

Hard-Shell® 480 PCR Plates

- 96-well
- 384-well

Premium Plates Specifically Designed for the Roche LightCycler 480

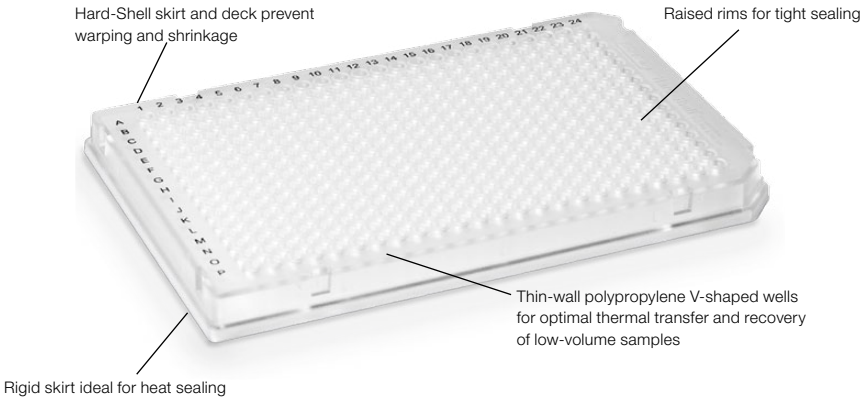
Hard-Shell PCR plates represent an innovative concept in PCR plate design. The skirt and deck are molded from a rigid thermostable polymer, preventing robotic handling problems due to plate warping. This hard shell prevents the distortion and shrinkage that may occur when regular single-component polypropylene PCR plates are exposed to the high temperatures of thermal cycling or heat sealing. The thin-wall wells are molded of virgin polypropylene selected for low DNA binding. The raised rims around each well allow tight sealing using a variety of procedures, including pressure-, adhesive-, and heat-sealing methods.



**Hard-Shell 96-Well
480 PCR Plate**



**Hard-Shell 384-Well
480 PCR Plate**



BIO-RAD

Hard-Shell 96-Well 480 PCR Plates and Seals

8- x 12-well array, for the LightCycler 480

HSR-9905	White well, clear shell, bar-coded on 1 side (Row A), 25 plates
HSR-9905K	White well, clear shell, bar-coded on 1 side (Row A), 100 plates and 100 Microseal® 'C' optical seals
HSR-9901	Clear well, clear shell, bar-coded on 1 side (Row A), 25 plates
HSR-9901K	Clear well, clear shell, bar-coded on 1 side (Row A), 100 plates and 100 Microseal 'C' optical seals

Hard-Shell 384-Well 480 PCR Plates and Seals

16- x 24-well array, for the LightCycler 480



HSR-4805	White well, clear shell, bar-coded on 1 side (Row A), 50 plates
HSR-4805K	White well, clear shell, bar-coded on 1 side (Row A), 100 plates and 100 Microseal 'C' optical seals
HSR-4801	Clear well, clear shell, bar-coded on 1 side (Row A), 50 plates
HSR-4801K	Clear well, clear shell, bar-coded on 1 side (Row A), 100 plates and 100 Microseal 'C' optical seals

Seals and Accessories

181-4000	PX1™ PCR Plate Sealer , semiautomated heat sealing instrument
181-4030	Optically Clear Heat Seal , for use with PX1 PCR plate sealer, 100 seals
MSB-1001	Microseal 'B' Adhesive Seals , optically clear, 100 seals
MSC-1001	Microseal 'C' Optical Seals , 100 adhesive seals
MSR-0001	Sealing Roller , for adhesive film seals
CHO-1411	Chill-out™ Liquid Wax , clear, optical grade, 100 ml

Please request bulletin 5496 for more information about Hard-Shell PCR plates, including dimension details, and bulletin 6090 for more information about seals and accessories.

Hard-Shell PCR Plates for the LightCycler 480

	96-Well 480 PCR Plates	384-Well 480 PCR Plates
		
Catalog number	HSR-99xx	HSR-48xx
Maximum well volume	200 µl	50 µl
ANSI/SBS standard dimensions	Yes	Yes
Autoclave specification	121°C, 20 min	
Centrifuge specification	Up to 2,250 x g in a swing-out rotor	
Application temperature	–80 to 121°C	
Storage	Room temperature; avoid sun and UV light	
Cleanroom production	Free of DNase, RNase, and human DNA	

Please request bulletin 5496 for more information about instrument compatibility with Hard-Shell 96-well and 384-well 480 PCR plates.



Microseal 'C' optical seals are optically clear adhesive films designed for optical assays, such as those performed in real-time PCR. They can be used for microplates with raised well rims.

Key Features

- Films provide a tight seal for PCR and qPCR
- Pressure- and heat-sensitive adhesive sticks to the plate, not your gloves
- Free from DNase, RNase, and human DNA contaminants

LightCycler is a trademark of Roche Diagnostics GmbH.

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 Numbers 7,347,977; 6,340,589; and 6,528,302.

Practice of the polymerase chain reaction (PCR) may require a license.



**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 11 5044 5699 **Canada** 905 364 3435 **China** 86 21 6169 8500
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 9532 220 **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 64 9 415 2280 **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 **Thailand** 800 88 22 88
United Kingdom 020 8328 2000