



Better microplates means better results

Microplates for storage

In the day-to-day hustle of today's lab it's easy to overlook the importance of proper compound storage. After all, you want to ensure your samples and compounds will be protected until you are ready to use them, and for that not just any microplate will do. Storage plates need to be resistant to the types of solvents that may be used to preserve chemical libraries.

Revvity storage plates are made of polypropylene which is both resistant to the solvents used in compound storage and suitable for use at temperatures down to -80 °C. Each StorPlate™ is DNase, RNase and pyrogen free with support certification available on request.



These plates are offered with a variety of options, including U-bottom and V-bottom wells, and deep-well plates for higher volumes.

Product	Features	Maximum volume without StorMat	Maximum volume with StorMat	Well number	Case of	Part number
StorPlate-96 U	U-shaped well bottom	450 µL	350 µL	96-well	50	6008190
					200	6008199
StorPlate-96 V	V-shaped well bottom	450 µL	300 µL	96-well	50	6008290
					200	6008299
Deep-well StorPlate - 96 U (1.1 mL)	Deep well, U-shaped bottom	1.1 mL	1.01 mL	96-well	50	6008390
Deep-well StorPlate - 96 U (2.0 mL)	Deep well, U-shaped bottom	2.2 mL	2.0 mL	96-well	50	6008490
StorPlate-384 V	Deep well, V-shaped bottom	120 µL	90 µL	384-well	50	6008590
					200	6008598
Deep-well StorPlate-384 V	Deep well, V-shaped bottom	240 µL	200 µL	384-well	50	6008690
StorMat-96	Silicone Microplate lid, Piercable with Self-sealing ports.	-	-	96-well	50	6008096
StorMat-384	Silicone Microplate lid, Piercable with Self-sealing ports.	-	-	384-well	50	6008384

Please ask a representative for more information or visit www.revvity.com.