

MatTek Dishes Price List 2025



Part Number	Description	Unit Size	Unit Price
Chambered Cell Culture Slides			
CCS-2	2-well Chambered Cell Culture Slide	Case of 12	\$101.00
CCS-4	4-well Chambered Cell Culture Slide	Case of 12	\$101.00
CCS-8	8-well Chambered Cell Culture Slide	Case of 12	\$101.00

Coverslip Kits

CSGK/F	Glass Coverslip Kit, No. 1.5 Coverslip, 22 X 22 mm, 35 mm Falcon Dish	Case (500)	\$377.00
CSGK/F-F	Glass Coverslip Kit, No. 1.5 Coverslip, 22 X 22 mm, 35 mm Falcon Dish w/ foam for shipping	Case (500)	\$387.00
CSGK/M	Glass Coverslip Kit, No. 1.5 Coverslip, 22 X 22 mm, 35 mm MatTek Dish	Case (500)	\$377.00

Accessories

P DCF OS 30	Glass Bottom Dishes and Multi-Well Plate Coverslip Removal Fluid	Bottle (50 mL)	\$30.00
-------------	------------------------------------------------------------------	----------------	---------

35 mm Glass Bottom Dishes

P35G-0.170-14-C	35 mm Dish, No. 1.5 High Tolerance Coverslip 170 +/- 5 um, 14 mm Glass Diameter	Case (75)	\$356.00
P35G-0.170-14-C.S	35 mm Dish, No. 1.5 High Tolerance Coverslip 170 +/- 5 um, 14 mm Glass Diameter	Sleeve (5)	\$62.00
P35G-0-10-C	35 mm Dish, No. 0 Coverslip, 10 mm Glass Diameter, Uncoated	Case (200)	\$407.00
P35G-0-10-C.S	35 mm Dish, No. 0 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (10)	\$51.00
P35G-0-10-C-HA	35 mm Dish, No. 0 Coverslip, 10 mm Glass Diameter, High Adherence	Case (200)	\$458.00
P35G-0-10-C-HA.S	35 mm Dish, No. 0 Coverslip, 10 mm Glass Diameter, High Adherence	Sleeve (10)	\$51.00
P35G-0-14-C	35 mm Dish, No. 0 Coverslip, 14 mm Glass Diameter, Uncoated	Case (200)	\$407.00
P35G-0-14-C.S	35 mm Dish, No. 0 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (10)	\$51.00
P35G-0-14-C-HA	35 mm Dish, No. 0 Coverslip, 14 mm Glass Diameter, High Adherence	Case (200)	\$458.00
P35G-0-14-C-HA.S	36 mm Dish, No. 0 Coverslip, 14 mm Glass Diameter, High Adherence	Sleeve (10)	\$51.00
P35G-0-20-C	35 mm Dish, No. 0 Coverslip, 20 mm Glass Diameter, Uncoated	Case (200)	\$540.00
P35G-0-20-C.S	35 mm Dish, No. 0 Coverslip, 20 mm Glass Diameter, Uncoated	Sleeve (10)	\$62.00
P35G-0-7-C	35 mm Dish, No. 0 Coverslip, 7 mm Glass Diameter, Uncoated	Case (200)	\$509.00
P35G-0-7-C.S	35 mm Dish, No. 0 Coverslip, 7 mm Glass Diameter, Uncoated	Sleeve (10)	\$62.00
P35G-1.0-14-C	35 mm Dish, No. 1.0 Coverslip, 14 mm Glass Diameter, Uncoated	Case (200)	\$407.00
P35G-1.0-14-C.S	35 mm Dish, No. 1.0 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (10)	\$51.00
P35G-1.0-20-C	35 mm Dish, No. 1.0 Coverslip, 20 mm Glass Diameter, Uncoated	Case (200)	\$540.00
P35G-1.0-20-C.S	35 mm Dish, No. 1.0 Coverslip, 20 mm Glass Diameter, Uncoated	Sleeve (10)	\$62.00
P35G-1.5-10-C	35 mm Dish, No. 1.5 Coverslip, 10 mm Glass Diameter, Uncoated	Case (200)	\$387.00
P35G-1.5-10-C.S	35 mm Dish, No. 1.5 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (10)	\$51.00
P35G-1.5-10-C-HA	35 mm Dish, No. 1.5 Coverslip, 10 mm Glass Diameter, High Adherence	Case (200)	\$438.00
P35G-1.5-10-C-HA.S	35 mm Dish, No. 1.5 Coverslip, 10 mm Glass Diameter, High Adherence	Sleeve (10)	\$51.00
P35G-1.5-14-C	35 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, Uncoated	Case (200)	\$387.00
P35G-1.5-14-C.S	35 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (10)	\$51.00
P35G-1.5-14-C-GRD	35 mm Dish, No. 1.5 Gridded Coverslip, 14 mm Glass Diameter (Cells and grid in same focal plane)	Case (75)	\$540.00
P35G-1.5-14-C-GRD.S	35 mm Dish, No. 1.5 Gridded Coverslip, 14 mm Glass Diameter (Cells and grid in same focal plane)	Sleeve (5)	\$62.00
P35G-1.5-14-C-GRD-DOWN*	36 mm Dish, No. 1.5 Gridded Coverslip, 14 mm Glass Diameter (Cells and grid in different focal plane)	Case (200)	\$560.00
P35G-1.5-14-C-HA	35 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, High Adherence	Case (200)	\$438.00
P35G-1.5-14-C-HA.S	35 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, High Adherence	Sleeve (10)	\$51.00

