

Competitor Cross Guide: Mettler XPR vs. Cubis® II



Model	Readability (mg)	Weighing capacity (g)	Linearity (mg)*	Corner load (mg)*	Typical Repeatability (mg) 5%/max*	Typical Stabilization time (s)*	Minimum Weight (mg)*	Pan size (mm)	Chamber size DWH or DH (mm)
Ultramicro Balances									
XPR2U	0.0001	2.1	0.001	0.0015	0.00015	<10	0.3	16	77 x 55
XPR3DUE	0.001 0.01	1.2 3.2	0.006		0.0008	6	1.6	32 x 32	
MCA2.7S-2S00-M	0.0001	2.1	0.0007	0.0005	0.00015	7	0.3	20	80 x 67
Micro Balances									
XPR6UD5	0.0005	6.1	0.003	0.002	0.0003	<8	0.6	27	77 x 55
XPR2	0.001	2.1	0.003	0.002	0.0005	<8	1	27	77 x 55
MCA6.6S-2S00-M-QP99	0.001	6.1	0.003	0.003	0.0005	5	0.82	30	80 x 67
XPR10	0.001	10.1	0.003	0.003	0.0004	<8	0.82	27	77 x 55
XPR16DUE	0.001 0.01	3.2 12	0.006		0.0012	6	2.4	32 x 32	
MCA10.6S-2S00-M-QP99	0.001	10.1	0.003	0.003	0.0005	5	0.82	30	80 x 67
High Capacity Micro Balances									
XPR36	0.001	32	0.005	0.006	0.0007	3.5	1.4	40 x 40	85 x 69 x 114
MCA36S-3S00-D-QP99-MDS	0.001	32	0.005	0.006	0.0007	3.5	1.4	50 90	185 x 159 x 234
XPR56	0.001	52	0.006	0.01	0.0007	3.5	1.4	40 x 40	85 x 69 x 114
MCA66S-3S00-D-QP99-MDS	0.001	61	0.005	0.01	0.0007	3.5	1.4	50 90	185 x 159 x 234
XPR26DR	0.001 0.01	5.1 22	0.006	0.008	0.001	3.5	2	-	-
XPR36DR	0.001 0.01	8.1 32	0.008	0.008	0.001	2.5	2	40 x 40	85 x 69 x 114
MCA36P-3S00-D-QP99-MDS	0.001 0.01	10 32	0.006	0.008	0.0007	3.5	1.4	50 90	185 x 159 x 234
XPR56DR	0.001 0.01	11 52	0.01	0.012	0.001	2.5	2	40 x 40	85 x 69 x 114
MCA66P-3S00-D-QP99-MDS	0.001 0.01	12 61	0.008	0.012	0.002	3.5	1.4	50 90	185 x 159 x 234
XPR106DUH	0.002 0.005	41 120	0.03	0.04	0.003	4	6	64 x 56	160 x 137 x 235
MCA116S-3S00-D-QP99-MDS	0.002	111	0.02	0.02	0.0025 0.01	3.5	5	50 90	185 x 159 x 234

