

# Crystallization Plates

## Overview

Cat.- No.	Plate name	Manufacturer	Max. reservoir volume	# protein wells per reservoir	Protein drop volume	Material / Treatment
<b>24 Well Plates</b>						
CPL-101(S/L)	Linbro Plate	MP Biomedicals	3500 µl	-	-	Polystyrene
CPL-119	ComboPlate™	Greiner Bio-One	3300 µl	-	-	Polystyrene
CPL-131	ComboPlate™, pregreased	Greiner Bio-One	3300 µl	-	-	Polystyrene
CPL-130	Crystalgen SuperClear™ Plates	Crystalgen	3500 µl	-	-	Polystyrene
CPL-132	Crystalgen SuperClear™ Plates, pregreased	Crystalgen	3500 µl	-	-	Polystyrene
CPL-137(S/L)	CrystalPlate 24	Axygen Biosciences	3500 µl			Polystyrene
CPL-102	CombiClover™ Plate	Emerald BioSystems	1500 µl	4	50 µl	Polypropylene
CPL-103	CompactClover™ Plate	Emerald BioSystems	250 µl	4	50 µl	Polypropylene
CPL-105	CombiClover™ Junior Plate	Emerald BioSystems	500 µl	4	10 µl	Polypropylene
<b>48 Well Plates</b>						
CPL-138(S/L)	CrystalPlate 48	Axygen Biosciences	1500 µl	-	-	Polystyrene
<b>96 Well Plates</b>						
CPL-118(S/L)	CrystalQuick™ - 609101	Greiner Bio-One	320 µl	3	4,1 µl	Polystyrene
CPL-107(S/L)	CrystalQuick™ - 609120	Greiner Bio-One	320 µl	3	1,9 µl	Polystyrene
CPL-108(S/L)	CrystalQuick™ - 609171	Greiner Bio-One	140 µl	1	3,9 µl	Polystyrene
CPL-120(S/L)	CrystalQuick™ - 609801	Greiner Bio-One	320 µl	3	4,1 µl	LBR
CPL-121(S/L)	CrystalQuick™ - 609820	Greiner Bio-One	320 µl	3	1,9 µl	LBR
CPL-122(S/L)	CrystalQuick™ - 609871	Greiner Bio-One	140 µl	1	3,9 µl	LBR
CPL-123(S/L)	CrystalQuick™ - 609130	Greiner Bio-One	320 µl	3	4,1 µl	Hydrophobic Surface
CPL-124(S/L)	CrystalQuick™ - 609830	Greiner Bio-One	320 µl	3	4,1 µl	LBR & Hydrophobic S.
CPL-125(S/L)	CrystalQuick™ - 609180	Greiner Bio-One	140 µl	1	3,9 µl	Hydrophobic Surface
CPL-126(S/L)	Corning® Plate 3773	Corning Life Science	210 µl	1	10 µl	Advanced Polymer
CPL-127(S/L)	Corning® Plate 3785	Corning Life Science	210 µl	1	7 µl	Advanced Polymer
CPL-109(S/L)	Corning® Plate 3556	Corning Life Science	100 µl	1	4 µl	Hydrophobic COC
CPL-110(S/L)	Corning® Plate 3551	Corning Life Science	100 µl	1	4 µl	Hydrophobic COC
CPL-111(S/L)	Corning® Plate 3552	Corning Life Science	100 µl	3	2 µl	PZero Polymer
CPL-112(S/L)	Corning® Plate 3553	Corning Life Science	100 µl	3	2 µl	PZero Polymer
CPL-113(S/L)	Corning® Plate 3554	Corning Life Science	100 µl	3	2 µl	PZero Polymer
CPL-114(S/L)	Corning® Plate 3555	Corning Life Science	100 µl	3	2 µl	PZero Polymer
CPL-115(S/L)	Corning® Plate 3557	Corning Life Science	100 µl	5	1 µl	PZero Polymer
CPL-116(S/L)	Corning® Plate 3550	Corning Life Science	100 µl	5	1 µl	PZero Polymer
CPL-133	CrystalClear 'Duo'	Douglas Instruments	100 µl	2	2 µl	Polystyrene
CPL-134	CrystalClear 'P'	Douglas Instruments	100 µl	1	2 µl	Polystyrene
CPL-135(S/L)	AxyGem™ Crystallography Plate	Axygen Biosciences	250 µl	1	4,3 µl	Polystyrene
CPL-139	Innovaplate™ SD-2	Innovadyne	80 µl		5 µl	New Patented Polymer
CPL-106	CompactClover™ Junior Plate	Emerald BioSystems	300 µl	1	6 µl	Polypropylene
CPL-141	Clover 384™ Plate	Emerald BioSystems	70 µl	4	1 µl	Polystyrene
CPL-142	Clover 384™ Plate, LBR	Emerald BioSystems	70 µl	4	1 µl	LBR
<b>Microbatch</b>						
CPL-117(S/L)	Imp@ct™ Plate - 673170	Greiner Bio-One			8 µl	Polystyrene
CPL-144(S/L)	Imp@ct™ Plate – 673096, black, 96 well	Greiner Bio-One	-	-	8 µl	Polystyrene
CPL-145(S/L)	Imp@ct™ Plate – 790801, LBR, 1536 well	Greiner Bio-One	-	-	10,1 µl	Polystyrene
CPL-146	Terasaki – 653102, 60 well	Greiner Bio-One	-	-	11,45 µl	Polystyrene
CPL-147	Terasaki – 654102, 72 well	Greiner Bio-One	-	-	11,45 µl	Polystyrene
CPL-128(S/L)	Vapor Batch Plate, hydrophobic	Douglas Instruments	-	-	9,3 µl	Hydrophobic Surface
CPL-129(S/L)	Vapor Batch Plate, hydrophilic	Douglas Instruments	-	-	9,3 µl	Polystyrene