



revvity

Better microplates means better results.

Microplates for research

We know you are working hard to produce the next big breakthrough and the last thing you need to worry about is sub-par quality in something so commonly used, yet often overlooked, as a microplate. That's why our microplates are designed to give you optimal performance, whether you are working with fluorescence, luminescence or absorbancebased assays.

What sets our microplates apart from others? The manufacturing process. Our microplates are made using quality plastic injected into a mold in a clean room. This, together with our strict QC tolerance testing, ensures you are getting the best quality microplate possible. In addition, we offer a selection of treatments and coatings that fit all your research needs.

Try our OptiPlates™, AlphaPlate™, or ProxiPlates™, available in black, gray and white, for your next fluorescence, luminescence, or absorbance assay.

Pair our microplates with our range of detection reagents, plate readers, and automation technology for optimal assay performance and to simplify your workflow from plating to detection.



Microplates									
Application	AlphaPlate [®] ™	CulturPlate™	DELFLIA®	IsoPlate™	OptiPlate™	ProxiPlate™	SpectraPlate™	ViewPlate™	VisiPlate™
Fluorescence Assays		x	x	x	x	x		x	x
Luminescence Assays	x	x		x	x	x		x	x
Absorbance/Colorimetrics Assays				x			x	x	

OptiPlate						
Application	Color	Feature	Well number	Case of 50	Case of 200	
Fluorescence Luminescence	White	Standard untreated	96-well	6055290	6055299	
			384-well	6007290	6007299	
			1536-well	6004290		
		½Area, untreated	96-well	6002290	6002299	
			Low binding	96-well	6065420	
				384-well	6057480	
		½Area, low binding	96-well	6052480		
			High binding	96-well	6055240	6055249
		384-well		6005620	6005629	
		1536-well		6004620		
		Black	Standard untreated	96-well	6055270	6055279
				384-well	6007270	6007279
	1536-well			6004270		
	½Area, untreated		96-well	6002270	6002279	
			Low binding	96-well	6065400	
	384-well			6057260		
	½Area, low binding		96-well	6052260		
			High binding	96-well	6055220	6055229
	384-well	6005520		6005529		
	1536-well	6004520				
	Gray	High sensitivity	1536-well	6005540		
			1536-well	6004360		

CulturPlate						
Application	Color	Feature	Well number	Case of 50	Case of 160	Case of 200
Fluorescence Luminescence	White/opaque	Tissue-culture treated	96-well	6055680	6055688	6055689
			384-well	6007680	6007688	6007689
			1536-well	6004680		
	Black/opaque	Tissue-culture treated	96-well	6055660	6055668	6055669
			384-well	6007660	6007668	
			1536-well	6004660		

ProxiPlate						
Application	Color	Feature	Well number	Case of 50	Case of 160	Case of 200
Fluorescence Luminescence	White	Standard untreated	96-well	6006290		6006290
			384-well	6008280		6008289
		Low binding	384-well	6059480		
		Tissue-culture treated	384-well	6008230	6008238	6008239
	Black	Standard	96-well	6006270		6006279
			384-well	6008260		6008269
		Low binding	384-well	6059260		
		Tissue-culture treated	384-well	6008210		6008219

AlphaPlate						
Application	Color	Feature	Well Number	Case of 5	Case of 50	Case of 200
Luminescence	Light gray	Standard untreated	384-well		6005350	6005359
			1536-well		6004350	
		Shallow well	384-well	6008350		6008359
		Low binding	384-well		6057340	
		Low binding, shallow well	384-well		6059340	
		Low binding, ½Area	96-well		6052340	

SpectraPlate							
Application	Color	Feature	Well number	Case of 50	Case of 160	Case of 200	
Absorbance	Clear	Standard untreated	96-well	6005640		6005649	
			384-well	6007640		6007649	
			1536-well	6004640			
		½ Area	96-well	6002640		6002649	
		Shallow well	384-well	6008640		6008649	
		Low binding	96-well	6055630			
			384-well	6057630			
		Low binding, ½Area	96-well	6052630			
		Low binding, shallow well	384-well	6059630			
		High binding	96-well	6005600			6005609
			384-well	6007500			6007509
			1536-well	6004500			
		Tissue-culture treated	96-well	6005650			6005659
			384-well	6007650	6007658		6007659
			1536-well	6004650			
		Tissue-culture treated, shallow well	384-well	6008650			

ViewPlate										
Application	Color (frame/bottom)	Feature	Well Number	Case of 8	Case of 10	Case of 40	Case of 50	Case of 60	Case of 160	
Fluorescence Luminescence Absorbance	White/clear	Standard untreated	384-well				6007490			
			1536-well				6004490			
		Tissue-culture treated	96-well					6005181		
			384-well					6007480		
			1536-well					6004480		
			Tissue-culture treated, ½Area	96-well			6005760		6004470	6005768
	Black/clear	Standard untreated	384-well					6007470		
			1536-well							
		Tissue-culture treated	96-well					6005182		
			384-well				6007460			
			1536-well				6004460			
			Tissue-culture treated, glass bottom	96-well				6005430		
		PDL coated	96-well					6005710		
			384-well						6007710	6007718
			1536-well			6004710				
			PDL coated, glass bottom	96-well	6005530					
		Collagen coated	96-well					6005810		6005818
			384-well					6007810		6007818
			1536-well			6004810				
			Collagen coated, glass bottom	96-well	6005720					

1. Two sleeves of 25

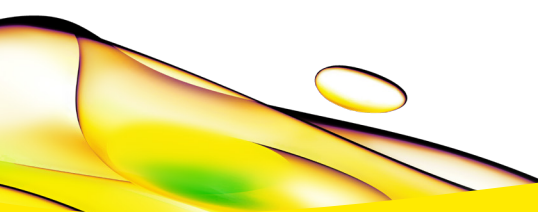
IsoPlate							
Application	Color (frame/well)	Feature	Well number	Case of 50	Case of 160	Case of 200	
Fluorescence Luminescence Absorbance	White/clear	Standard untreated	96-well	6005040		6005049	
		High binding	96-well	6005590		6005599	
		Tissue-culture treated	96-well	6005070	6005078		
	Black/clear	Standard untreated	96-well	6005020			6005029
		High binding	96-well	6005570			6005579
		Tissue-culture treated	96-well	6005050			
	Black/white	Standard untreated	96-well	6005030			6005039
		High binding	96-well	6005580			6005589
		Tissue-culture treated	96-well	6005060	6005068		

VisiPlate						
Application	Color (frame/bottom)	Feature	Well Number	Case of 14	Case of 56	Case of 68
Fluorescence Luminescence Absorbance	White/clear	Standard untreated	24-well			1450-602
		Tissue-culture treated	24-well	1450-603	1450-604	
	Black/clear	Tissue-culture treated	24-well	1450-605	1450-606	

Better microplates means better results

DELFLA					
Application	Color	Feature	Well number	Case of 10	Case of 60
Fluorescence	Yellow	Uncoated	96-well		AAAND-0001
		Streptavidin coated	96-well	AAAND-0005	
		Anti-rabbit coated	96-well	AAAND-0004	
	Clear	Streptavidin coated	8 x 12-well strips	4009-0010	
		Anti-rabbit coated	8 x 12-well strips	4008-0010	
		Anti-mouse coated	8 x 12-well strips	4007-0010	
	White	Streptavidin coated	384-well	CC11-H10	

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



revvity