of 5–125 μ l (350 μ l maximum)
- Full-height (20.75 mm) wells that fit most thermal cyclers, real-time PCR detection systems, and DNA sequencers
- Warp-free half-height skirt for improved robotic handling and labeling surface

Compatible Instruments

C1000, S1000, DNA Engine family, PTC-100, iCycler[®], iQ[™]5, iCycler iQ[®], MyiQ[™], MyiQ[™]2, Chromo4, Applied Biosystems 0.2 ml tube cyclers, standard real-time systems, and DNA sequencers (3130, 7300, 7500, 9700, Veriti, etc.), Eppendorf Mastercycler series, Stratagene (Agilent) Mx series, Idaho Technology LightScanner



Hard-Shell 96-Well Semi-Skirted PCR Plate

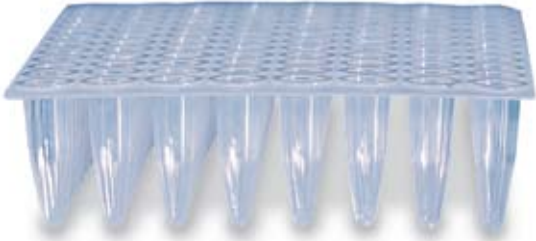
Ordering Information

Description	Clear Wells	White Wells	Black Wells
Hard-Shell Full-Height 96-Well Semi-Skirted PCR Plates, 25			
Clear Shell	HSS-9601	—	—
Green Shell	HSS-9641	—	—
Black Shell	—	HSS-9665	—
Clear Shell, bar-coded	HSS-9901	—	—

Multiplate™ 96-Well PCR Plates

Multiplate 96-Well Unskirted PCR Plates, Full-Height (High-Profile)

The single-component polypropylene construction of Multiplate PCR plates confers very low protein binding and excellent retention of sample. When less than a full plate is needed, these plates are easily cut with scissors to the required size. The plates are designed for compatibility with a wide variety of sealing methods for oil-free cycling, and are suitable for PCR volumes of 5–125 µl.



Multiplate Full-Height 96-Well PCR Plate

Compatible Instruments

C1000, S1000, DNA Engine family, PTC-100, MyCycler™, iCycler, iQ5, iCycler iQ, MyiQ, MyiQ2, Chromo4, Applied Biosystems 0.2 ml tube cyclers, standard real-time systems, and DNA sequencers, Eppendorf Mastercycler series, Stratagene (Agilent) Mx series

Ordering Information	
Catalog #	Description
Multiplate 96-Well Unskirted PCR Plates	
MLP-9601	Multiplate 96-Well Unskirted PCR Plates , natural, 25 plates
MLP-9651	Multiplate 96-Well Unskirted PCR Plates , white, 25 plates
MLP-9611	Multiplate 96-Well Unskirted PCR Plates , pink, 25 plates
MLP-9631	Multiplate 96-Well Unskirted PCR Plates , blue, 25 plates

96-Well PCR Plates

www.bio-rad.com/pcrplastics/

Multiplate 96-Well Unskirted PCR Plates, Low-Profile

Multiplate low-profile PCR plates combine the unskirted feature of the original Multiplate plate and the 5 mm lower overall height. The lower height (15.50 mm) reduces the potential for condensate formation and offers advantages for fast PCR, low-volume reactions, and light capture in fluorescence assays, such as those used in real-time PCR. A rigid top surface provides firm handling, while still allowing the plate to be cut for use in other formats.

Compatible Instruments

C1000, S1000, DNA Engine family, PTC-100, CFX96, Chromo4, Opticon, Applied Biosystems 0.1 ml tube cyclers and fast real-time systems, Eppendorf Mastercycler series, Idaho Technology LightScanner



Multiplate Low-Profile 96-Well PCR Plate

Ordering Information

Catalog #	Description
-----------	-------------

Multiplate Low-Profile 96-Well Unskirted PCR Plates

MLL-9601	Multiplate Low-Profile 96-Well Unskirted PCR Plates, natural, 25 plates
MLL-9651	Multiplate Low-Profile 96-Well Unskirted PCR Plates, white, 25 plates

iQ™ 96-Well Semi-Skirted Real-Time PCR Plates, Full-Height

These semi-skirted, full-height PCR plates are optimized for iQ5, iCycler iQ, MyiQ2, and MyiQ real-time PCR detection systems. The semi-skirted design adds stiffness and a labeling surface. Plates are perforated every three columns for easy setup of triplicate reactions.

Compatible Instruments

C1000, S1000, DNA Engine family, PTC-100, MyCycler, iCycler, MyiQ, iCycler iQ, iQ5, MyiQ2, Chromo4, Applied Biosystems 0.2 ml tube cyclers and standard real-time systems, Eppendorf Mastercycler series, Stratagene (Agilent) Mx series



Ordering Information

Catalog #	Description
-----------	-------------

iQ 96-Well PCR Plates

223-9441	iQ 96-Well PCR Plates, 25
----------	---------------------------

Microseal® 96-Well PCR Plates

Microseal 96-Well Semi-Skirted PCR Plates, Full-Height (High-Profile)

These PCR plates are designed for Applied Biosystems 0.2 ml tube cyclers, real-time PCR systems, and DNA sequencers. The semi-skirted design adds stiffness and a surface for labeling. They are not recommended for use in Bio-Rad cyclers because the raised ridges around the plate prevent proper sealing in these instruments.

Compatible Instruments

Applied Biosystems 0.2 ml tube cyclers, standard real-time systems, and DNA sequencers

Microseal 96-Well Skirted PCR Plates, Low-Profile

These PCR plates feature single-component construction and a skirted design that is suitable for high-throughput plate handling. The robot-friendly design features low-binding polypropylene construction, locator holes, and flat vertical sidewalls for secure handling and easy bar coding. Raised rims around wells provide an excellent surface for tight sealing with a variety of sealing methods, and allow easy release of the sealer from the plate. Bar-coded plates are also available.



