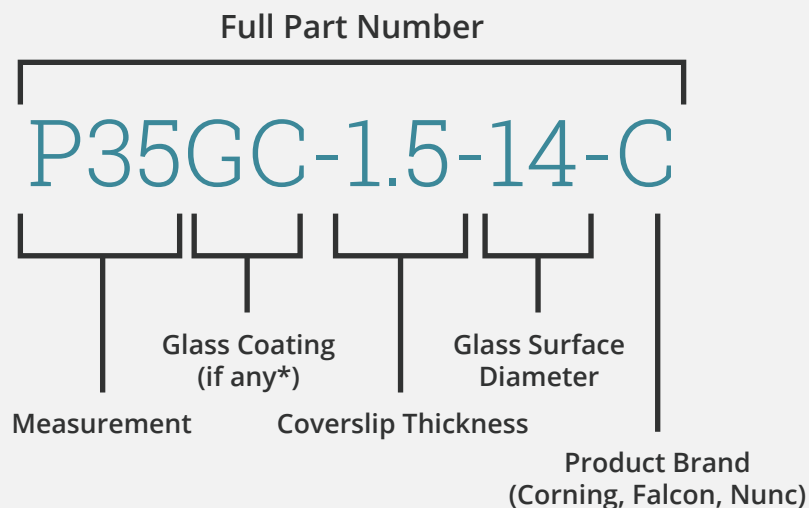


Free sample sleeves are available for the following products

ANATOMY OF A MATTEK PART NUMBER



* Coatings are denoted with C for Poly-D-Lysine and COL for Collagen

PRODUCT	DISH DIAMETER/ PLATE TYPE	COATING	GLASS THICKNESS	GLASS DIAMETER	SLEEVE QTY
---------	------------------------------	---------	--------------------	-------------------	---------------

Coverslip Kits

CSGK/C-HT	35 mm	Uncoated	No. 1.5	18 mm	10
CSGK/F	35 mm	Uncoated	No. 1.5	22 mm	20

Glass Bottom Multi-well Plates

P06G-0-20-F	6-well	Uncoated	No. 0	20 mm	1
P06G-1.0-20-F	6-well	Uncoated	No. 1.0	20 mm	1
P06G-1.5-14-F	6-well	Uncoated	No. 1.5	14 mm	1
P06G-1.5-20-F	6-well	Uncoated	No. 1.5	20 mm	1
P12G-0-14-F	12-well	Uncoated	No. 0	14 mm	1
P12G-1.0-14-F	12-well	Uncoated	No. 1.0	14 mm	1
P12G-1.5-10-F	12-well	Uncoated	No. 1.5	10 mm	1
P12G-1.5-14-F	12-well	Uncoated	No. 1.5	14 mm	1
P24G-0-13-F	24-well	Uncoated	No. 0	13 mm	1
P24G-1.0-13-F	24-well	Uncoated	No. 1.0	13 mm	1
P24G-1.5-13-F	24-well	Uncoated	No. 1.5	13 mm	1
P96G-1.5-5-F	96-well	Uncoated	No. 1.5	5 mm	1
P96GC-1.5-5-F	96-well	Uncoated grid	No. 1.5	5 mm	1

PRODUCT	DISH DIAMETER/ PLATE TYPE	COATING	GLASS THICKNESS	GLASS DIAMETER	SLEEVE QTY
Glass bottom dishes					
P35G-0.170-14-C	35 mm	Uncoated	No. 1.5	14 mm	5
P35G-0-10-C	35 mm	Uncoated	No. 0	10 mm	10
P35G-0-14-C	35 mm	Uncoated	No. 0	14 mm	10
P35G-0-20-C	35 mm	Uncoated	No. 0	20 mm	10
P35G-1.0-14-C	35 mm	Uncoated	No. 1.0	14 mm	10
P35G-1.0-20-C	35 mm	Uncoated	No. 1.0	20 mm	10
P35G-1.5-10-C	35 mm	Uncoated	No. 1.5	10 mm	10
P35G-1.5-14-C	35 mm	Uncoated	No. 1.5	14 mm	10
P35G-1.5-20-C	35 mm	Uncoated	No. 1.5	20 mm	10
P35G-2-14-CGRD	35 mm	Uncoated grid	No. 2.0	14 mm	5
P35GC-0-10-C	35 mm	Poly-D-Lysine	No. 0	10 mm	10
P35GC-0-14-C	35 mm	Poly-D-Lysine	No. 0	14 mm	10
P35GC-1.0-14-C	35 mm	Poly-D-Lysine	No. 1.0	14 mm	10
P35GC-1.5-10-C	35 mm	Poly-D-Lysine	No. 1.5	10 mm	10
P35GC-1.5-14-C	35 mm	Poly-D-Lysine	No. 1.5	14 mm	10
P35GCOL-0-10-C	35 mm	Collagen	No. 0	10 mm	10
P35GCOL-0-14-C	35 mm	Collagen	No. 0	14 mm	10
P35GCOL-1.0-14-C	35 mm	Collagen	No. 1.0	14 mm	10
P35GCOL-1.5-14-C	35 mm	Collagen	No. 1.5	14 mm	10
P50G-0-14-F	50 mm	Uncoated	No. 0	14 mm	10
P50G-1.5-14-F	50 mm	Uncoated	No. 1.5	14 mm	10
P50G-2-14-FGRD	50 mm	Uncoated grid	No. 2.0	14 mm	5
P50GC-14-F	50 mm	Poly-D-Lysine	No. 1.5	14 mm	10

Part Number Key

Dish diameter - Glass Bottom Dishes (35 or 50 mm)

P35 35 mm

Coating

G Uncoated

GC Poly-D-Lysine Coated

GCOL Collagen Coating

Coverslip Thickness

No. 0 0.085 - 0.13 mm

No.1.0 0.13 - 0.16 mm

No. 1.5 0.16 - 0.19 mm

No. 2.0 0.19 - 0.23 mm

Glass Diameter

10 Glass Diameter (10,14, 20 or 30 mm)

Other

GRD Gridded Coverslip

Number of wells - Multi-Well Culture Plates (6, 12, 24, or 96)

P06 6-well

Coating

G Uncoated

GC Poly-D-Lysine Coated (special order)

GCOL Collagen Coating (special order)

Coverslip Thickness

No. 0 0.085 - 0.13 mm

No.1.0 0.13 - 0.16 mm

No. 1.5 0.16 - 0.19 mm

No. 2.0 0.19 - 0.23 mm

Glass Diameter

10 Glass Diameter (10,13, 14 or 20 mm)