P35G-1.5-20-C	35 mm Dish, No. 1.5 Coverslip, 20 mm Glass Diameter, Uncoated	Case (200)	\$509.00
P35G-1.5-20-C.S	35 mm Dish, No. 1.5 Coverslip, 20 mm Glass Diameter, Uncoated	Sleeve (10)	\$62.00
P35G-1.5-20-C-HA	35 mm Dish, No. 1.5 Coverslip, 20 mm Glass Diameter, High Adherence	Case (200)	\$560.00
P35G-1.5-20-C-HA.S	35 mm Dish, No. 1.5 Coverslip, 20 mm Glass Diameter, High Adherence	Sleeve (10)	\$62.00
P35G-1.5-7-C	35 mm Dish, No. 1.5 Coverslip, 7 mm Glass Diameter, Uncoated	Case (200)	\$509.00
P35G-1.5-7-C.S	35 mm Dish, No. 1.5 Coverslip, 7 mm Glass Diameter, Uncoated	Sleeve (10)	\$62.00
P35GC-0-10-C	35 mm Dish, No. 0 Coverslip, 10 mm Glass Diameter, Poly-D-Lysine Coated	Case (200)	\$509.00
P35GC-0-10-C.S	35 mm Dish, No. 0 Coverslip, 10 mm Glass Diameter, Poly-D-Lysine Coated	Sleeve (10)	\$62.00
P35GC-0-14-C	35 mm Dish, No. 0 Coverslip, 14 mm Glass Diameter, Poly-D-Lysine Coated	Case (200)	\$509.00
P35GC-0-14-C.S	35 mm Dish, No. 0 Coverslip, 14 mm Glass Diameter, Poly-D-Lysine Coated	Sleeve (10)	\$62.00
P35GC-1.0-14-C	35 mm Dish, No. 1.0 Coverslip, 14 mm Glass Diameter, Poly-D-Lysine Coated	Case (200)	\$509.00
P35GC-1.0-14-C.S	35 mm Dish, No. 1.0 Coverslip, 14 mm Glass Diameter, Poly-D-Lysine Coated	Sleeve (10)	\$62.00
P35GC-1.5-10-C	35 mm Dish, No. 1.5 Coverslip, 10 mm Glass Diameter, Poly-D-Lysine Coated	Case (200)	\$489.00
P35GC-1.5-10-C.S	35 mm Dish, No. 1.5 Coverslip, 10 mm Glass Diameter, Poly-D-Lysine Coated	Sleeve (10)	\$62.00
P35GC-1.5-14-C	35 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, Poly-D-Lysine Coated	Case (200)	\$489.00
P35GC-1.5-14-C.S	35 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, Poly-D-Lysine Coated	Sleeve (10)	\$62.00
P35GC-1.5-14-C-HA	35 mm Dish, No. 1.5 Coverslip, 20 mm Glass Diameter, High Adherence, Poly-D-Lysine Coated	Case (200)	\$540.00
P35GC-1.5-14-C-HA.S	36 mm Dish, No. 1.5 Coverslip, 20 mm Glass Diameter, High Adherence, Poly-D-Lysine Coated	Sleeve (10)	\$62.00
P35GCOL-0-14-C	35 mm Dish, No. 0 Coverslip, 14 mm Glass Diameter, Collagen Coated	Case (200)	\$489.00
P35GCOL-0-14-C.S	35 mm Dish, No. 0 Coverslip, 14 mm Glass Diameter, Collagen Coated	Sleeve (10)	\$62.00
P35GCOL-1.0-14-C	35 mm Dish, No. 1.0 Coverslip, 14 mm Glass Diameter, Collagen Coated	Case (200)	\$489.00
P35GCOL-1.0-14-C.S	35 mm Dish, No. 1.0 Coverslip, 14 mm Glass Diameter, Collagen Coated	Sleeve (10)	\$62.00
P35GCOL-1.5-10-C	35 mm Dish, No. 1.5 Coverslip, 10 mm Glass Diameter, Collagen Coated	Case (200)	\$489.00
P35GCOL-1.5-10-C.S	35 mm Dish, No. 1.5 Coverslip, 10 mm Glass Diameter, Collagen Coated	Sleeve (10)	\$62.00
P35GCOL-1.5-14-C	35 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, Collagen Coated	Case (200)	\$489.00
P35GCOL-1.5-14-C.S	35 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, Collagen Coated	Sleeve (10)	\$62.00