Microseal Full-Height 96-Well Semi-Skirted PCR Plate



Microseal Low-Profile 96-Well Skirted PCR Plate

Compatible Instruments

C1000, S1000, DNA Engine family, PTC-100, CFX96, Chromo4, Opticon, Eppendorf Mastercycler series

Ordering Information

Catalog # Description

Microseal 96-Well Semi-Skirted PCR Plates, Full-Height

MSS-9601 Microseal 96-Well Semi-Skirted PCR Plates, natural, 25

Microseal 96-Well Skirted PCR Plates, Low-Profile

MSP-9601 Microseal 96-Well Skirted PCR Plates, natural, 50

MSP-9605 Microseal 96-Well Skirted PCR Plates, bar-coded, natural, 50

Concord™ 96-Well Skirted Polycarbonate PCR Plates, Low-Profile

Concord 96-well PCR plates have a thin-wall polycarbonate construction that allows excellent heat transfer. A minimum sample volume of 20 µl is recommended for oil-free cycling with these plates.



Ordering Information

Catalog # Description

CON-9601* Concord 96-Well Polycarbonate PCR Plates, low-profile, skirted, natural, 25

CVR-9601 Dust Covers, for Concord polycarbonate PCR wells, nonsealing, 25

* Not recommended for use with ³⁵S.

384-Well PCR Plates

Hard-Shell® 384-Well Skirted PCR Plates

Hard-Shell PCR plates are designed to withstand the stresses of thermal cycling and robotic handling. The patented two-component design provides superior stability and flatness, allowing precise positioning for automation. Features include:

- Reaction volumes of 1–30 µl (50 µl maximum)
- Extremely uniform wells that reduce well-to-well variability in optical assays, such as those used in real-time PCR
- Compatibility with most 384-well thermal cyclers, real-time PCR detection systems, and DNA sequencers
- White-well option for increased fluorescent signal strength
- Color-coded skirts with clear or white wells
- Warp-free skirt and deck for improved robotic handling
- Footprint and well spacing that match ANSI/SBS standard dimensions
- Low-cost, user-readable bar code option for convenient database tracking



Compatible Instruments

C1000, S1000, DNA Engine family, CFX384™, Applied Biosystems cyclers, real-time systems, and DNA sequencers, Eppendorf Mastercycler series, Idaho Technology LightScanner

Ordering Information

Description	Clear Wells	White Wells	Black Wells
Hard-Shell Thin-Wall 384-Well Skirted PCR Plates			
Clear Shell, 50	HSP-3801	HSP-3805	—
White Shell, 50	HSP-3851	—	—
Red Shell, 50	HSP-3811	—	—
Yellow Shell, 50	HSP-3821	—	—
Blue Shell, 50	HSP-3831	—	—
Green Shell, 50	HSP-3841	—	—
Black Shell, 50	—	HSP-3865	HSP-3866
Clear Shell, bar-coded, 50	HSP-3901	HSP-3905	—
Clear Shell, bulk pack of 500 plates	HSP-3801B	—	—

384-Well PCR Plates

www.bio-rad.com/pcrplastics/

Microseal® 384-Well Skirted PCR Plates

Microseal 384-well skirted PCR plates feature single-component construction and a skirted design that is suitable for high-throughput plate handling. The plates are available in clear, opaque white, and black for luminescence- and fluorescence-based assays. Bar-coded plates are also available.

Compatible Instruments

C1000, S1000, DNA Engine family, iCycler, CFX384, Applied Biosystems cyclers, real-time systems, and DNA sequencers, Eppendorf Mastercycler series, Idaho Technology LightScanner

Ordering Information

Description	50 Plates	Bar-Coded, 50 Plates
Microseal 384-Well Skirted PCR Plates		
Natural	MSP-3842	MSP-3846
White	MSP-3852	—
Black	MSP-3862	—

PCR Plate Accessories

Microseal® 384 Plate Positioner

To achieve consistent results in automated processes such as dispensing and retrieving small volumes (<5 µl) in a PCR plate, precise positioning and uniform well depth are required. Placing a Microseal 384-well PCR plate onto the Microseal 384 plate positioner secures the PCR plate and flattens it if needed. The positioner footprint (127.90 x 85.85 mm) is designed to fit securely into the plate nest of commonly used automated liquid handlers.

Heat Sealer Support Plate

This thin aluminum support plate ensures optimal sealing of Hard-Shell 384-well PCR plates when using automated heat sealers, so less fine-tuning of heat seal conditions is needed.










Microseal 384 Plate Positioner










Ordering Information	
Catalog #	Description
ADR-3841	Microseal 384 Plate Positioner , for automated liquid handling
ADR-3001	Heat Sealer Support Plate , for supporting Hard-Shell 384-well plates in automated heat sealers

Selection Guide

www.bio-rad.com/pcrplastics/

Instrument Compatibility of PCR Plastic Consumables








	0.2 ml Tubes			48- and 96-			
	Individual Full-Height	Strips Full-Height	Strips Low-Profile	Hard-Shell Semi-Skirted Full-Height	Hard-Shell Skirted Low-Profile	Multiplate Unskirted Full-Height	Multiplate Unskirted Low-Profile
							
Catalog #	pages 4, 5	pages 6, 7	pages 6, 7	pages 10, 12	pages 10, 11	page 13	page 14
Thermal Cycler							
Bio-Rad C1000 and S1000	•	•	•	•	•	•	•
Bio-Rad DNA Engine line, PTC-100	•	•	•	•	•	•	•
Bio-Rad MyCycler	•	•				•	
Bio-Rad iCycler	•	•		•		•	
Bio-Rad MJ Mini	•	•	•			•	•
Applied Biosystems 0.2 ml tube cyclers (2720, 9700, Veriti)	•	•		•		•	
Applied Biosystems 0.1 ml tube cyclers (9800 fast, Veriti fast)			•				•
Applied Biosystems 384-well cyclers (9700, Veriti)							
Eppendorf Mastercycler series	•	•	•	•	•	•	•

Well Plates				384-Well Plates		Sealers and Caps		
iQ Semi-Skirted Full-Height	Concord Skirted Low-Profile	Microseal Semi-Skirted Full-Height	Microseal Skirted Low-Profile	Hard-Shell 384	Microseal 384	Microseal 'B' Seals	Optical Cap Strips	Domed Cap Strips
								
page 15	page 17	page 16	page 16	pages 18, 19	page 20	pages 28–30	pages 6, 7	pages 6, 7
•	•		•	•	•	•	•	•
•	•		•	•	•	•	•	•
•						•	•	•
•						•	•	•
•		•				•	•	•
						•	•	•
				•	•	•		
•	•		•	•	•	•	•	•









Selection Guide

www.bio-rad.com/pcrplastics/

Instrument Compatibility of PCR Plastic Consumables (cont.)