Model	Readability (mg)	Weighing capacity (g)	Linearity (mg)*	Corner load (mg)*	Typical Repeatability (mg) 5%/max*	Typical Stabilization time (s)*	Minimum Weight (mg)*	Pan size (mm)	Chamber size DWH or DH (mm)
Semi-Micro Balances									
XPR226DR	0.01	120	0.03	0.04	0.005	2	10	78 x 73	160 x 137 x 235
MCA226S-3S00-D-QP99-QAPP1002	0.005	220	0.03	0.04	0.004 0.015	1.5	8	50	185 x 159 x 234
XPR105	0.001	120	0.03	0.04	0.007	2	14	78 x 73	160 x 137 x 235
MCA125S-3S00-D-QP99-QAPP1002	0.001	120	0.03	0.04	0.007 0.015	1.5	13	90	185 x 137 x 234
XPR205	0.001	220	0.03	0.06	0.007	2	14	78 x 73	160 x 137 x 235
MCA225S-3S00-D-QP99-QAPP1002	0.001	220	0.03	0.05	0.007 0.015	1.5	13	90	160 x 137 x 235
XPR105DR	0.001 0.01	41 121	0.05	0.06	0.007	1.5	14	78 x 73	160 x 137 x 235
XPR105DUE	0.01 0.1	42 120	0.06		0.01	1.5	20	78 x 73	
MCA125P-2S00-A-QP99	0.001 0.01	60 120	0.065	0.1	0.01 0.02	2	13	85 x 85	192 x 154 x 250
XPR205DR	0.001 0.01	81 220	0.05	0.08	0.007	1.5	14	78 x 73	160 x 137 x 235
XPR205DU	0.001 0.01	81 220	0.06	0.1	0.01	1.5	20	78 x 73	160 x 137 x 235
XPR225DR	0.001 0.01	121 220	0.03	0.08	0.007	1.5	14	78 x 73	160 x 137 x 235
XPR205DUE	0.001 0.01	81 220	0.06		0.01	1.5	20	78 x 73	
XPR225DUE	0.001 0.01		0.06		0.01	1.5	20	78 x 73	
XPR305DR	0.001 0.01	121 320			0.02	1.5	40	78 x 73	160 x 137 x 235
XPR225DU	0.001 0.01	121 220	0.05	0.1	0.007	1.5	14	78 x 73	160 x 137 x 235
XPR205D5	0.05	220	0.06	0.08	0.02	1.5	41	78 x 73	160 x 137 x 235
MCA225P-3S00-D-QP99-QAPP1002	0.001 0.01	120 220	0.103	0.06	0.007 0.02	1.5	13	90	185 x 159 x 234
Analytical Balances									
XPR104DUE	0.1 1	81 120	0.06		0.04	1.5	82	78 x 73	
XPR204E	0.1	220	0.06		0.04	1.5	82	78 x 73	
XPR204DUE	0.1 1	121 220	0.06		0.04	1.5	82	78 x 73	
XPR204	0.1	220	0.06	0.1	0.04	1.5	82	78 x 73	160 x 137 x 235
MCA224S-2S00-A-QP99	0.1	220	0.13	0.12	0.05 0.05	1	100	85 x 85	192 x 154 x 250
XPR204S	0.1	210	0.2	0.1	0.12	2	240	68 x 68	179.5 x 155 x 237
MCA224S-2S00-U-QP99	0.1	220	0.13	0.12	0.05 0.05	1	100	85 x 85	192 x 154 x 250
XPR304E	0.1	320	0.1		0.04	1.5	82	78 x 73	
XPR404S	0.1	410	0.2	0.1	0.06	2	120	68 x 68	179.5 x 155 x 237
XPR504S	0.1	510	0.3	0.1	0.07	2	140	68 x 68	179.5 x 155 x 237
MCA524S-2S00-U-QP99	0.1	520	0.2	0.2	0.04 0.05	1	82	85 x 85	192 x 154 x 250
XPR504SDR	0.1 1	101 510	0.4	0.8	0.08 0.4	2	160	68 x 68	179.5 x 155 x 237
MCA524P-2S00-A-QP99	0.1 0.2 0.5	120 240 520	0.2	0.2	0.04 0.05	1	82	85 x 85	192 x 154 x 250
Precision Balances									
XPR303SN	1	310	2	1	0.5	1.5		127 x 127	
MCA323S-2S00-R-QP99	1	320	0.6	0.1	0.4 0.5	0.8		140 x 140	
XPR303S	1	310	2	1	0.5	1.5		127 x 127	179.5 x 155 x 236
MCA323S-2S00-E-QP99	1	320	0.6	1	0.4 0.5	0.8		140 x 140	
XPR603SN	1	610	2	1	0.5	1.5		127 x 127	
MCA623S-2S00-R-QP99	1	620	0.6	1	0.4 0.5	0.8		140 x 140	
XPR603S/A	1	610	2	1	0.5	1.5		127 x 127	179.5 x 155 x 236

