

Better imaging microplates means better results.

Phenoplates™ for high-content screening

To get the best results from your high-content assays you need the best imaging microplates.

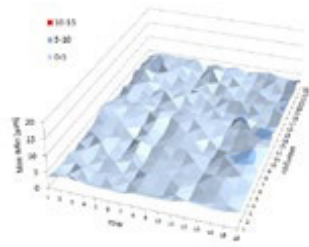
Drawing on many years of experience in high-content screening, Revvity's expert team have developed and validated the Phenoplate™ line of microplates.*

Our Phenoplates have been engineered to deliver the **highest performance** and **superior image quality** for high-content applications.



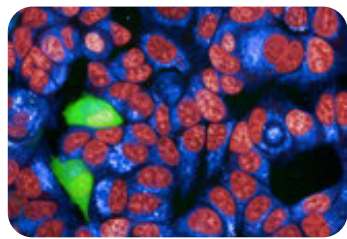
Optimal clarity and fast auto-focusing

Due to flatness of the plate bottom



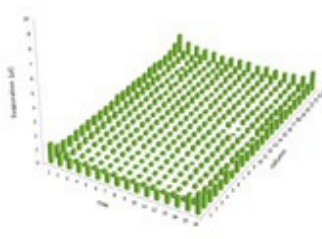
Superior image quality

From high optical quality of cyclic olefin imaging surface



Reduced evaporation

From newly designed, low-profile polystyrene lid



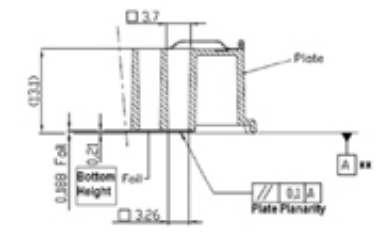
Corner spacers

Avoid damaging the imaging surface when stacking with unique corner spacers



Better well access

Ultra-low plate bottom provides access to all wells when using high NA water immersion objectives



* Phenoplates were formerly known as CellCarrier Ultra™ microplates.

96-well PhenoPlates	Part numbers			
Description	Case of 10	Case of 40	Case of 50	Case of 160
PhenoPlate 96-well, black, optically clear flat-bottom, tissue-culture treated, lids		6055302	6055300	6055308
PhenoPlate 96-well, black, optically clear flat-bottom, poly-D-lysine-coated, lids		6055500		6055508
PhenoPlate 96-well, black, optically clear flat-bottom, fibronectin-coated, lids	6055600	6055602		
PhenoPlate 96-well, black, optically clear flat-bottom, collagen-coated, lids		6055700		6055708
PhenoPlate 96-well, black, optically clear flat-bottom, ultra-low-attachment-coated, lids	6055800	6055802		

384-well PhenoPlates	Part numbers			
Description	Case of 10	Case of 40	Case of 50	Case of 160
PhenoPlate 384-well, black, optically clear flat-bottom, tissue-culture treated, lids		6057302	6057300	6057308
PhenoPlate 384-well, black, optically clear flat-bottom, tissue-culture treated, non-irradiated, lids				6057328
PhenoPlate 384-well, black, optically clear flat-bottom, poly-D-lysine-coated, lids		6057500		6057508
PhenoPlate 384-well, black, optically clear flat-bottom, fibronectin-coated, lids	6057600	6057602		
PhenoPlate 384-well, black, optically clear flat-bottom, collagen-coated, lids		6057700		6057708
PhenoPlate 384-well, black, optically clear flat-bottom, ultra-low-attachment-coated, lids	6057800	6057802		

Custom microplate services	
Coatings for culturing cells	Description
Tissue culture (TC)	Allows for cell attachment and binding to the bottom surface of the microplate (for adherent cells)
Poly-D-Lysine (PDL)	Enhances cell attachment and binding for cells that are difficult to attach or when wash steps are needed
Collage (COL)	Promotes cell attachment and proliferation when keratinocytes and hepatocytes are used
Fibronectin	Provides a substrate that mimics the natural cellular environment and enhances cell adhesion (ideal for imaging of stem cells)
Ultra-low attachment (ULA)	For further reduction of nonspecific binding, used in cell-colony high-content screening assays, 3D cultures, and other imaging applications
Immunological coatings	
High binding (HB)	Allows for capturing proteins and antibodies to the microplate, for washed-based assays
Low binding (LB)	Results in reduced binding to proteins and nucleic acids for biochemical assays
Streptavidin coating	Perfect for time-resolved fluorescence applications
Custom barcoding	
Our microplates can be supplied with high-quality barcode labels. Waterproof, scratchproof, and DMSO-resistant, these plastic labels withstand temperatures up to -80 °C and are available in different label formats and barcode types. We can label all microplate sides, with multiple labels per plate, which can include custom sequences and information.	

Please ask a representative for more information or for a free sample pack. Alternatively, visit www.revivity.com

