



## Cell culture microplates



### Better microplates means better results.

In research and drug discovery today, there is increasing demand for more physiologically relevant assays. As the need for more physiologically relevant assays has increased, so has the need for multi-well microplates designed to support the growth, attachment, and differentiation of cells, whether for ongoing culture or temporarily before cell-based assay. Growing and assaying cells on a single plate eliminates the need for a detachment step and therefore deems the assay more physiologically relevant.

Revvity's CulturPlates™ and PhenoPlates™ are specifically designed to promote cell attachment and growth. These microplates are appropriate for not only cell culture, but also imaging (PhenoPlates, formerly known as CellCarrier™ Ultra plates) or fluorescence and luminescence assays (CulturPlate).



Additionally, Revvity offers a variety of coatings that can be applied to a multitude of our plates to promote cell growth.

- Tissue-culture (TC) treatment allows for cell attachment and binding to the bottom surface of the microplate and is used for assays that use adherent cell lines.
- Poly-D-Lysine (PDL) treatment enhances cell attachment and binding and 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 and is used with working with keratinocytes and hepatocytes, specifically.
- Ultra-low attachment (ULA) treatment allows for further reduction of non-specific binding and is used in cell colony high content screening assays, 3D cultures and other imaging applications.

PhenoPlate										
Color	Feature	Lid	Well Number	Case of 10	Case of 20	Case of 40	Case of 50	Case of 80	Case of 160	Case of 200
Black	Collagen-coated	Yes	96-well			6055700			6055708	
			384-well			6057700			6057708	
	PDL-coated	Yes	96-well			6055500			6055508	
			384-well			6057500			6057508	
	Tissue culture-treated	Yes	96-well			6055302	6055300		6055308	
			384-well			6057302	6057300		6057308	
	ULA-coated	Yes	96-well	6055800		6055802				
			384-well	6057800		6057802				
	Non-irradiated tissue culture-treated	Yes	384-well						6057328	

CellCarrier										
Color	Feature	Lid	Well Number	Case of 10	Case of 20	Case of 40	Case of 50	Case of 80	Case of 160	Case of 200
Black	Tissue culture-treated	Yes	1536-well		6004550			6004558		

CellCarrier Spheroid										
Color	Feature	Lid	Well Number	Case of 10	Case of 20	Case of 40	Case of 50	Case of 80	Case of 160	Case of 200
Clear	Ultra Low Attachment, round bottom	Yes	96-well	6055330		6055334			6057328	

CulturPlate										
Color	Feature	Lid	Well Number	Case of 10	Case of 20	Case of 40	Case of 50	Case of 80	Case of 160	Case of 200
White	Tissue-culture treated	Yes*	24-well				6005168			
			96-well				6055680		6055688	6055689
			384-well				6007680		6007688	6007689
			1536-well				6004680			
Black	Tissue-culture treated	Yes*	96-well				6055660		6055668	6055669
			384-well				6007660		6007668	
			1536-well				6004660			

IsoPlate										
Color	Feature	Lid	Well Number	Case of 10	Case of 20	Case of 40	Case of 50	Case of 80	Case of 160	Case of 200
White	Tissue-culture treated, optically clear bottom	Yes	96-well				6005070		6005078	
Black	Tissue-culture treated, optically clear bottom	Yes	96-well				6005050			
White/Black	Tissue-culture treated, black frame, white bottom	Yes	96-well				6005060		6005068	

ProxiPlate										
Color	Feature	Lid	Well Number	Case of 10	Case of 20	Case of 40	Case of 50	Case of 80	Case of 160	Case of 200
White	Tissue-culture treated	Yes*	384-well				6008230		6008238	6008239
Black	Tissue-culture treated	Yes*	384-well				6008210			6008219

ScintiPlate										
Color	Feature	Lid	Well Number	Case of 10	Case of 20	Case of 40	Case of 50	Case of 80	Case of 160	Case of 200
White	Tissue-culture treated	Yes	96-well				6005390			

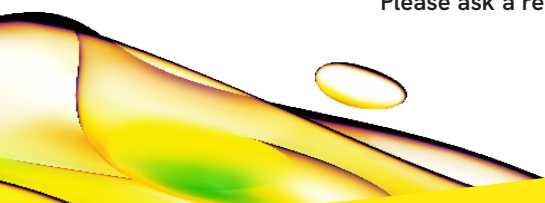
SpectraPlate										
Color	Feature	Lid	Well Number	Case of 10	Case of 20	Case of 40	Case of 50	Case of 80	Case of 160	Case of 200
Clear	Tissue-culture treated	Yes*	96-well				6055650		6055658	
			384-well				6007650		6007658	6007659
			1536-well				6004650			
	Tissue-culture treated, shallow well	Yes	384-well				6008650			

ViewPlate										
Color	Feature	Lid	Well Number	Case of 10	Case of 20	Case of 40	Case of 50	Case of 80	Case of 160	Case of 200
White	Tissue cultured treated, optically clear bottom	Yes	96-well				6005181			
			384-well			6007480				
			1536-well			6004480				
	Tissue cultured treated, optically clear bottom, ½Area	Yes	96-well			6005760			6005768	
Black	Tissue cultured treated, optically clear bottom	Yes	96-well				6005182	6005225 (2 sleeves of 25)		
			384-well			6007460				
			1536-well			6004460				
	Tissue-culture treated, glass bottom	Yes	96-well			6005430				
	Tissue-culture treated, PDL coated, optically clear bottom	Yes	384-well			6007710			6007718	
			1536-well	6004710						
	Tissue-culture treated, PDL coated, glassbottom	Yes	96-well	6005530 (case of 8)						
	Tissue-culture treated, Collagen coated, optically clear bottom	Yes	384-well			6007810			6007818	
			1536-well	6004810						
	Tissue-culture treated, Collagen coated, glassbottom	Yes	96-well	6005720 (case of 8)						

VisiPlate										
Color	Feature	Lid	Well Number	Case of 10	Case of 20	Case of 40	Case of 50	Case of 80	Case of 160	Case of 200
White	Tissue-culture treated, optically clear bottom	Yes	24-well	1450-603 (case of 14)			1450-604 (case of 56)			
Black	Tissue-culture treated, optically clear bottom	Yes	24-well	1450-605 (case fo 14)			450-606 (case of 56)			

\*Case of 200 does not include lids. Lids for 96-well plates are part number 6005619 and lids for 384-well plates are part number 6007619.

Please ask a representative for more information or for a free sample pack.



revvity