SARTURIUS

Quintix® Pro



Benefits

- Enhance the user experience with new premium features
- Offering a wide range of use in different environments and applications
- Energy efficiency, reduced carbon footprint, and sustainability, minimizing environmental impact from the manufacturing plant to the bench

Product Information

It's easy to overlook the standard lab balance. To expect the ordinary. Until something new comes along that redefines what a standard balance can be.

Quintix® Pro Laboratory Balances deliver the performance you need, with added premium features that raise the bar in usability, flexibility, and eco-conscience design.

Technical Specifications

General Technical Data

Ambient Conditions		Value
Installation site	For indoor use only, max. height above sea level	3000 m
Temperature	Environment (metrological data)* Environment Storage and transport	+10 - +30° C +5 - +40° C -20 - +60° C
Relative humidity**	At temperatures up to 31° C, non-condensing, then linear decrease from max. 80% at 31° C to max. 50% at 40° C	80%
No heat from heating systems or direct sunlight		
No electromagnetic fields		
Power Supply Device		Value
Input voltage		15 V _{DC} (±10%)
Power consumption, max		4 W
Only by Sartorius power supply unit YEPS01-15V0\	W with interchangeable country-specific plug-in AC adaptors	
Power Supply Unit		Value
Type: Sartorius power supply unit YEPS01-15V0W		
Primary	Voltage Frequency Current consumption, maximum	100 - 240 V _{AC} (±10%) 50 - 60 Hz 0.2 A
Secondary	Voltage Current, maximum	15 V _{DC} (±5%) 0.53 A
Short-circuit protection		Electronic
Protection class according to IEC 60950-1		11
Pollution level according to IEC 61010-1		2
Overvoltage category according to IEC 60664-1		11
Other data: See label on the power supply unit		
Electromagnetic Compatibility		
Interference resistance: Suitable for use in industria	areas	
Transient emissions	Class B Suitable for use in residential areas and areas that are connected to a low voltage network that also supplies residential buildings	
Materials		
Housing	Polybutylene terephthalate (PBT)	
Control module	Glass	
Draft shield	Glass polybutylene terephthalate (PBT)	
Weighing pan	Stainless steel	
Warm-up Time		Value
Device, approx		2 hr

 $^{{}^{\}star} \, \text{For conformity-assessed (verified) balances in accordance with EU requirements, refer to the information on the balance.} \\$

^{**} For conformity-assessed (verified) balances in accordance with EU requirements, the legal regulations apply.

Interfaces			
Specifications of the RS232 Interface			
Type of interface	Serial interface		
Interface operation	Full duplex		
Level	RS232		
Connection	D-sub connector, 9-pin		
Maximum cable length	10 m		
Pin assignment		Die 1. Net endered	Die / Met essiene el
, mossignment	$ \begin{bmatrix} 1 & 0 & 0 & 0 & 0 & 5 \\ 6 & 0 & 0 & 0 & 0 & 9 \end{bmatrix} $	Pin 1: Not assigned Pin 2: Data output (T×D) Pin 3: Data input (R×D) Pin 4: Not assigned Pin 5: Internal ground	Pin 6: Not assigned Pin 7: Clear to Send (CTS) Pin 8: Request to Send (RTS) Pin 9: Not assigned
Specifications for the USB-C Interface (back)			
Communication	USB UTL		
Connectable devices	Sartorius printers, Sartorius secor	nd display, FTDI cable or US	B memory stick
Specifications for the PC-USB Interface (back)			
Communication	USB Device		
Connectable devices	PC		
Specifications for the USB-C Interface (front)			
Communication	USB UTL		
Connectable devices	Sartorius printers, Sartorius secor	nd display, FTDI cable or US	B memory stick
Calibration	·		`
 Internal calibration isoCAL 	3-point calibration		
Internal calibration	External calibration		
Selectable Weight Units*			
Gram, kilogram, carat, pound, ounce, troy ounce, milligram, parts per pound, China tael, mommes, A			
Display			
Intuitive graphic touch technology			
Built-in Applications			
Weighing Dosing	Mixing	 Animal weighing 	3
Mass unit conversion	Components	■ Peakhold	
Counting Descriptions	Density determination	■ Pipette smart ch	
Percentage weighingCalculation	StatisticsCheckweighing	Differential weigUnderfloor weig	
Languages			·····9
Chinese, Czech, Dutch, English, French, German, Russian, Spanish, Turkish	Hungarian, Italian, Japanese, Korean,	Polish, Portuguese,	
Protection			
Chemical resistant housing parts	 Dust cover for balances (also av 		
Glass parts of the draft shield are coated to	■ In-use-cover for housing (also a	-,	
reduce electrostatic influences	In-use-cover for weighing pan ((available as accessory)	
Display foil (available as assessory)			
Anti-theft Lock			
Kensington lock and lockdown capability for cable	e or chain		