50 mm Glass Bottom Dishes

P50G-0-14-F	50 mm Dish, No. 0 Coverslip, 14 mm Glass Diameter, Uncoated	Case (200)	\$438.00
P50G-0-14-F.S	50 mm Dish, No. 0 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (10)	\$51.00
P50G-0-30-F	50 mm Dish, No. 0 Coverslip, 30 mm Glass Diameter, Uncoated	Case (200)	\$591.00
P50G-0-30-F.S	50 mm Dish, No. 0 Coverslip, 30 mm Glass Diameter, Uncoated	Sleeve (10)	\$71.00
P50G-1.5-14-F	50 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, Uncoated	Case (200)	\$407.00
P50G-1.5-14-F.S	50 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (10)	\$51.00
P50G-1.5-30-F	50 mm Dish, No. 1.5 Coverslip, 30 mm Glass Diameter, Uncoated	Case (200)	\$591.00
P50G-1.5-30-F.S	50 mm Dish, No. 1.5 Coverslip, 30 mm Glass Diameter, Uncoated	Sleeve (10)	\$71.00
P50GC-1.5-14-F	50 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, Poly-D-Lysine Coated	Case (200)	\$509.00
P50GC-1.5-14-F.S	50 mm Dish, No. 1.5 Coverslip, 14 mm Glass Diameter, Poly-D-Lysine Coated	Sleeve (10)	\$62.00

60 mm Glass Bottom Dishes

P60G-1.5-20-F	60 mm Dish, No. 1.5 Coverslip, 20 mm Glass Diameter, Uncoated	Case (100)	\$591.00
P60G-1.5-20-F.S	60 mm Dish, No. 1.5 Coverslip, 20 mm Glass Diameter, Uncoated	Sleeve (5)	\$82.00
P60G-1.5-30-F	60 mm Dish, No. 1.5 Coverslip, 30 mm Glass Diameter, Uncoated	Case (100)	\$591.00
P60G-1.5-30-F.S	60 mm Dish, No. 1.5 Coverslip, 30 mm Glass Diameter, Uncoated	Sleeve (5)	\$82.00

100 mm Glass Bottom Dishes

P100G-1.5-30-F	100 mm Dish, No. 1.5 Coverslip, 30 mm Glass Diameter, Uncoated	Case (75)	\$540.00
P100G-1.5-30-F.S	100 mm Dish, No. 1.5 Coverslip, 30 mm Glass Diameter, Uncoated	Sleeve (5)	\$82.00

