

# Polypropylene Reservoir Microplates

## For automated liquid handlers and handheld pipettors

All microplates are molded using chemical-resistant polymers to ANSI/SLAS standard footprints, with optimized base geometries to minimize dead volumes. They are available with one to 24 cavities per plate, with container heights between 19 and 44 mm, and with clear or black lids. Recommended for method scale-up and scale-down applications, and for creating design of experiment (DoE) arrays.

To add items to your "My Favorites"\* list at the [Agilent online store](#), simply click the part number. Then, enter the quantities for the products you need.



The polypropylene reservoir microplates listed below are organized by the number of cavities in the reservoir and the volume in each cavity.

Description	Cavity Volume (mL)	Part Number
Single cavity polypropylene, 96 pyramids base geometry, 44 mm height, case of 25	300	<a href="#">201244-100</a>
Single cavity polypropylene, 96 pyramids base geometry, 44 mm height, case of 5, sample pack of 201244-100	300	<a href="#">204756-100</a>
Single cavity black polypropylene, 96 pyramids base geometry, 44 mm height, case of 25	300	<a href="#">201246-100</a>
Single cavity irradiated polypropylene, 96 pyramids base geometry, 44 mm height, case of 25	300	<a href="#">204017-100</a>
Single cavity polypropylene, 12-column base geometry, 44 mm height, case of 25	290	<a href="#">201250-100</a>
Single cavity irradiated polypropylene, 12-column base geometry, 44 mm height, case of 25	290	<a href="#">204093-100</a>
Single cavity polypropylene, 8-row base geometry, 44 mm height, case of 25	290	<a href="#">201252-100</a>
Single cavity polypropylene, 8-row base geometry, 44 mm height, irradiated, case of 25	290	<a href="#">204504-100</a>
Single cavity polypropylene, 384 pyramid base geometry, 44 mm height, case of 25	282	<a href="#">201264-100</a>
Single cavity polypropylene, 384 pyramid base geometry, 30.6 mm height, case of 25	170	<a href="#">201302-100</a>
Single cavity polypropylene, 384 pyramid base geometry, 30.6 mm height, irradiated, case of 25	170	<a href="#">204377-100</a>
Single cavity polypropylene, 12-column base geometry, 30.6 mm height, case of 25	170	<a href="#">200686-100</a>
Single cavity polypropylene, 12-column base geometry, 30.6 mm height, irradiated, case of 25	170	<a href="#">204363-100</a>

Description	Cavity Volume (mL)	Part Number
Single cavity polypropylene, 384 pyramid base geometry, 19 mm height, case of 25	92	201266-100
Single cavity polypropylene, 384 pyramid base geometry, 19 mm height, irradiated, case of 25	92	204612-100
Single cavity polypropylene, flat base geometry, 19 mm height, case of 25	90	201270-100
Single cavity polypropylene, 96 pyramids base geometry, 19 mm height, case of 25	86	201254-100
Single cavity polypropylene, 96 pyramids base geometry, 19 mm height, irradiated, case of 25	86	204484-100
2-column polypropylene, 292 mL max, pyramid base geometries, 44 mm height, case of 25	146	203852-100
2-column polypropylene, 292 mL max, pyramid base geometries, 44 mm height, irradiated, case of 25	146	204359-100
3 column polypropylene, 285 mL max, pyramid base geometries, 44 mm height, case of 25	95	204249-100
3 column polypropylene, 285 mL max, pyramid base geometries, 44 mm height, irradiated, case of 25	95	204361-100
4 column polypropylene, 292 mL max, pyramid base geometries, 44 mm height, case of 25	73	201308-100
4 column polypropylene, 292 mL max, pyramid base geometries, 44 mm height, irradiated, case of 25	73	204282-100
6 column polypropylene, 282 mL max, 24 pyramid base geometries, 44 mm height, case of 25	47	201284-100
6 column polypropylene, 282 mL max, 24 pyramid base geometries, 44 mm height, irradiated, case of 25	47	204284-100
12 column polypropylene, 252 mL max, pyramid base geometries, 44 mm height, case of 25	21	201256-100
12 column polypropylene, 252 mL max, pyramid base geometries, 44 mm height, case of 5, sample pack of 201256-100	21	204758-100
12 column irradiated polypropylene, 252 mL max, pyramid base geometries, 44 mm height, case of 25	21	204095-100
12 column polypropylene, 84 mL max, pyramid base geometries, 19 mm height, case of 25	7	201280-100
12 column polypropylene, 84 mL max, pyramid base geometries, 19 mm height, irradiated, case of 25	7	204365-100
24 column polypropylene, 78 mL max, pyramid base geometries, 19 mm height, case of 25	3.25	201296-100
24 column polypropylene, 78 mL max, pyramid base geometries, 19 mm height, irradiated, case of 25	3.25	204373-100
4 row polypropylene, 292 mL max, pyramid base geometries, 44 mm height, case of 25	73	201312-100
4 row polypropylene, 292 mL max, pyramid base geometries, 44 mm height, irradiated, case of 25	73	204280-100
8 row polypropylene, 256 mL max, pyramid base geometries, 44 mm height, case of 25	32	201260-100
8 row irradiated polypropylene, 256 mL max, pyramid base geometries, 44 mm height, case of 25	32	204019-100
8 row polypropylene, 85 mL max, pyramid base geometries, 19 mm height, case of 25	10.6	201282-100
8 row polypropylene, 85 mL max, pyramid base geometries, 19 mm height, irradiated, case of 25	10.6	204367-100
16 row polypropylene, 78 mL max, pyramid base geometries, 19 mm height, case of 25	4.9	201294-100
16 row polypropylene, 78 mL max, pyramid base geometries, 19 mm height, irradiated, case of 25	4.9	204371-100
Unique 3 cavity polypropylene, 2 control wells (7 mL col 24 rows A-L and 2 mL col 24 rows M-P) and 273 mL main, 384 pyramid base geometries, 44 mm height, case of 25		201286-100
Unique 5 cavity polypropylene, 4 control wells (cols 23-24 rows A-H, I-P) of 2.6 mL each and 145 mL main, 384 pyramid base geometries, 30.6 mm height, case of 25		201300-100
Unique 5 cavity polypropylene, 4 control wells (cols 1,2,11,12) of 3.25 mL each and 86.3 mL main, 24 pyramid base geometries, 19 mm height, case of 25		201278-100
Unique 24 cavity polypropylene, 2 rows x 12 columns, 3.5 ml/cavity, 84 mL max, pyramid base geometries, 19 mm height, case of 25		201298-100
Unique 24 cavity polypropylene, 2 rows x 12 columns, 3.5 ml/cavity, 84 mL max, pyramid base geometries, 19 mm height, irradiated, case of 25		204375-100

\*First time using "My Favorites?" You will be asked to enter your email address for account verification. If you have an existing Agilent account, you will be able to log in. However, if you don't have a registered Agilent account, you will need to [register](#) for one. This feature is valid only in regions that are e-commerce enabled. All items also can be ordered through your regular sales and distributor channels.

See the difference for yourself. Request your free sample:

[explore.agilent.com/microplate-sample](https://www.agilent.com/microplate-sample)

Find microplates for any workflow:

[www.agilent.com/chem/microplates](https://www.agilent.com/chem/microplates)

DE44280.4374074074

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021  
 Printed in the USA, September 26, 2021  
 5994-3224EN

