

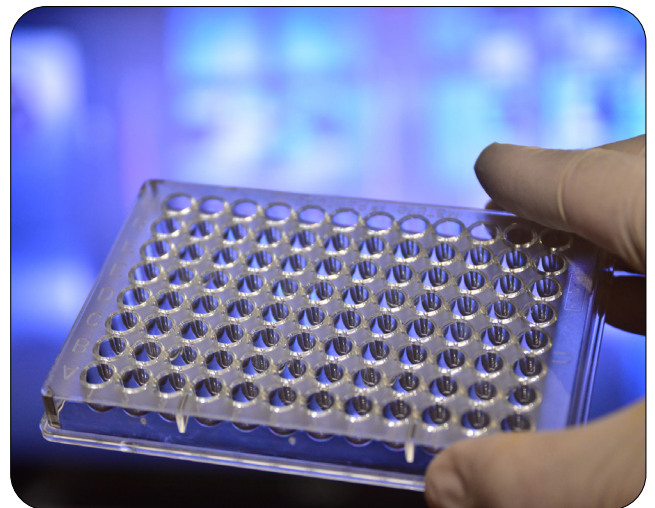


Discover better.

Choosing the right microplate is a critical, often overlooked, part of an assay. The right microplate helps provides valuable data, whereas the wrong microplate can lead to missed or inaccurate data leading to missed project timelines and ultimately higher costs.

New technologies for drug development are advancing at record speed. These technologies have given rise to a variety of applications that require their own unique consumables. At Revvity, we provide microplates that help you discover better. Covering a vast array of application areas, Revvity microplates are available in different colors, well format, plastic type, as well as microplate coatings.

Our microplates have been engineered to deliver the highest quality data for applications ranging from high throughput screening to imaging.



Storage

Microplates for storing reagents, biomolecules or other samples



Imaging

Microplates for microscopy and high content screening offering the lowest plate bottom for superior image acquisition



Research

Microplates for TR-FRET, HTRF, Alpha, Radiometric, luminescence, fluorescence, and others for general research



Cell culture

Microplates designed to improve consistent and reliable cell growth for general cell culture as well as 3D cell culture



NGS

Microplates for Next Generation Sequencing used in sample preparation workflows

Microplates														
Application	1/2 Area Plate	Alpha Plate*	Pheno Plate†	CellCarrier Spheroid	Cultur Plate	DELTA	Hardshell PCR	Iso Plate	Opti Plate	Proxi Plate	Spectra Plate	Stor Plate	View Plate	Visi Plate
Storage												✓		
Imaging			✓	✓									✓	
Fluorescence Assays	✓				✓	✓		✓	✓	✓			✓	✓
Luminescence Assays	✓	✓			✓			✓	✓	✓			✓	✓
Absorbance/ Colorimetrics Assays	✓							✓			✓		✓	
Cell Culture			✓	✓	✓			✓**			✓**		✓**	✓**
NGS							✓							

* Select plates available in 1/2Area

** Tissue-culture treated only

† Formerly known as CellCarrier Ultra Plates

Available Coatings

At the surface of a microplate well, interaction between a sample and the microplate takes place. It's important to choose the correct surface coating to maximize your success. Revvity's reputation for high quality precision microplates extends to expertise in a variety of surface chemistries. Depending on your application, you can choose microplates pre-coated with popular surface chemistries or Revvity's technical support staff can assist you in the selection and qualification of the right microplate coating. We also offer custom microplate services which provide services for custom coatings as well as microplate barcoding.

Coatings for Culturing Cells

Revvity offers a variety of options for general cell culture needs. Our microplates undergo special physical surface treatment which results in improved and consistent cell attachment.

Tissue-culture (TC) treatment allows for cell attachment and binding to the bottom surface of the microplate. It is used for assays that use adherent cell lines.

Poly-D-lysine (PDL) treatment enhances cell attachment and binding. It is used when working with cells that are difficult to attach and when wash steps are needed.

Collagen (COL) treatment enhances cell attachment and proliferation. It is used with working with keratinocytes and hepatocytes.

Ultra-low attachment (ULA) treatment allows for further reduction of non-specific binding. It is used in cell colony high content screening assays, 3D cultures and other imaging applications.

Immunological Coatings

Generally, for immunology-based applications, immobilization of biomolecules to the well surface of a microplate is required. Revvity offers both a low and high binding surface. The high binding surface features a relatively high number of polar groups, whereas the number of polar groups is limited on the low binding surface.

High binding (HB) treatment allows for capturing proteins and antibodies to the microplate. It is used for washed-based assays such as ELISA and DELFIA.

Low binding (LB) treatment results in reduced binding to proteins and nucleic acids. It is used for biochemical assays to reduce non-specific binding.

Bar coding and other custom services available.

Contact us at <https://www.revvity.com/contact-us#sales> for more information.

TopSeal

Revvity's TopSeal™ is a range of plate seals that are applied to the top surface of the plate and are used to prevent evaporation or radioactive contamination during assay incubation steps and/or plate reading measurements. It is available as either a press-on adhesive or heat-activated seal and can be used for a wide array of applications in place of a lid.

Microplates for Radiometric Detection

Revvity offers the complete solution for your radiometric assay needs. Assays using radiochemicals can be run in various high-throughput formats, including filtration, liquid scintillation counting and scintillant coated- plate assays, and proximity assays. Whether you are reading samples using any of our plate based radiometric instruments or for ultra-HTS of SPA assays, we offer the ideal microplate for your radiometric assays.

For more information visit us at www.revvity.com

The Revvity logo is displayed in a lowercase, sans-serif font. The letters are black and have a clean, modern appearance. The 'v' and 'y' have a slight curve to them. The logo is positioned in the lower right quadrant of the page, above a large yellow wavy graphic that spans the bottom of the page.