6-well Multiwell Plates

P06G-0-10-F	6-Well Plate, No. 0 Coverslip, 10 mm Glass Diameter, Uncoated	Case (10)	\$234.00
P06G-0-10-F.S	6-Well Plate, No. 0 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (1)	\$41.00
P06G-0-14-F	6-Well Plate, No. 0 Coverslip, 14 mm Glass Diameter, Uncoated	Case (10)	\$234.00
P06G-0-14-F.S	6-Well Plate, No. 0 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (1)	\$41.00
P06G-0-20-F	6-Well Plate, No. 0 Coverslip, 20 mm Glass Diameter, Uncoated	Case (10)	\$234.00
P06G-0-20-F.S	6-Well Plate, No. 0 Coverslip, 20 mm Glass Diameter, Uncoated	Sleeve (1)	\$41.00
P06G-1.0-10-F	6-Well Plate, No. 1.0 Coverslip, 10 mm Glass Diameter, Uncoated	Case (10)	\$234.00
P06G-1.0-10-F.S	6-Well Plate, No. 1.0 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (1)	\$41.00
P06G-1.0-14-F	6-Well Plate, No. 1.0 Coverslip, 14 mm Glass Diameter, Uncoated	Case (10)	\$234.00
P06G-1.0-14-F.S	6-Well Plate, No. 1.0 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (1)	\$41.00
P06G-1.0-20-F	6-Well Plate, No. 1.0 Coverslip, 20 mm Glass Diameter, Uncoated	Case (10)	\$234.00
P06G-1.0-20-F.S	6-Well Plate, No. 1.0 Coverslip, 20 mm Glass Diameter, Uncoated	Sleeve (1)	\$41.00
P06G-1.5-10-F	6-Well Plate, No. 1.5 Coverslip, 10 mm Glass Diameter, Uncoated	Case (10)	\$234.00
P06G-1.5-10-F.S	6-Well Plate, No. 1.5 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (1)	\$41.00
P06G-1.5-14-F	6-Well Plate, No. 1.5 Coverslip, 14 mm Glass Diameter, Uncoated	Case (10)	\$234.00
P06G-1.5-14-F.S	6-Well Plate, No. 1.5 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (1)	\$41.00
P06G-1.5-20-F	6-Well Plate, No. 1.5 Coverslip, 20 mm Glass Diameter, Uncoated	Case (10)	\$234.00
P06G-1.5-20-F.S	6-Well Plate, No. 1.5 Coverslip, 20 mm Glass Diameter, Uncoated	Sleeve (1)	\$41.00

12-well Multiwell Plates

P12G-0-10-F	12-Well Plate, No. 0 Coverslip, 10 mm Glass Diameter, Uncoated	Case (10)	\$336.00
P12G-0-10-F.S	12-Well Plate, No. 0 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (1)	\$51.00
P12G-0-14-F	12-Well Plate, No. 0 Coverslip, 14 mm Glass Diameter, Uncoated	Case (10)	\$336.00
P12G-0-14-F.S	12-Well Plate, No. 0 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (1)	\$51.00
P12G-1.0-10-F	12-Well Plate, No. 1.0 Coverslip, 10 mm Glass Diameter, Uncoated	Case (10)	\$336.00
P12G-1.0-10-F.S	12-Well Plate, No. 1.0 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (1)	\$51.00
P12G-1.0-14-F	12-Well Plate, No. 1.0 Coverslip, 14 mm Glass Diameter, Uncoated	Case (10)	\$336.00
P12G-1.0-14-F.S	12-Well Plate, No. 1.0 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (1)	\$51.00
P12G-1.5-10-F	12-Well Plate, No. 1.5 Coverslip, 10 mm Glass Diameter, Uncoated	Case (10)	\$336.00
P12G-1.5-10-F.S	12-Well Plate, No. 1.5 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (1)	\$51.00
P12G-1.5-14-F	12-Well Plate, No. 1.5 Coverslip, 14 mm Glass Diameter, Uncoated	Case (10)	\$336.00
P12G-1.5-14-F.S	12-Well Plate, No. 1.5 Coverslip, 14 mm Glass Diameter, Uncoated	Sleeve (1)	\$51.00

24-well Multiwell Plates

P24G-0-10-F	24-Well Plate, No. 0 Coverslip, 10 mm Glass Diameter, Uncoated	Case (10)	\$387.00
P24G-0-10-F.S	24-Well Plate, No. 0 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (1)	\$62.00
P24G-0-13-F	24-Well Plate, No. 0 Coverslip, 13 mm Glass Diameter, Uncoated	Case (10)	\$387.00
P24G-0-13-F.S	24-Well Plate, No. 0 Coverslip, 13 mm Glass Diameter, Uncoated	Sleeve (1)	\$62.00
P24G-1.0-10-F	24-Well Plate, No. 1.0 Coverslip, 10 mm Glass Diameter, Uncoated	Case (10)	\$387.00
P24G-1.0-10-F.S	24-Well Plate, No. 1.0 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (1)	\$62.00
P24G-1.0-13-F	24-Well Plate, No. 1.0 Coverslip, 13 mm Glass Diameter, Uncoated	Case (10)	\$387.00
P24G-1.0-13-F.S	24-Well Plate, No. 1.0 Coverslip, 13 mm Glass Diameter, Uncoated	Sleeve (1)	\$62.00

