

Competitor Cross Guide



Quintix® Pro

Quintix® Pro Models with Motorized Level Feet

Sartorius Quintix® Pro	Sartorius Secura®	Mettler Toledo® MX Series	Ohaus Explorer™ EXP	Precisa Series 390 HM
-	-	MX105DU	EXP125D/AD	HM 125SM-FR
SECURA125-1x	SECURA125-1x	MX105	EXP125/AD	HM 125SM
SECURA225D-1x	SECURA225D-1x	MX205DU	EXP225D/AD	HM 225SM-DR
-	-	-	EXP225/AD	-
QTX124IMU-1x	SECURA124-1S	MX104	EXP124/AD	HM 120A
QTX224IMU-1x	SECURA224-1S	MX204	EXP224/AD	HM 220A
QTX324IMU-1x	SECURA324-1S	MX304	EXP324/AD	HM 320A
-	-	-	-	HM 520A
-	SECURA213-1S	-	EXP223/AD	-
QTX323IMU-1x	SECURA313-1S	MX303	-	-
QTX423IMU-1x	SECURA513-1S	-	EXP423/AD	-
QTX623IMU-1x	SECURA613-1S	MX603	EXP623/AD	-
QTX1203IMU-1x	SECURA1103-1S	MX1203	EXP1203/AD	-
QTX1503IMU-1x	-	-	-	-
-	SECURA612-1S	-	-	-
QTX1202IMO-1x	SECURA1102-1S	-	-	-
QTX2202IMO-1x	SECURA2102-1S	MX2002	EXP2202	-
QTX3202IMO-1x	SECURA3102-1S	-	-	-
QTX4202IMO-1x	SECURA5102-1S	MX4002	EXP4202	-
QTX6202IMO-1x	SECURA6102-1S	MX6002	EXP6202	-
-	-	-	EXP8202	-
-	-	MX12002	EXP12202	-
QTX3201IMO-1x	SECURA3101-1S	-	-	-
QTX6201IMO-1x	SECURA6101-1S	MX6001	EXP6201	-
QTX8201IMO-1x	-	MX8001	EXP8201	-
QTX10201IMO-1x	-	-	EXP10201	-
QTX12201IMO-1x	-	MX12001L	-	-

Quintix® Pro Models with LevelControl

Sartorius Quintix® Pro	Sartorius Quintix®	Mettler Toledo® MR Series	Ohaus Explorer™ EXR	A&D BM Series	Precisa Series 360 EP	Radwag® AS PS.5Y Series
QUINTIX125D-1x	QUINTIX125D-1x		EXR125D		-	AS 60/220.5Y
QUINTIX65-1x	QUINTIX65-1x		EXR125		EP 125SM	AS 120.5Y
-	-		EXR225D		EP 225SM-DR	AS 82/220.5Y
-	-		EXR225		-	-
-	QUINTIX64-1S	-	-	-	-	-
QTX124IRU-1x	QUINTIX124-1S	MR104	EXR124	-	EP 120A	AS 160.5Y
QTX224IRU-1x	QUINTIX224-1S	MR204	EXR224	BM-200	EP 220A	AS 220.5Y
QTX324IRU-1x	-	MR304	EXR324	BM-300	EP 320A	AS 310.5Y
-	-	-	-	-	EP 420A	-
-	-	-	-	BM-500	EP 520A	AS 520.5Y
-	QUINTIX213-1S	MR203	EXR223	-	-	PS 210.5Y
QTX323IRU-1x	QUINTIX313-1S	MR303	-	-	EP 320M	PS 360.5Y
QTX423IRU-1x	QUINTIX513-1S	MR503	EXR423	-	-	-
QTX623IRU-1x	QUINTIX613-1S	MR603	EXR623	-	EP 620M	PS 600.5Y
-	-	-	-	-	EP 920M	PS 750.5Y
QTX1203IRU-1x	-	-	EXR1203	-	EP 1220M	PS 1000.5Y
QTX1503IRU-1x	-	-	-	-	EP 2200M	PS 3000.5Y
-	QUINTIX412-1S	-	-	-	-	-
-	QUINTIX612-1S	-	-	-	-	-
QTX1202IRO-1x	QUINTIX1102-1S	MR1002	-	-	EP 1200C	-
QTX2202IRO-1x	QUINTIX2102-1S	MR2002	EXR2202	-	EP 2200C	PS 2100.5Y.M
QTX3202IRO-1x	QUINTIX3102-1S	MR3002	-	-	-	PS 3500.5Y.M
QTX4202IRO-1x	QUINTIX5102-1S	MR4002	EXR4202	-	EP 4200C	PS 4500.5Y.M
QTX6202IRO-1x	QUINTIX6102-1S	MR6002	EXR6202	-	EP 6200C	PS 6100.5Y.M
-	-	-	EXR8202	-	EP 8200C	PS 8100.5Y.M
-	-	-	EXR12202	-	-	PS 10100.5Y.M
QTX3201IRO-1x	QUINTIX2101-1S	-	-	-	-	-
-	QUINTIX5101-1S	-	-	-	-	-
QTX6201IRO-1x	QUINTIX6101-1S	MR6001	EXR6201	-	EP 6200D	-
QTX8201IRO-1x	-	-	EXR8201	-	EP 8200D	-
QTX10201IRO-1x	-	-	EXR10201	-	-	-
QTX12201IRO-1x	-	-	-	-	EP 12200D	-

x = term for country-specific models

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600

 For further information, visit
www.sartorius.com