	0.2 ml Tubes						48- and 96-
	Individual Full-Height	Strips Full-Height	Strips Low-Profile	Hard-Shell Semi-Skirted Full-Height	Hard-Shell Skirted Low-Profile	Multiplate Unskirted Full-Height	Multiplate Unskirted Low-Profile
							
Catalog #	pages 4, 5	pages 6, 7	pages 6, 7	pages 10, 12	pages 10, 11	page 13	page 14
Real-Time PCR System							
Bio-Rad CFX96 or CFX384*			•		•		•
Bio-Rad MyiQ2, iQ5, iCycler iQ, MyiQ		•		•		•	
Bio-Rad Chromo4		•	•	•	•	•	•
Bio-Rad Opticon, Opticon 2			•		•		•
Bio-Rad MiniOpticon*			•				•
Applied Biosystems standard systems (7300, 7500, 7900HT)		•		•		• (except 7900HT)	
Applied Biosystems fast systems (7500 fast, StepOne, StepOnePlus)			•				•
Eppendorf Mastercycler ep <i>realplex</i>		•	•	•	•	•	•
Stratagene (Agilent) Mx series		•		•		•	
Corbett (QIAGEN) Rotor-Gene	•						
Other Instruments							
Applied Biosystems DNA sequencers (3100, 3700, 3730)				•		•	
Idaho Technology LightScanner				•	•	•	•

* CFX384 and MiniOpticon real-time PCR detection systems are factory calibrated for white tubes and white-well plates. White plastics are recommended due to their superior performance.

Well Plates				384-Well Plates		Sealers and Caps	
iQ Semi-Skirted Full-Height	Concord Skirted Low-Profile	Microseal Semi-Skirted Full-Height	Microseal Skirted Low-Profile	Hard-Shell 384	Microseal 384	Microseal 'B' Seals	Optical Cap Strips
							
page 15	page 17	page 16	page 16	pages 18, 19	page 20	pages 28–30	pages 6, 7
			•	•	•	•	•
•						•	•
•			•			•	•
			•			•	•
• (except 7900HT)		•		•	•	•	•
						•	•
•			•			•	•
•						•	•
		•		•	•		
			•	•	•		

or signal-to-noise ratio. Using clear tubes or clear-well plates on these instruments will require user calibration.

Selection Guide

www.bio-rad.com/pcrplastics/

Comparison of PCR Plates*

Feature	Multiplate 48- or 96-Well Unskirted**		Hard-Shell 96-Well	Hard-Shell Skirted		iQ 96-Well	Microseal			Concord Polycarbonate 96-Well
	Full-Height	Low-Profile	Semi-Skirted	96-Well	384-Well	Semi-Skirted	96-Well	96-Well	384-Well	96-Well
Thermal cycling volume	5–125 µl	5–125 µl	5–125 µl	5–125 µl	1–30 µl	5–125 µl	5–125 µl	5–125 µl	5–25 µl	20–150 µl
Maximum well volume	330 µl	200 µl	330 µl	250 µl	50 µl	330 µl	330 µl	220 µl	40 µl	210 µl
Automation/ bar coding			•	•	•			•	•	
Rigid, warp-free skirt and deck			•	•	•					
Availability of white or black wells	•	•	•	•	•				•	
Color-coded skirts with clear wells			•	•	•					
Well profile	High	Low	High	Low		High	High	Low		Low

* See sealer selection chart, page 27.

** Multiplate unskirted plates can be cut to the size needed when running small numbers of reactions.

Sealers

Effective sealing is essential for PCR and qPCR reactions. Besides cap strips (pages 6, 7), Bio-Rad offers many sealing options to best fit your needs.

Sealer Selection Guide

Feature or Use	Microseal				96-Well	Microseal		Auto-Sealing	Chill-out™
	'A' Film	'B' Seal	Sealing Kit	'F' Foil	Sealing Mats	'P' Pads*	'P+' Pads*	Lids	Liquid Wax
	pages 28–30				pages 29, 30	pages 31, 32		pages 31, 32	page 33
Thermal cycling volume									
96-well	≥10 µl	≥10 µl	≥5 µl	≥25 µl		≥15 µl	≥5 µl	≥3 µl	≥2 µl
384-well	≥5 µl	≥5 µl		≥5 µl		≥10 µl	≥4 µl	≥1 µl	≥2 µl
Recommended for real-time systems		•	•						•
Automated sealing						•	•	•	•
Reusable					•	•	•	•	
Cold storage		•	•	•				•	•
Pierceable seal				•					•
Adhesive-based seal		•	•	•					
Sealing 96-well plates	•	•	•	•	•	•	•	•	•
Sealing 384-well plates	•	•		•		•	•	•	•
Sealing Concord plates	•			•					•
Suggested for cyclers without heated lids									•

* Reaction volume data based on use with a motorized, high-pressure lid.

PCR Plate Sealers

Microseal® 'A' Film

Microseal 'A' film quickly and effectively seals the full range of Bio-Rad PCR plates and tubes. The pliant inner layer is designed to seal tightly during cycling, yet release smoothly to minimize the risk of aerosol formation and cross-contamination of samples. This film is easily cut for use with fewer than 96 wells; a peel-away release liner protects the sealing surface from contamination (not suitable for qPCR). An optional sealing roller provides a quick way to firmly seat Microseal 'A' film on an entire array of wells.