P24G-1.5-10-F	24-Well Plate, No. 1.5 Coverslip, 10 mm Glass Diameter, Uncoated	Case (10)	\$387.00
P24G-1.5-10-F.S	24-Well Plate, No. 1.5 Coverslip, 10 mm Glass Diameter, Uncoated	Sleeve (1)	\$62.00
P24G-1.5-13-F	24-Well Plate, No. 1.5 Coverslip, 13 mm Glass Diameter, Uncoated	Case (10)	\$387.00
P24G-1.5-13-F.S	24-Well Plate, No. 1.5 Coverslip, 13 mm Glass Diameter, Uncoated	Sleeve (1)	\$62.00
P24GTOP-1.5-F	Glass Coverslip Lid For DIC In P06, P12 and P24 Multi-Well Products	Case (10)	\$458.00
P24GTOP-1.5-F.S	Glass Coverslip Lid For DIC In P06, P12 and P24 Multi-Well Products	Sleeve (1)	\$72.00

48-well Multiwell Plates

P48G-1.5-6-F	48-Well Plate, No. 1.5 Coverslip, 6 mm Glass Diameter, Uncoated	Case (10)	\$407.00
P48G-1.5-6-F.S	48-Well Plate, No. 1.5 Coverslip, 6 mm Glass Diameter, Uncoated	Sleeve (1)	\$72.00

96-well Multiwell Plates

P96G-1.5-5-F	96-Well Plate, No. 1.5 Coverslip, 5 mm Glass Diameter, Uncoated	Case (5)	\$234.00
P96G-1.5-5-F.S	96-Well Plate, No. 1.5 Coverslip, 5 mm Glass Diameter, Uncoated	Sleeve (1)	\$72.00
P96GC-1.5-5-F	96-Well Plate, No. 1.5 Coverslip, 5 mm Glass Diameter, Poly-D-Lysine Coated	Case (5)	\$305.00
P96GC-1.5-5-F.S	96-Well Plate, No. 1.5 Coverslip, 5 mm Glass Diameter, Poly-D-Lysine Coated	Sleeve (1)	\$92.00
PBK96G-1.5-5-F	96-Well Black Plate, No. 1.5 Coverslip, 5 mm Glass Diameter, Uncoated	Case (5)	\$254.00
PBK96G-1.5-5-F.S	96-Well Black Plate, No. 1.5 Coverslip, 5 mm Glass Diameter, Uncoated	Sleeve (1)	\$72.00

384-well Multiwell Plates

P384G-1.5-10872-C	Single Specimen Plate, No. 1.5 Coverslip, 98 mm x 67 mm Viewing Area, Uncoated	Case (5)	\$285.00
P384G-1.5-10872-C.S	Single Specimen Plate, No. 1.5 Coverslip, 98 mm x 67 mm Viewing Area, Uncoated	Sleeve (1)	\$72.00
PBK384G-1.5-C	384-Well Black Plate, No. 1.5 Coverslip, Uncoated	Case (5)	\$305.00
PBK384G-1.5-C.S	384-Well Black Plate, No. 1.5 Coverslip, Uncoated	Sleeve (1)	\$82.00

Cell Culture Inserts

CCI24-PC-0.4	PermaCell Cell Culture Insert, 24-well, PC, 0.4 μ m	Inserts (24)	\$123.00
CCI24-PC-8	PermaCell Cell Culture Insert, 24-well, PC, 8 μ m	Inserts (24)	\$123.00
CCI24-PET-0.4	PermaCell Cell Culture Insert, 24-well, PET, 0.4 μ m	Inserts (24)	\$123.00
CCI24-PTFE-0.4	PermaCell Cell Culture Insert, 24-well, PTFE, 0.4 μ m	Inserts (24)	\$133.00
CCI12-HANGTOP	12-Well Plate with Specialized Lid for Air-lifting CCI24 PermaCell Cell Culture Inserts	1 Plate	\$51.00
CCI24-HANGTOP	24-Well Plate with Specialized Lid for Air-lifting CCI24 PermaCell Cell Culture Inserts	1 Plate	\$51.00