 $^{{}^*\}text{The availability of units depends on national legislation and is therefore country-specific.}\\$

Model specific data

For models with internal adjustment feature

Model QTX	Unit	324lxy-1z ¹	224lxy-1z ¹	124lxy-1z ¹	
Readability Scale interval (d)	mg	0.1	0.1	0.1	
Maximum capacity (Max)	g	320	220	120	
Weighing system		EMC	EMC	EMC	
Repeatability					
At 5% load, typical value	± mg	0.08	0.08	0.08	
At approx. maximum load, typical value	± mg	0.1	0.1	0.1	
Linearity deviation					
Limits	± mg	0.2	0.2	0.2	
Typical value	± mg	0.06	0.06	0.06	
Eccentricity deviation (Deviation when load is	off-center, positior	ns according to OIML R76)		
Test weight	g	200	100	50	
Tolerance	± mg	0.4	0.4	0.4	
Typical value	± mg	0.2	0.12	0.1	
Sensitivity drift between +10° C and +30° C	± ppm/K	1	1	1	
Tare maximum capacity (subtractive)		<100%	of maximum capacity		
isoCAL:					
Temperature change	К	1.5	1.5	1.5	
Time interval	h	4	4	4	
For models with approval:					
Accuracy class		I	1	ĺ	
Туре		BC-QA	BC-QB	BC-QB	
Verification scale interval (e)	mg	1	1	1	
Minimum load (Min)	mg	10	10	10	
Minimum initial weighing according to USP (U	nited States Pharm	nacopeia), Chap. 41			
Optimum minimum initial weighing	g	0.082	0.082	0.082	
Typical minimum initial weighing	g	0.16	0.16	0.16	
Typical measurement time	S	≤2.0	≤2.0	≤2.0	
Typical stabilization time	S	≤1.5	≤1.5	≤1.5	
Recommended calibration weight					
External calibrated test weight	g	200	200	100	
Accuracy class in accordance with OIML R1	11-1	E2	E2	E2	
Weighing pan size	mm	Ø 90	Ø 90	Ø 90	
Weighing chamber height*	mm	240	240	240	
Dimension (W x D x H)	mm	220×377×346	220×377×346	220×377×346	
Net weight, approx.	kg	6.70	6.70	6.70	
Gross weight, approx.	kg	8.65	8.65	8.65	

 $^{^{\}star}$ upper edge of the weighing pan to the lower edge of the upper draft shield panel

¹ Levelling, x =

x = R: Real-time level support

x = M: Motorized, automatic level feet

¹ Draft shield, y = y = U : Manual, glass y = O : No draft shield

¹ Country-specific marking in model, z =

z = S: Standard balances without country-specific additions

 $z = \mathsf{SAR}: \mathsf{Standard} \ \mathsf{balances} \ \mathsf{with} \ \mathsf{country-specific} \ \mathsf{additions} \ \mathsf{for} \ \mathsf{Argentina}$