Microseal 'A' Film

Sealing Roller

Microseal® 'B' Adhesive Seals, Optically Clear

Microseal 'B' seals provide an adhesive-based sealing option for thermal cycling using thin-wall PCR plates. The strong adhesive layer ensures secure sample storage (as low as -20°C and up to 120°C) or transport before or after cycling, as well as tight sealing during thermal cycling when supplemental pressure is applied by a heated lid. This clear polyester film allows easy inspection of sample wells and effective light transmission for optical assays. The sealing surface is protected from contamination by a peel-away release liner. Perforated end-tabs allow removal of overhanging film for automation and other applications.



Microseal 'B' Adhesive Seals

Microseal 'B' adhesive seal accessories include:

- **Optical compression pad (96-well)** — enhances the seal integrity of Microseal 'B' clear seals when used in real-time PCR detection systems
- **Optical film sealing kit** — contains 100 Microseal 'B' clear seals and an optical compression pad

Microseal® 'F' Foil

These aluminized foil seals act as a barrier against evaporation from -80°C to 150°C . They are thin enough to pierce with a pipet tip for recovery of samples from individual wells, and they are suitable for use with automated systems, such as the ABI 3730 DNA analyzer. In addition to cold storage applications, they can be used for thermal cycling for sample volumes of $25\ \mu\text{l}$ (96-well) or $5\ \mu\text{l}$ (384-well).

96-Well PCR Plate Sealing Mats

These reusable mats are convenient for sealing 96-well PCR plates; they are not suitable for qPCR.

Pressure Pad

This foam pad with magnet distributes lid pressure uniformly on sealing film on plates used in thermal cyclers.



Optical Compression Pad



Microseal 'F' Foil



96-Well PCR Plate Sealing Mat

www.bio-rad.com/pcrplastics/

Optical Sealing Tape

An adhesive-based sealing option for optical assays demanding a low-fluorescent background. Recommended for the Idaho Technology LightScanner system.

Ordering Information

Catalog #	Description
MSA-5001	Microseal 'A' Film , package of 50 seals
MSR-0001	Sealing Roller , for film seals
MSB-1001	Microseal 'B' Adhesive Seals , optically clear, 100
ADR-3296	Optical Compression Pad , for improved film sealing of 96-well plates in DNA Engine Opticon 2 and Chromo4 systems
MSO-1001	Optical Film Sealing Kit , for 96-well plates, includes optical compression pad, 100 Microseal 'B' clear adhesive seals
MSF-1001	Microseal 'F' Foil , package of 100 seals
223-9442	96-Well PCR Plate Sealing Mats , 5
ADR-5001	Pressure Pad , uniformly distributes lid pressure for sealing film
223-9444	Optical Sealing Tape , package of 100 sheets

Sealing Pads for Automation

Microseal® 'P' and 'P+' Sealing Pads

These reusable sealing pads are designed to adhere to a motorized heated lid. They are ideal for applications such as cycle sequencing, in which several successive runs may be sealed with the same pad. Use 'P' pads with Power Bonnet™ lids, and 'P+' pads with Moto Alpha™ units. Microseal 'P+' pads provide improved sealing of low-volume reactions — as low as 5 µl in 96-well plates and 1 µl in 384-well plates. Each pad may be used for approximately 25 runs. Microseal 'P' and 'P+' pads are intended for use exclusively with the higher sealing pressures provided by a motorized lid.

Auto-Sealing Lids for PCR Plates

These lids are reusable automation-friendly sealers that prevent evaporation and contamination during reaction assembly and seal tightly for thermal cycling when the cycler lid is closed. They can be reused up to 50 times. The lids are constructed of metal with an attached compressible pad, and they can be applied by robotic gripper devices.



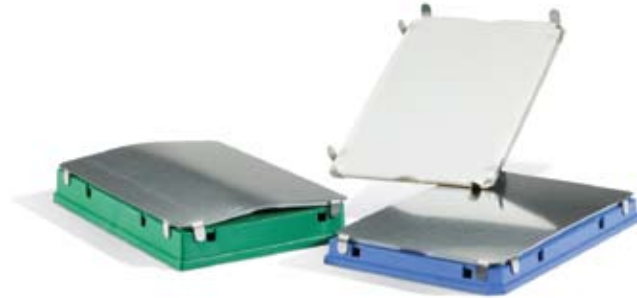
Microseal 'P+' Sealing Pad

Sealers

www.bio-rad.com/pcrplastics/

Three varieties of auto-sealing lids are available:

- **Flat lid** — allows low-volume cycling (down to 1 μ l) in 384-well plates using a Moto Alpha unit
- **Arched lid** — self-releasing; can also be removed by robotic devices
- **Arched lid with wide tabs** — especially well suited for gripping with robotic grippers



Auto-Sealing Lids

Ordering Information

Catalog #	Description
-----------	-------------

Microseal 'P' and 'P+' Pads

MSP-1001	Microseal 'P' Pads , reusable, for Power Bonnet lids, 10
MSP-1002	Microseal 'P+' Pads , reusable, for Moto Alpha unit lids, 10
MSP-1003	Microseal 'P' Replacement Pads , for MSL-2032, reusable, 10

Auto-Sealing Lids for PCR Plates

MSL-2012	Flat Auto-Sealing PCR Plate Lids , reusable, 4
MSL-2022	Arched Auto-Sealing PCR Plate Lids , reusable, 4
MSL-2032	Arched Auto-Sealing PCR Plate Lids With Wide Tabs , reusable, 4

Chill-out™ Liquid Wax

Chill-out liquid wax provides an excellent vapor barrier that may be used instead of mineral oil in thermal cyclers without heated lids. After cycling, the tubes are chilled below 10°C to harden the wax. The solid layer protects samples from spills or aerosol formation, but is easy to penetrate with a pipet tip for sample retrieval. Chill-out liquid wax is available in a clear formulation for use in fluorescence assays, such as those used in real-time PCR, and a bright red formulation, easily visible when recovering reaction products (not recommended for sealing Microseal 384-well PCR plates).