Coverslips

PCS-0-10872	Coverslips, No. 0, German Glass, 108 X 72 mm Rectangle, Non-Sterile	Case (10)	\$86.00
PCS-0-15	Coverslips, No. 0, German Glass, 15 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-0-17	Coverslips, No. 0, German Glass, 17 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-0-18	Coverslips, No. 0, German Glass, 18 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-0-1818	Coverslips, No. 0, German Glass, 18 x 18 mm Square, Non-Sterile	Case (100)	\$61.00
PCS-0-25	Coverslips, No. 0, German Glass, 25 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-0-35	Coverslips, No. 0, German Glass, 35 mm Round, Non-Sterile	Case (100)	\$71.00
PCS-1.0-15	Coverslips, No. 1.0, German Glass, 15 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-1.0-17	Coverslips, No. 1.0, German Glass, 17 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-1.0-18	Coverslips, No. 1.0, German Glass, 18 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-1.0-1818	Coverslips, No. 1.0, German Glass, 18 x 18 mm Square, Non-Sterile	Case (100)	\$61.00
PCS-1.0-25	Coverslips, No. 1.0, German Glass, 25 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-1.5-10	Coverslips, No. 1.5, German Glass, 10 mm Round, Non-Sterile	Case (100)	\$61.00

PCS-1.5-10872	Coverslips, No. 1.5, German Glass, 108 X 72 mm Rectangle, Non-Sterile	Case (10)	\$86.00
PCS-1.5-15	Coverslips, No. 1.5, German Glass, 15 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-1.5-17	Coverslips, No. 1.5, German Glass, 17 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-1.5-18	Coverslips, No. 1.5, German Glass, 18 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-1.5-1818	Coverslips, No. 1.5, German Glass, 18 x 18 mm Square, Non-Sterile	Case (100)	\$61.00
PCS-1.5-1818-GRD	Coverslips, No. 1.5, Gridded German Glass, 18 x 18 mm Square, Non-Sterile	Case (25)	\$127.00
PCS-1.5-2222	Coverslips, No. 1.5, German Glass, 22 x 22 mm Square, Non-Sterile	Case (100)	\$61.00
PCS-1.5-25	Coverslips, No. 1.5, German Glass, 25 mm Round, Non-Sterile	Case (100)	\$61.00
PCS-1.5-35	Coverslips, No. 1.5, German Glass, 35 mm Round, Non-Sterile	Case (100)	\$71.00
PCS-1.5-5024	Coverslips, No. 1.5, German Glass, 50 x 24 mm Rectangle, Non-Sterile	Case (100)	\$71.00
PCS-170-1818	Coverslips, No. 1.5, High Tolerance 170 +/- 5 um, German Glass, 18 x 18 mm Square	Case (100)	\$81.00
PCS-1.5-11677*	Coverslips, No. 1.5, German Glass , Non-Sterile, 116 x 77 mm Rectangle, Non-sterile	Case (1000)	\$2,648.00

Organ-on-a-Chip Systems by AIM Biotech

CBLK	Cooling Block	Each (1)	\$220.00
DAX-1/ IDTX3	idenTx3 3D Cell Culture Chip, Pack of 25 Chips	Each (25)	\$484.00
HOL-2	idenTx Holder, sterile AIM Chip idenTx Holder	Each (10)	\$163.35
IDTX40	idenTx 40 3D Microfluidics Plate, 40 individual experiments, Box of 5	Box (5)	\$1,925.00
IDTX9	idenTx 9 3D Microfluidics Plate, 9 individual experiments, Box of 8	Box (8)	\$595.32
LUC-1	Luer Connectors with 1 sheet of 20 pre-cut Inlet Seals	Each (36)	\$121.00
OGNXI	organiX Insert Individually Packaged, Sterile organiX Inserts, pack of 24	Inserts (24)	\$361.90
OGNXP-1EA	organiX Plate 24-well, 3D Tissues Culture Plate, Pack of 1	Pack (1)	\$350.90
OGNXP-3EA	organiX Plate 24-well, 3D Tissues Culture Plate, Pack of 3	Pack (3)	\$950.40

Spheroid Culture Plates

F202001	BIOFLOAT 96 well Plate	Each (1)	\$23.00
F202003	BIOFLOAT 96 well Plate 4 Pack	4 Pack	\$82.00
F224384	BIOFLOAT 384 well Plate 4PCS	4 Pack	\$215.00