

# Multiplate™ 96- and 48-Well Unskirted PCR Plates

Multiplate PCR plates are constructed of thin-wall polypropylene for low protein binding and excellent sample retention. When fewer than 96 wells are needed, a Multiplate PCR plate can easily be cut with scissors to the required size. They are designed for compatibility with a wide variety of sealing systems for oil-free cycling of 5–125 µl reactions. They are free of DNase, RNase, and human DNA.

## High-Profile (Full-Height) PCR Plates

The versatile unskirted design with full-height (20.70 mm) wells makes these plates compatible with many instruments.

- PCR volumes of 5–125 µl are recommended
- Maximum well volume of 300 µl

### 96-well, 8 x 12 well array, 25 plates/pkg

MLP-9601 Natural

### 48-well, 6 x 8 well array, 50 plates/pkg

MLP-4801 Natural

## Low-Profile PCR Plates

The lower overall height (15.50 mm) of low-profile plates 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.

- PCR volumes of 5–125 µl are recommended
- Maximum well volume of 200 µl
- Available in natural and opaque white colors

### 96-well, 8 x 12 well array, 25 plates/pkg

MLL-9601 Natural  
MLL-9651 White

### 48-well, 6 x 8 well array, 50 plates/pkg

MLL-4801 Natural  
MLL-4851 White

## Recommended Sealing Options for Unskirted Plates

### Strip Caps

TCS-0803 **Optical Flat 8-Cap Strips**, ultraclear, 120  
TCS-0801 **Domed 8-Cap Strips**, natural, 120  
TCS-1201 **Domed 12-Cap Strips**, natural, 200

### Sealing Films

MSA-5001 **Microseal® 'A' Film**, 50  
MSB-1001 **Microseal 'B' Adhesive Seals**,  
optically clear, 100  
MSF-1001 **Microseal 'F' Foil**, 100

### Other Sealing Options

223-9442 **96-Well Plate Sealing Mats**, reusable, 5  
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

### Other Accessories

TRC-0501 **96-Place Racks**, with covers, 5  
ECT-1000 **Easy Cap™ Tool**, ensures tight seal for  
0.2 ml tubes or 96-well plates  
ECT-2000 **Strip Cap Tool**, for seating 8- and  
12-cap strips  
MSR-0001 **Sealing Roller**, for film seals



## Instrument Compatibility

|   | High-Profile PCR Plates | Low-Profile PCR Plates |
|---|-------------------------|------------------------|
| Catalog number  | MLP-xxxx                | MLL-xxxx               |
| Well height   | 20.70 mm                | 15.50 mm               |
| PCR reaction volume   | 5–125 µl                | 5–125 µl               |
| Maximum well volume   | 300 µl                  | 200 µl                 |
| <b>Thermal Cycler Compatibility</b>   |                         |                        |
| Bio-Rad C1000™/S1000™   | •                       | •                      |
| Bio-Rad MJ Mini™  | •                       | •                      |
| Bio-Rad DNA Engine®, DNA Engine Tetrad® and Tetrad 2, DNA Engine Dyad®, and Dyad Disciple™      | •                       | •                      |
| Bio-Rad iCycler® and MyCycler™  | •                       |                        |
| Bio-Rad PTC-100®  | •                       | •                      |
| Life Technologies cyclers with regular 0.2 ml blocks (2720, 9700, Veriti, etc.)                 | •                       |                        |
| Life Technologies cyclers with fast 0.1 ml blocks (9800 fast, Veriti fast, etc.)                |                         | •                      |
| Eppendorf Mastercycler series   | •                       | •                      |
| <b>Real-Time PCR Detection System Compatibility</b>   |                         |                        |
| Bio-Rad CFX96™  |                         | •                      |
| Bio-Rad MiniOpticon™  |                         | •                      |
| Bio-Rad iQ™5, iCycler iQ®, MyiQ™, and MyiQ™2  | •                       |                        |
| Bio-Rad Chromo4™  | •                       | •                      |
| Bio-Rad DNA Engine Opticon® and Opticon 2   |                         | •                      |
| Life Technologies systems with regular blocks (7300, 7500, 7900HT, etc.)                        | •                       |                        |
| Life Technologies systems with fast blocks (7500 fast, 7900HT fast, StepOne, StepOnePlus, etc.) |                         | •                      |
| Eppendorf Mastercycler ep <i>realplex</i>   | •                       | •                      |
| Stratagene Mx series  | •                       |                        |
| <b>Other Instrument Compatibility</b>   |                         |                        |
| Life Technologies DNA sequencers (3700, 3730, etc.)   | •                       |                        |
| Idaho Technology LightScanner   | •                       | •                      |



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.

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's 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 7,074,367.

LightScanner is a trademark of Idaho Technology Inc. Mastercycler is a trademark of Eppendorf AG. Mx is a trademark of Stratagene Corporation. StepOne, StepOnePlus, and Veriti are trademarks of Applied Biosystems.