**Specifications**

Reaction Volume	Chill-out Wax Volume	Reaction Volume	Chill-out Wax Volume
2–20 μ l	10 μ l	60–80 μ l	25 μ l
20–40 μ l	15 μ l	>80 μ l	30 μ l
40–60 μ l	20 μ l		

Ordering Information

Catalog #	Description
CHO-1401	Chill-out Liquid Wax , red, 100 ml
CHO-1404	Chill-out Liquid Wax , red, 1 L
CHO-1411	Chill-out Liquid Wax , clear, optical grade, 100 ml
CHO-1414	Chill-out Liquid Wax , clear, optical grade, 1 L

Frame-Seal™ Incubation Chambers

Frame-Seal incubation chambers are easy-to-use, strongly adhesive hybridization chambers with flexible plastic coverslips. They provide vapor-tight sealing for FISH, colonies, in situ PCR, and PRINS and allow samples to be recovered easily. The seal withstands temperatures up to 97°C. Frame-Seal chambers should be used with plain (unprinted) glass slides. Slides with highly hydrophobic ink patterns (such as Teflon) are not recommended for use with Frame-Seal chambers.

- UV-treatable for inactivation of contaminating DNA
- Compatible with Bio-Rad and other horizontal-format thermal cyclers that accommodate slides



Ordering Information

Catalog #	Description
SLF-0201	Frame-Seal Incubation Chambers , 9 x 9 mm, 25 µl capacity, coverslips included, 100
SLF-0601	Frame-Seal Incubation Chambers , 15 x 15 mm, 65 µl capacity
SLF-1201	Frame-Seal Incubation Chambers , 17 x 28 mm, 125 µl capacity
SLF-3001	Frame-Seal Incubation Chambers , 19 x 60 mm, 300 µl capacity

Liquid Handling Solutions

For liquid-handling applications, Bio-Rad offers pipet tips, reagent reservoirs, and test tubes.



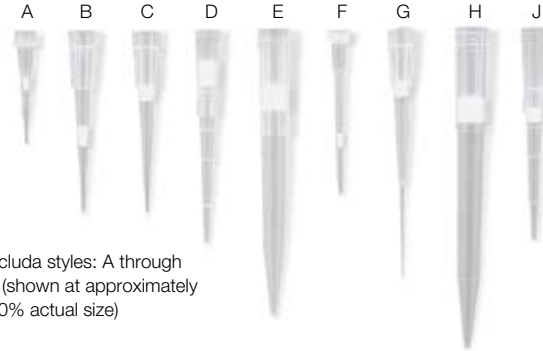
Pipet Tips

Bio-Rad's pipet tips are made from virgin polypropylene and have been accurately molded for an airtight fit. The finest molds and manufacturing methods are used, so the tips are guaranteed to have a smooth interior surface, which is essential for precision pipetting.

- Tips and racks are autoclavable at 120°C at 15 lb of pressure for 15 minutes; tips presterilized by E-beam irradiation are also available
- All materials used in both natural and colored pipet tips have been formulated without heavy metals
- Xcluda™ pipet tips have been independently tested and certified to be free of DNase, RNase, and pyrogens
- Pipet tip racks are manufactured from polypropylene, with a plastic content code of 5, and are free of labels for convenient recycling

Xcluda™ Aerosol Barrier Pipet Tips

Xcluda aerosol barrier pipet tips guard against aerosol contamination of samples, a feature of particular importance in PCR experiments. The barrier will not seal on contact with liquid if unintentional overpipetting occurs, which protects samples from accidental loss. Offered in nine different styles, the tips fit a variety of pipets. They are available presterilized in fully enclosed racks and are independently tested and certified to be free of DNase, RNase, and pyrogens.



Xcluda styles: A through J (shown at approximately 60% actual size)

Ordering Information

Catalog #	Description		
211-2001	Xcluda Style A , 0.5–10 µl, 960	211-2026	Xcluda Style F , 0.5–10 µl, 1,000
211-2006	Xcluda Style B , 2–20 µl, 960	211-2031	Xcluda Style G , 10–100 µl, 1,000
211-2011	Xcluda Style C , 10–100 µl, 960	211-2036	Xcluda Style H , 100–1,000 µl, 1,000
211-2016	Xcluda Style D , 20–200 µl, 960	211-2041	Xcluda Style J , 10–100 µl, 960
211-2021	Xcluda Style E , 100–1,000 µl, 960		

Pipet Tips for Gel Loading

Seque/Pro™ Capillary Pipet Tips

Seque/Pro capillary pipet tips have an average OD of <0.3 mm, making them ideal for loading sequencing gels or IEF tube gels. Seque/Pro tips perform best when used with Eppendorf ultra micro pipets (0.5–10 µl), and can also be used with Rainin Pipetman 10 and 20 µl pipets.



Prot/Elec™ Pipet Tips

Prot/Elec tips fit easily within a gap of 0.75 mm between vertical slab gel plates while maintaining a large bore for fast sample flow. The 200 µl capacity tips are molded to fit Rainin Pipetman 20, 100, and 200 pipets and Eppendorf 20 and 100 µl pipets. Prot/Elec tips are also available with an aerosol barrier to guard against airborne contaminants and cross-contamination of samples (see Xcluda style G).



