



PhenoPlates for high-content screening

High-content screening has the capability to generate valuable, information-rich image data that bring greater insights into cellular biology and drug discovery screening.

Out of focus images and inaccurate data collection are common issues caused by microplates that are not carefully designed for the unique challenges of high-content screening. This can result in missed data points, sample re-screening, missed project timelines, and ultimately higher costs.

Drawing on many years of experience Revvity expert team has developed the PhenoPlate line of microplates (formerly known as CellCarrier™ Ultra plates). These microplates have been engineered to deliver superior images and the highest quality data for high-content applications.

- Optimal clarity and fast autofocus from excellent flatness of the plate bottom
- Superior image quality from high optical quality of cyclic olefin imaging surface

- Better well access when using water immersion and high NA objectives with ultra-low plate bottom
- Reduced evaporation from new design, low-profile polystyrene lid
- Avoid damaging the imaging surface when stacking with unique corner spacers
- Available with different coatings to suit your application

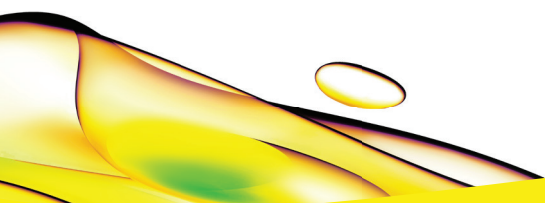
PhenoPlates are part of a range of microplates for high content applications, and belong to our complete solution for high content imaging that includes the Operetta®CLS™ high-content analysis and Opera Phenix®Plus high-content screening systems, software, and automation.

By using PhenoPlates, specifically engineered for high content applications, you can be confident that you'll acquire high quality images providing more information-rich data for easier analysis through unmatched image clarity.



Description	Part numbers				
	Case of 10	Case of 40	Case of 50	Case of 160	Case of 200
96-well PhenoPlates					
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, collagen-coated, lids		6055700		6055708	
PhenoPlate 96-well, black, optically clear flat-bottom, fibronectin-coated, lids	6055600	6055602			
PhenoPlate 96-well, black, optically clear flat-bottom, ultra-low-attachment-coated, lids	6055800	6055802			
384-well PhenoPlates	Case of 10	Case of 40	Case of 50	Case of 160	Case of 200
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, collagen-coated, lids		6057700		6057708	
PhenoPlate 384-well, black, optically clear flat-bottom, fibronectin-coated, lids	6057600	6057602			
PhenoPlate 384-well, black, optically clear flat-bottom, ultra-low-attachment-coated, lids	6057800	6057802			
PhenoPlate lids					
Lid, clear, for PhenoPlates, sterile					6057328

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



revvity