Model	Readability (mg)	Weighing capacity (g)	Linearity (mg)*	Corner load (mg)*	Typical Repeatability (mg) 5%/max*	Typical Stabilization time (s)*	Pan size (mm)	Chamber size DWH or DH (mm)
Precision Balances (continued)								
XPR603S	1	610	2	1	0.5	1.5	127 x 127	179.5 x 155 x 236
MCA623S-2S00-E-QP99	1	620	0.6	1	0.4 0.5	0.8	140 x 140	
XPR1203S	1	1,210	2	1	0.4	1.5	127 x 127	179.5 x 155 x 236
MCA1203S-2S00-E-QP99	1	1,200	1	1	0.5 0.6	1	140 x 140	
XPR3003S	1	3,100	6	3	0.6	2	127 x 127	179.5 x 155 x 236
MCA3203S-2S00-E-QP99	1	3,200	2	1	0.6 0.6	1	140 x 140	
XPR5003S	1	5,100	5	3	1	2	127 x 127	179.5 x 155 x 236
MCA5203S-2S00-E-QP99	1	5,200	2	1	0.6 0.6	1	140 x 140	
XPR3003SD5	5	3,100	6	5	3	2	205 x 172	
MCA3203S-2S00-R-QP99	1	3,200	2	1	0.6 0.6	1	140 x 140	
XPR6003SD5	5	6,100	7	10	3	2	205 x 172	
XPR603SNDR	1 10	120 610	6	0.8	0.8 4	1.5	127 x 127	
MCA623P-2S00-R-QP99	1 2 5	150 300 620	1.5	3	0.4 0.5	0.8	140 x 140	
XPR603SDR/A	1 10	120 610	6	1.5	0.8 4	1.5	127 x 127	179.5 x 155 x 236
XPR603SDR	1 10	120 610	6	1.5	0.8 4	1.5	127 x 127	179.5 x 155 x 236
MCA623P-2S00-E-QP99	1 2 5	150 300 620	1.5	3	0.4 0.5	0.8	140 x 140	
XPR1202S	10	1,210	6	6	4	1.2	205 x 172	
MCA1202S-2S00-O-QP99	10	1,200	6	10	4 4	0.8	206 x 206	
XPR2002S	10	2,100	6	10	4	1.2	205 x 172	
MCA2202S-2S00-O-QP99	10	2,200	6	10	4 4	0.8	206 x 206	
XPR4002S	10	4,100	6	10	4	1.2	205 x 172	
MCA4202S-2S00-O-QP99	10	4,200	6	10	4 4	1	206 x 206	
XPR6002S	10	6,100	6	10	4	1.2	205 x 172	
MCA6202S-2S00-O-QP99	10	6,200	6	10	4	1	206 x 206	
XPR8002S	10	8,100	6	12	4	1.5	205 x 172	
MCA8202S-2S00-O-QP99	10	8,200	6	10	4 4	1	206 x 206	
XPR10002S	10	10,100	6	12	4	1.5	205 x 172	
MCA10202S-2S00-O-QP99	10	10,200	6	10	5 5	0.8	206 x 206	
XPR15002L	10	15.100	6	25	10	1.5	203 x 170	
MCA14202S-2S00-O-QP99	10	14.200	10	10	5 5	0.8	206 x 206	
XPR6002SDR	10 100	1,200 6,100	20	30	5 0	1.2	205 x 172	
MCA6202P-2S00-O-QP99	10 20 50	1,500 3,000 6,200	20	30	4 15	1	206 x 206	
XPR20002LDR	10 100	4.200 20,100	20	40	18 60	1.5	172 x 205	
MCA20201S-2S00-O-QP99	100	20,200	60	200	20 20	1.5	400 x 300	
XPR4001S	100	4,100	20	60	40	0.8	190 x 223	
XPR6001S	100	6,100	20	60	40	0.8	190 x 223	
MCA5201S-2S00-O-QP99	100	5,200	20	100	20 20	0.8	206 x 206	
XPR8001S	100	8,100	30	60	40	1	190 x 223	
MCA8201S-2S00-O-QP99	100	8,200	30	100	20 20	0.8	206 x 206	
High Capacity Balances								
XPR10001S	100	10,100	30	60	40	1	190 x 223	
XPR10001L	100	10,100	60	100	40	1.5	280 x 360	
MCA11201S-2S00-O-QP99	100	11,200	30	100	20 20	0.8	206 x 206	

Model	Readability (mg)	Weighing capacity (g)	Linearity (mg)*	Corner load (mg)*	Typical Repeatability (mg) 5%/max*	Typical Stabilization time (s)*	Pan size (mm)	Chamber size DWH or DH (mm)
High Capacity Balances (continued)								
XPR16001L	100	16,100	60	100	40	1.5	280 x 360	
MCA20201S-2S00-O-QP99	100	20,200	60	200	20 20	1.5	400 x 300	
XPR32001L	100	32,100	100	100	40	1.5	280 x 360	
MCA36201S-2S00-O-QP99	100	36,200	100	200	20 50	1.5	400 x 300	
XPR64001L	100	64,100	150	150	40	1.8	282 x 362	
MCA70201S-2S00-O-QP99	100	70,200	150	300	40 40	1.5	400 x 300	
XPR32000L	1,000	32,100	200	300	400	1.2	280 x 360	
MCA36200S-2S00-O-QP99	1,000	36,200	200	500	200 200	1	400 x 300	
XPR64000L	1,000	64,100	200	300	400	1.5	282 x 362	
MCA70200S-2S00-O-QP99	1,000	70,200	200	600	200 200	1	400 x 300	

* Typical values

Specifications subject to change without notice.
Copyright Sartorius Lab Instruments GmbH & Co. KG.
Status: 03 | 2024