Ordering Information

Catalog #	Description
223-9911	Seque/Pro Capillary Pipet Tips , in enclosed autoclavable rack, 0.5–10 µl, 200
223-9912	Seque/Pro Capillary Pipet Tips , sterilized in enclosed rack, 0.5–10 µl, 200
223-9915	Prot/Elec Pipet Tips , bulk pack, plastic bag in dust-free box, 1–200 µl, 1,000
223-9917	Prot/Elec Pipet Tips , racked, 12 x 17 format, 204 per rack with a cover on each rack, 1–200 µl, 1,020
223-9916	Prot/Elec Pipet Tips , racked, 8 x 12 format, 96 per rack with a cover on each rack, 1–200 µl, 960

Tip Selection Guide

Pipet Type	Tip Type
Costar (8-Pette, 12-Pette)	
20–200 µl	28
25–200 µl	28
Eppendorf	
0.5–10 µl	Seque/Pro, Xcluda F, 31
2–20 µl	Prot/Elec, Xcluda B, 14, 35, 37–39
10–100 µl	Prot/Elec, Xcluda C and G, 35, 37–39
50–250 µl	39, 40, 41
100–1,000 µl	Xcluda H, 41
Excalibur	
1–200 µl	Prot/Elec, 35, 37–39
200–1,000 µl	40, 41
Thermo Labsystems Finnpiquette	
5–40 µl	Prot/Elec, Xcluda B, C, G, 35, 37–39
40–200 µl	Prot/Elec, Xcluda C and G, 35, 37–39
200–1,000 µl	Xcluda E
Oxford Benchmate	
0.5–10 µl	Seque/Pro, Xcluda A, 14
10–50 µl	Xcluda J, 35, 37–39
40–200 µl	Prot/Elec, Xcluda C, D, G, 35, 37–39
200–1,000 µl	Xcluda E

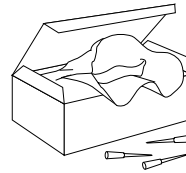
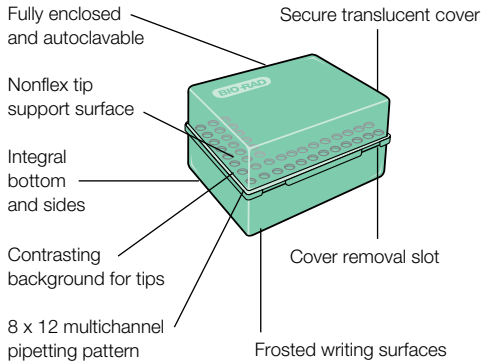
Pipet Type	Tip Type
Rainin Pipetman and EDP Series	
0.1–10 µl	Seque/Pro, Xcluda F, 31
2–20 µl	Prot/Elec, Seque/Pro, Xcluda B, 31, 35, 37–39
10–100 µl	Prot/Elec, Xcluda C and G, 35, 37–39
20–200 µl	Prot/Elec, Xcluda D, 35, 37–39
100–1,000 µl	Xcluda E, 40, 41
Socorex	
0.5–10 µl	Seque/Pro
1–200 µl	Prot/Elec, 35, 37–39
200–1,000 µl	40, 41
Titertek Flow	
5–200 µl	26
5–300 µl	Xcluda D
Volac	
1–20 µl	Prot/Elec, 31, 35, 37–39
1–200 µl	Prot/Elec, 35, 37–39
200–1,000 µl	40, 41

Standard Pipet Tips

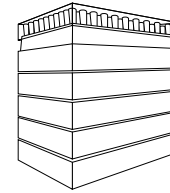


MTP

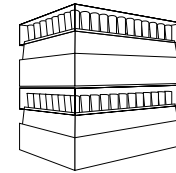
8 x 12 format, enclosed, nonflex racks are suitable for repeated reloading and autoclaving. Racked tips are also available presterilized.



BR
Bulk tips, plastic bag in a dust-free box.



RBR
Racked tips, one cover on a stack of nested racks.



TBR
Ideal for autoclaving, each rack has its own lid and covered bottom for full enclosure.

Tip Packaging Options

Packaging Option	Standard Tip Type Number									
	14	26	28	31	35	37	38	39	40	41
BR	•		•	•	•	•	•	•	•	•
RBR					•					
TBR	•				•				•	•
MTP		•			•	•	•			
MTP-S (sterilized)			•		•	•	•			

Ordering Information

Catalog # Description

MTP 8 x 12 Format, Enclosed Racks

223-9301 **MTP-26 Tips**, natural, 5–200 µl, 960
 223-9303 **MTP-35 Tips**, yellow, 1–200 µl, 960
 223-9304 **MTP-37 Tips**, natural, 1–200 µl, 960
 223-9313 **MTP-38 Tips**, yellow, graduated, beveled, 1–200 µl, 960

MTP-S 8 x 12 Format, Enclosed Racks, Presterilized*

223-9307 **MTP-28-S Tips**, natural, 25–200 µl, 960
 223-9308 **MTP-35-S Tips**, yellow, 1–200 µl, 960
 223-9309 **MTP-37-S Tips**, natural, 1–200 µl, 960
 223-9318 **MTP-38-S Tips**, yellow, graduated, beveled, 1–200 µl, 960

continues

Ordering Information

Catalog #	Description
BR Bulk Tips	
223-9014	BR-14 Tips , natural, 0.1–10 µl, 1,000
223-9028	BR-28 Tips , natural, 25–200 µl, 1,000
223-9031	BR-31 Tips , natural, 0.5–10 µl, 1,000
223-9035	BR-35 Tips , yellow, 1–200 µl, 1,000
223-9037	BR-37 Tips , natural, 1–200 µl, 1,000
223-9038	BR-38 Tips , yellow, graduated, beveled, 1–200 µl, 1,000
223-9039	BR-39 Tips , natural, graduated, beveled, 1–200 µl, 1,000
223-9040	BR-40 Tips , blue, 100–1,000 µl, 500
223-9041	BR-41 Tips , natural, 100–1,000 µl, 500
RBR Racked Tips, One Cover on Nested Racks	
223-9135	RBR-35 Tips , yellow, 1–200 µl, 1,000
TBR Racked Tips, Cover on Each Rack	
223-9354	TBR-14 Tips , natural, 0.1–10 µl, 1,000
223-9347	TBR-35 Tips , yellow, 1–200 µl, 1,000
223-9350	TBR-40 Tips , blue, 100–1,000 µl, 1,000
223-9351	TBR-41 Tips , natural, 100–1,000 µl, 1,000

BR Bulk Tips**RBR Racked Tips, One Cover on Nested Racks****TBR Racked Tips, Cover on Each Rack**

* Presterilized racks and tips are not suitable for autoclaving. Order nonsterile racks for reloading tips and autoclaving.

