

Plate Name	Cat.-No.	Manufacturer	Wells/ Plate	Reservoir Well Volume	Protein Wells/ Reservoir	Protein Well Volume	Material
Hanging Drop							
Crystalgen SuperClear™ Plate	CPL-130	Crystalgen	24	3.500 µl	–	–	Polystyrene
Crystalgen SuperClear™ Plate, pregreased	CPL-132	Crystalgen	24	3.500 µl	–	–	Polystyrene
Linbro Plate Space Saver	CPL-166	MP Biomedicals	24	3.500 µl	–	–	Polystyrene
HDP Hanging Drop Crystallization Plate	CPL-151	Swissci	96	200 µl	–	5 µl	UVP
In-situ-1 Crystallization Plate	CPL-160	MiTeGen	96	40 µl	1	droplets 50 nL - 2 µl	COC
Sitting Drop							
MRC Maxi Plate	CPL-154	Swissci	48	200 µl	1	10 µl	UVP
MRC Maxi Plate	CPL-155	Swissci	48	200 µl	1	10 µl	Polystyrene
In-situ-1 Crystallization Plate	CPL-160	MiTeGen	96	40 µl	1	droplets 50 nL - 2 µl	COC
CrystalDirect™ Plate	CPL-167	MiTeGen	96	50 µl	1	droplets up to 1 µl	COC
MRC 2 Well Crystallization Plate	CPL-139	Swissci	96	85 µl	2	5 µl	UVP
MRC 2 Well Crystallization Plate	CPL-153	Swissci	96	85 µl	2	5 µl	Polystyrene
MRC 2 Well Crystallization Plate	CPL-163	Swissci	96	85 µl	2	5 µl	UVXPO
3 Well Crystallization Plate	CPL-149	Swissci	96	30 µl	3	5 µl	UVP
3 Well Crystallization Plate	CPL-150	Swissci	96	30 µl	3	5 µl	Polystyrene
3 Well Crystallization Plate	CPL-168	Swissci	96	30 µl	3	5 µl	UVXPO
3 Well Crystallization Plate - Low Profile	CPL-164	Swissci	96	30 µl	3	5 µl	UVP
3 Well Crystallization Plate - Low Profile	CPL-165	Swissci	96	30 µl	3	5 µl	Polystyrene
3 Well Crystallization Plate - Low Profile	CPL-169	Swissci	96	30 µl	3	5 µl	UVXPO
Lipidic Cubic Phase Plates							
IMISX™ Plate Kit	CPL-162	MiTeGen	96	600 nL + 100 nL LCP	–	–	COC, Glass
LCP Glass Sandwich Set	CPL-159	Marienfeld	96	1 µl + 50 nL LCP	–	–	Glass
Lipidic Cubic Phase Screening Kit	CPL-156	Swissci	96	1 µl + 50 nL LCP	–	–	UVP
Microbatch Plates							
MRC Under Oil Crystallization Plate	CPL-152	Swissci	96	20 µl	–	–	UVP
Dialysis							
Diaplate Kit	X-W82010	Swissci	96	–	–	3.2 µl	Polystyrene

Product name	Cat.-No.	Manufacturer
Sealing Tapes		
HD Clear™ Sealing Tape - 3 inches wide	CTP-101M	Duck
HD Clear™ Sealing Tape - 1.88 inches wide	CTP-100	Duck
Greiner VIEWseal™	CTP-102	Greiner Bio-One
UVXPO Sealing Sheets	CTP-106	Swissci
96 Well Masterblocks & Cap Mats		
Greiner 780271 – Masterblock 2 ml	CMB-102	Greiner Bio-One
AxyMat Silicone Sealing Mat	CCM-103	Axygen

