

Competitor Cross Guide

Quintix[®] Pro

Models with Motorized Level Feet

Sartorius Quintix [®] Pro	Sartorius Secura [®]	Mettler Toledo MX	Precisa HM Series
QTX124IMU-1x	SECURA124-1S	MX104	HM120A
QTX224IMU-1x	SECURA224-1S	MX204	HM220A
QTX324IMU-1x	SECURA324-1S	MX304	HM320A
-	-	-	HM520A
-	SECURA213-1S	-	-
QTX323IMU-1x	SECURA313-1S	MX303	-
QTX423IMU-1x	SECURA513-1S	-	-
QTX623IMU-1x	SECURA613-1S	MX603	-
QTX1203IMU-1x	SECURA1103-1S	MX1203	-
QTX1503IMU-1x	-	-	-
-	SECURA612-1S	-	-
QTX1202IMO-1x	SECURA1102-1S	-	-
QTX2202IMO-1x	SECURA2102-1S	MX2002	-
QTX3202IMO-1x	SECURA3102-1S	-	-
QTX4202IMO-1x	SECURA5102-1S	MX4002	-
QTX6202IMO-1x	SECURA6102-1S	MX6002	-
-	-	MX12002	-
QTX3201IMO-1x	SECURA3101-1S	-	-
QTX6201IMO-1x	SECURA6101-1S	MX6001	-
QTX8201IMO-1x	-	MX8001	-
QTX10201IMO-1x	-	-	-
QTX12201IMO-1x	-	MX12001L	-

Models with LevelControl

Sartorius Quintix® Pro	Sartorius Quintix®	Mettler Toledo MR	Ohaus® Explorer® EX	A&D BM Series	Precisa EP Series	Radwag® AS PS.5Y Series
-	QUINTIX64-1S	-	-	-	-	-
QTX124IRU-1x	QUINTIX124-1S	MR104	EX124	-	EP120A	AS160.5Y
QTX224IRU-1x	QUINTIX224-1S	MR204	EX224	BM-200	EP220A	AS220.5Y
QTX324IRU-1x	-	MR304	EX324	BM-300	EP320A	AS310.5Y
-	-	-	-	-	EP420A	-
-	-	-	-	BM-500	EP520A	AS520.5Y
-	QUINTIX213-1S	MR203	EX223	-	-	PS210.5Y
QTX323IRU-1x	QUINTIX313-1S	MR303	-	-	EP320M	PS360.5Y
QTX423IRU-1x	QUINTIX513-1S	MR503	EX423	-	-	-
QTX623IRU-1x	QUINTIX613-1S	MR603	EX623	-	EP620M	PS600.5Y
-	-	-	-	-	EP920M	PS750.5Y
QTX1203IRU-1x	-	-	EX1103	-	EP1220M	PS1000.5Y
QTX1503IRU-1x	-	-	-	-	EP2200M	PS3000.5Y
-	QUINTIX412-1S	-	-	-	-	-
-	QUINTIX612-1S	-	-	-	-	-
QTX1202IRO-1x	QUINTIX1102-1S	MR1002	-	-	EP1200C	-
QTX2202IRO-1x	QUINTIX2102-1S	MR2002	EX2202	-	EP2200C	PS2100.5Y.M
QTX3202IRO-1x	QUINTIX3102-1S	MR3002	-	-	-	PS3500.5Y.M
QTX4202IRO-1x	QUINTIX5102-1S	MR4002	EX4202	-	EP4200C	PS4500.5Y.M
QTX6202IRO-1x	QUINTIX6102-1S	MR6002	EX6202	-	EP6200C	PS6100.5Y.M
-	-	-	-	-	EP8200C	PS8100.5Y.M
-	-	-	EX10202	-	-	PS10100.5Y.M
QTX3201IRO-1x	QUINTIX2101-1S	-	-	-	-	-
-	QUINTIX5101-1S	-	-	-	-	-
QTX6201IRO-1x	QUINTIX6101-1S	MR6001	EX6201	-	EP6200D	-
QTX8201IRO-1x	-	-	-	-	EP8200D	-
QTX10201IRO-1x	-	-	EX10201	-	-	-
QTX12201IRO-1x	-	-	EX12001	-	EP12200D	-

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
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906



For further information, visit
www.sartorius.com