Reservoirs

www.bio-rad.com/liquidhandling/

Reagent Reservoirs

Disposable reagent reservoirs are ideal for dispensing reagent with multichannel pipets. Their sloping design makes it easy to pick up liquid to the last drop. The 50 ml capacity is graduated to 25 ml in 5 ml increments. These presterilized reagent reservoirs are not autoclavable.



Reagent Reservoir

Ordering Information

Catalog #	Description
224-4872	Sterilized Reagent Reservoirs , graduated, polystyrene, 5 per package, box of 200

Titertube® Micro Test Tubes

Titertube 1 ml micro test tubes allow efficient preparation and storage of large numbers of samples in an 8 x 12 format.

Tubes and racks are made of polypropylene; they are autoclavable to 120°C and freezable to –80°C. The polyethylene plugs, which seal tightly for storage, are not designed for repeated use and cannot be autoclaved. The tubes are offered in sterile and unsterilized strips of eight, which may be cut apart to separate the tubes. Titertube micro test tubes have:

- A 1 ml capacity, 8.8 x 45 mm
- Spacing that matches 96-well plates
- Enclosed racks that stack to store 5,800 samples/ft³
- Racks with a grid for identification of contents



Ordering Information

Catalog #	Description
223-9390	Titertube Micro Test Tubes , unsterilized, 10 racks of 96 tubes
223-9395	Titertube Micro Test Tubes , presterilized,* 10 racks of 96 tubes
223-9391	Titertube Micro Test Tubes , unsterilized, bulk, 1,000
223-9393	Titertube Plugs , unsterilized, 960 in 120 strips of 8
223-9392	Titertube Plugs , presterilized,* 960 in 120 strips of 8
223-9394	Titertube Rack , empty, holds 96 Titertube micro test tubes, 10

* Presterilized racks and polyethylene plugs are not suitable for autoclaving; order unsterilized racks to reload tubes and autoclave.

EZ Micro™ and Standard Micro Test Tubes

EZ Micro graduated polypropylene test tubes have a frosted marking area and a flat top that is pierceable by a 19 gauge or thicker needle. Standard polypropylene micro test tubes are recommended for solvent extraction and heating, when a very tight cap seal is required. Because of their tighter cap fit, standard tubes may be more difficult to manipulate in applications that require repeated opening and closing of the tubes. Separate polyethylene caps for the 1.5 ml capless tubes have a knurled top for easy handling.

The micro test tubes are:

- Suitable for general-purpose benchtop centrifuge use; sturdy uniform walls easily withstand up to 13,000 x g
- Autoclavable to 120°C; freezable to -80°C
- Available in 2 ml, 1.5 ml, and 500 µl sizes



Ordering Information

Catalog #	Description
223-9503	EZ Micro Test Tubes , 500 μ l, natural, 1,000
223-9480	EZ Micro Test Tubes , 1.5 ml, natural, 500
223-9430	EZ Micro Test Tubes , 2 ml,* natural, 500
223-9501	Micro Test Tubes , standard, 1.5 ml, natural, 500
223-9500	Micro Test Tubes , capless, 1.5 ml, polypropylene, natural, graduated, 500
223-9490	Separate Caps , for capless micro test tubes, white, 1,000

* To fit centrifuges, 2 ml tubes have thinner walls than standard tubes and are intended for short runs at less than 13,000 x g. Heavy samples, high g forces, long runs, or centrifuges that radiate excess heat may damage these tubes.

Screwcap Micro Test Tubes

Screwcap micro test tubes and caps are made of polypropylene. O-rings are made of a blend of polyethylene and polypropylene. Features include:

- O-ring operating range from -55 to 150°C
- Uniform walls for uniform heat transfer
- Conical bottoms for pellet collection
- Knurled caps for easy handling



0.5 ml Conical 0.5 ml Skirted 1.5 ml Conical 2.0 ml Skirted Screwcap

Ordering Information

Catalog #	Description
224-0165	0.5 ml Conical Tubes , with installed O-ring screwcaps, sterilized, 500
224-0185	0.5 ml Skirted Tubes , with installed O-ring screwcaps, sterilized, 500
224-0100	1.5 ml Conical Tubes , with separate O-ring screwcaps, unsterilized, 500
224-0110	1.5 ml Conical Tubes , with installed O-ring screwcaps, sterilized, 500
224-0130	2.0 ml Skirted Tubes , with separate O-ring screwcaps, unsterilized, 500
224-0140	2.0 ml Skirted Tubes , with installed O-ring screwcaps, sterilized, 500

A

accessories, 8, 21, 35–48
 autoclaving, 41–43, 46, 47
 automation, products for, 10–12,
 16–20, 27–32
 auto-sealing lids, 27, 31, 32

B

bar-coded plates, 11, 12, 16–20, 26

C

cap strips, 4, 6, 7, 22–25
 capping tools, 8
 Chill-out™ liquid wax, 27, 33
 Concord™ 96-well PCR plates, 17,
 22–26
 contamination, 2, 5–7, 28–32, 34,
 37–39

E

Easy Cap™ tool, 8
 EZ Micro™ test tubes, 46, 47

F

Frame-Seal™ chambers, 34

H

Hard-Shell® 96-well PCR plates,
 10–12, 22–26

Hard-Shell 384-well PCR plates, 18,
 19, 22–26
 heat sealer support plate, 21

I

in situ PCR, *See* sealers, for slides
 iQ™ 96-well PCR plates, 15, 22–26

M

Microseal® 'A' film, 27, 28, 30
 Microseal 'B' seals, 27–30
 Microseal 'F' foil, 27, 29, 30
 Microseal 'P' and 'P+' pads, 27, 31, 32
 Microseal 96-well PCR plates,
 16, 17, 22–26
 Microseal 384 plate positioner, 21
 Microseal 384-well PCR plates,
 20, 22–26
 Multiplate™ 48-well PCR plates,
 9, 22–26
 Multiplate 96-well PCR plates,
 13, 22–26

