

Research use only. Not for use in diagnostic procedures.

96-Well Assay Microplates: OptiPlate-96™ / CulturPlate-96™ / SpectraPlate-96™

Part numbers:

Plate Family	Well Density	Color	Treatment	Case of 50	Case of 160	Case of 200
OptiPlate™	96	White	Untreated	6055290		6055299
	96	Black	Untreated	6055270		6055279
	96	White	High Binding	6055240		6055249
	96	Black	High Binding	6055220		6055229
	96	White	Low Binding	6065420		
	96	Black	Low Binding	6065400		
CulturPlate™	96	White	TC-treated	6055680	6055688	6055689
	96	Black	TC-treated	6055660	6055668	6055669
SpectraPlate™	96	Clear	Medium Binding	6055640		6055649
	96	Clear	TC-Treated	6055650	6055658	
	96	Clear	High Binding	6055620		6055629
	96	Clear	Low Binding	6065360		

Material and treatment:

Black, white, or transparent polystyrene microplate with 96 flat-bottom wells. Plate is available either untreated, with high protein binding affinity, low protein binding affinity, or tissue-culture (TC) treated.

Packaging:

Case of	Number of Bags in a Case	Number of Plates in a Bag	Lids Included?
50*	2	25	No, except for TC-treated plates which do include lids
160	8	20	Yes
200	8	25	No

* Case of 50 TC-treated are supplied in 50 individual pouches with lids

Plate Specification Overview:

Parameter	Specification
Material	Polystyrene
Well Density	96-well
Well Shape	Round
Well-Bottom shape	Flat
Plate Length (mm)	127.76 ± 0.2
Plate Width (mm)	85.48 ± 0.2
Plate Height (mm)	14.35 ± 0.2
A1 row offset (mm)	11.24 ± 0.1
A1 column offset (mm)	14.38 ± 0.1
Well-to-well spacing (mm)	9.0
Flange/Skirt height (mm)	2.6 ± 0.05
Well Depth (mm)	10.8 ± 0.1
Well Diameter (mm)	7.15 ± 0.05
Maximum Well Volume	412 µL
Working Volume	80 µL - 350 µL

Technical Drawing:

