

Assay, Collection, and Storage Microplates

For reaction chemistries, biological assays, low-temperature storage, and filtrates

All microplates are molded with medical-grade, chemical-resistant polymers and are available in 24-, 48-, 96-, or 384-well configurations of variable volumes. Mat seals and lids are also available.

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 assay, collection and storage microplates listed below are organized by well number and well volume.

Description	Well Volume (mL)	Part Number
24-well polypropylene, square well, pyramid bottoms, 44 mm height, case of 25	10	201272-100
24-well polypropylene, square well, round bottoms, 44 mm height, irradiated, case of 25	10	202061-100
24-well polypropylene, square well, round bottoms, 44 mm height, case of 25	10	202061-300
24-well irradiated polypropylene, square well, pyramid bottoms, 44 mm height, case of 25	10	204023-100
24-well polypropylene, square well, round bottoms, 44 mm height, case of 5, sample pack of 202061-100	10	204754-100
48-well (6 rows × 8 columns) polypropylene, square well, pyramid bottoms, 44 mm height, case of 25	4.8	201306-100
48-well (8 rows × 6 columns) polypropylene, rectangle well, pyramid bottoms, 44 mm height, case of 25	5	201238-100
48-well (8 rows × 6 columns) metallocene polypropylene, rectangle well, pyramid bottoms, 44 mm height, case of 25	5	203903-100
48-well (8 rows × 6 columns) polypropylene, rectangle well, pyramid bottoms, 44 mm height, irradiated, case of 25	5	204351-100
48-well (8 rows × 6 columns) polypropylene, rectangle well, pyramid bottoms, 68 mm height, case of 15	7.5	201236-100
96-well polypropylene, round well, round bottom, 14.3 mm height, case of 50	0.5	203942-100
96-well polypropylene, round well, round bottom, 14.3 mm height, case of 25	0.5	204600-100
96-well polypropylene, round well, round bottom, 14.3 mm height, irradiated, case of 25	0.5	204601-100

Description	Well Volume (mL)	Part Number
96-well polypropylene, round well, round bottom, 14.3 mm height, irradiated, case of 50	0.5	204602-100
96-well polypropylene, square well, flat bottoms, 19 mm height, case of 50	0.7	201242-100
96-well polypropylene, round well, round bottoms, 32 mm height, case of 25	1	203426-100
96-well, polypropylene, round well, round bottoms, 32 mm height, case of 5, sample pack of 203426-100	1	204755-100
96-well polypropylene, round well, round bottoms, 32 mm height, irradiated, case of 25	1	204357-100
96-well polypropylene, round well, round bottoms, 32 mm height, white, case of 25	1	204392-100
96-well polypropylene, square well, V bottoms, 30.6 mm height, case of 25	1.1	201276-100
96-well polypropylene, square well, V bottoms, 30.6 mm height, irradiated, case of 25	1.1	204355-100
96-well, ultra-high purity polypropylene, square well, V bottoms, 31 mm height, case of 5, sample pack of 201276-100	1	204753-100
96-well polypropylene, square well, pyramid bottoms, 44 mm height, case of 25	2	201240-100
96-well, polypropylene, square well, pyramid bottoms, 44 mm height, case of 5, sample pack of 201240-100	2	204751-100
96-well ultrahigh purity polypropylene, square well, pyramid bottoms, 44 mm height, case of 25	2	201379-100
96-well polypropylene, square well, pyramid bottoms, 44 mm height, irradiated, case of 25	2	204353-100
96-well ultrahigh purity polypropylene, square well, pyramid bottoms, 44 mm height, irradiated, case of 25	2	204379-100
384-well polypropylene, square well, spherical bottoms, 14.5 mm height, case of 60	0.058	201288-100
384-well black polypropylene, square well, spherical bottoms, 14.5 mm height, case of 60	0.058	201290-100
384-well polypropylene, square well, spherical bottoms, 19 mm height, case of 10	0.2	204586-100
384-well polypropylene, square well, spherical bottoms, 19 mm height, irradiated, case of 10	0.2	204587-100

*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

Find microplates for any workflow:

www.agilent.com/chem/microplates

DE44280.4346527778

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021
 Printed in the USA, September 27, 2021
 5994-3222EN