O

optical assays, sealers for, 4, 6, 7,
 22–25, 33

P

PCR plates
 48-well, 9, 22–26

96-well, 10–17, 22–26
 384-well, 18–20, 22–26
 bar-coded, 10–12, 18, 19, 26
 full-height (high-profile), 12, 13,
 15–17, 22–26
 low-profile, 9, 11, 14, 16, 17, 22–26
 polycarbonate, 17, 22–26
 semi-skirted, 10, 12, 15–17, 22–26
 skirted, 10, 11, 16–19, 22–26
 unskirted, 9, 10, 13, 14, 22–26

pipet tips, 35–43
 process sampling, 2
 Prot/Elec™ tips, 38, 39

R

racks, 8, 38, 39, 41–43, 45
See also accessories
 reaction vessels, *See* tubes,
See also PCR plates
 reagent reservoirs, 44

S

sealers
 caps, 4–7
 film, 27, 28, 30
 foil, 27, 29, 30
 for slides, 34
 pressure pad, 29, 30

reusable
 auto-sealing lids, 27, 31, 32
 mats, 27, 29, 30
 pads for motorized lids, 27, 31, 32
 tape, 30
 wax, 27, 33
 selection charts, 4, 22–27, 39, 42
 Seque/Pro™ tips, 38, 39
 slide sealers, *See* sealers, for slides
 storage, 6, 8, 27–33, 45

T

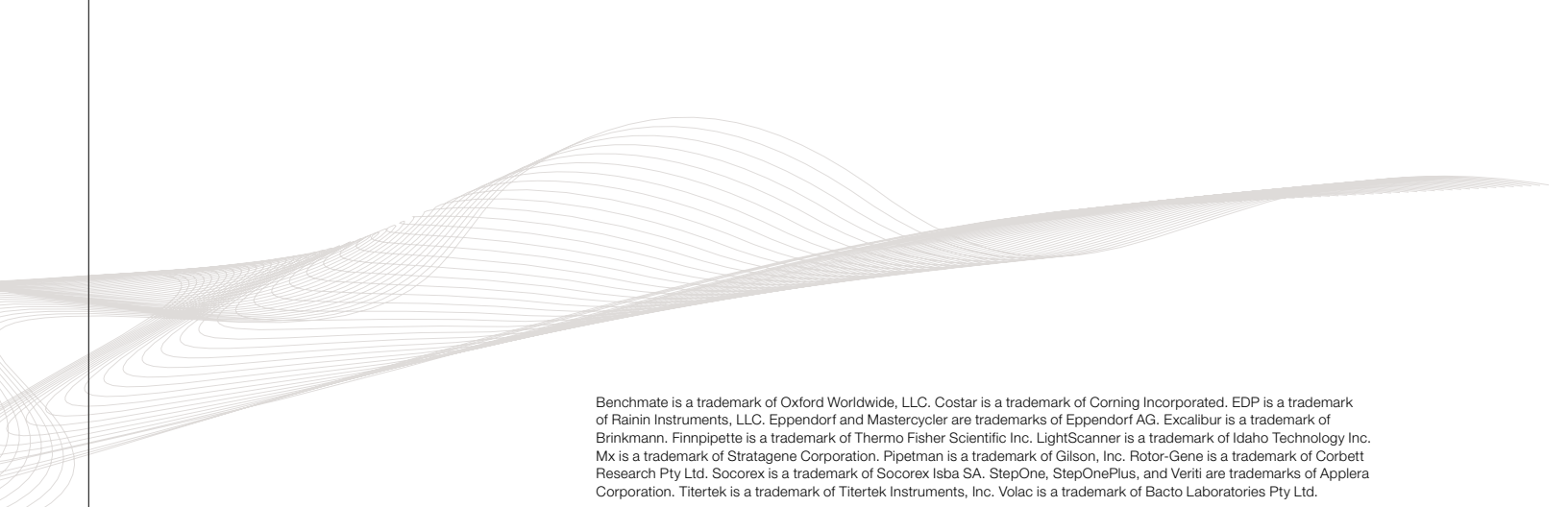
tips, pipet, *See* pipet tips
 Titertube® tubes, 45
 tubes

0.2 ml
 high-profile, 4–7
 individual, 5
 low-profile, 4, 6, 7
 screwcap, 48
 standard, 46, 47
 strips, 4, 6, 7, 22, 24
 0.5 ml, 4, 5
 micro test tubes, 45–48

X

Xcluda™ tips, 37, 39





Benchmate is a trademark of Oxford Worldwide, LLC. Costar is a trademark of Corning Incorporated. EDP is a trademark of Rainin Instruments, LLC. Eppendorf and Mastercycler are trademarks of Eppendorf AG. Excalibur is a trademark of Brinkmann. Finnpiquette is a trademark of Thermo Fisher Scientific Inc. LightScanner is a trademark of Idaho Technology Inc. Mx is a trademark of Stratagene Corporation. Pipetman is a trademark of Gilson, Inc. Rotor-Gene is a trademark of Corbett Research Pty Ltd. Socorex is a trademark of Socorex Isba SA. StepOne, StepOnePlus, and Veriti are trademarks of Applied Biosystems. Titertek is a trademark of Titertek Instruments, Inc. Volac is a trademark of Bacto Laboratories Pty Ltd.

Notice regarding Bio-Rad thermal cyclers and real-time systems:

Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in human in vitro diagnostics and all other applied fields under one or more of U.S. Patents 5,656,493, 5,333,675, 5,475,610 (claims 1, 44, 158, 160-163 and 167 only), and 6,703,236 (claims 1-7 only), or corresponding claims in their non-U.S. counterparts, owned by Applied Biosystems. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied Biosystems' United States Patent 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

Bio-Rad's thermal cyclers and real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Nos. 6,767,512 and 6,074,367.

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 Nos. 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 31 3689 6600
Canada 905 364 3435 **China** 86 20 8732 2339 **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** 30 210 777 4396 **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** 0508 805 500
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 **United Kingdom** 020 8328 2000