 $z = \mathsf{SJP} : \mathsf{Standard} \ \mathsf{balances} \ \mathsf{with} \ \mathsf{country-specific} \ \mathsf{additions} \ \mathsf{for} \ \mathsf{Japan}$ $z = \mathsf{SKR} : \mathsf{Standard} \ \mathsf{balances} \ \mathsf{with} \ \mathsf{country-specific} \ \mathsf{additions} \ \mathsf{for} \ \mathsf{South} \ \mathsf{Korea}$

z = CEU : Conformity-assessed balances with EU type examination

certificate without country-specific additions
z = CFR: Conformity-assessed balances with EU type examination certificate only for France

z = NUS : Balances with approval for Canada and USA

z = OBR : Balances with approval for Brazil

z = OCN: Balances with approval for China

z = OIN : Balances with approval for India z = OJP : Balances with approval for Japan

z = ORU : Balances with approval for Russia

1503lxy-1z ¹	1203lxy-1z ¹	623lxy-1z ¹	4231xy-1z ¹	3231xy-1z ¹
1	1	1	1	1
1.500	1,200	620	420	320
EMC	EMC	EMC	EMC	EMC
0.5	0.5	0.5	0.5	0.5
1	1	1	1	1
2	2	2	2	2
0.6	0.6	0.6	0.6	0.6
500	500	200	200	200
2	2	2	2	2
1	1	1	1	1
1,5	1,5	2	2	2
		<100% of maximum ca	apacity	
1.5	1.5	2	2	2
4	4	6	6	6
 I	I	II	II	II
BC-QC	BC-QC	BC-QD	BC-QD	BC-QD
10	10	10	10	10
100	100	20	20	20
0.82	0.82	0.82	0.82	0.82
1	1	1	1	1
≤1.5	≤1.5	≤1.5	≤1.5	≤1.5
≤1.0	≤1.0	≤1.0	≤1.0	≤1.0
1,000	1,000	500	200	200
E2	E2	F1	F1	F1
Ø 120	Ø 120	Ø 120	Ø 120	Ø 120
240	240	240	240	240
220×377×346	220×377×346	220×377×346	220×377×346	220×377×346
8.10	8.10	6.70	6.70	6.70
 10.03	10.03	8.65	8.65	8.65

Model QTX	Unit	6202lxy-1z ¹	4202lxy-1z ¹	3202lxy-1z ¹	2202lxy-1z ¹	
Readability Scale interval (d)	mg	10	10	10	10	
Maximum capacity (Max)	g	6,200	4,200	3,200	2,200	
Weighing system		EMC	EMC	EMC	EMC	
Repeatability						
At 5% load, typical value	± mg	5	5	5	5	
At approx. maximum load, typical value	± mg	10	10	10	10	
Linearity deviation						
Limits	± mg	20	20	20	20	
Typical value	± mg	6	6	6	6	
Eccentricity deviation (Deviation when load is	off-center, positior	ns according to OIML R	76)			
Test weight	g	2,000	2,000	2,000	1,000	
Tolerance	± mg	20	20	20	20	
Typical value	± mg	10	10	10	10	
Sensitivity drift between +10° C and +30° C	± ppm/K	2	2	2	2	
Tare maximum capacity (subtractive)		<1	L00% of maximum cap	acity		
isoCAL:						
Temperature change	K	2	2	2	2	
Time interval	h	6	6	6	6	
For models with approval:						
Accuracy class		II	II	II	II	
Туре		BC-QE	BC-QE	BC-QE	BC-QE	
Verification scale interval (e)	mg	100	100	100	100	
Minimum load (Min)	mg	500	500	500	500	
Minimum initial weighing according to USP (U	nited States Pharm	nacopeia), Chap. 41				
Optimum minimum initial weighing	g	8.2	8.2	8.2	8.2	
Typical minimum initial weighing	g	10	10	10	10	
Typical measurement time	S	≤1.0	≤1.0	≤1.0	≤1.0	
Typical stabilization time	S	≤0.9	≤0.9	≤0.9	≤0.9	
Recommended calibration weight						
External calibrated test weight	g	5,000	2,000	2,000	2,000	
Accuracy class in accordance with OIML R1	11-1	F1	F1	F1	F1	
Weighing pan size	mm	182×182	182×182	182×182	182×182	
Weighing chamber height*	mm	-	-	-	-	
Dimension (W x D x H)	mm	215×377×95	215×377×95	215×377×95	215×377×95	
Net weight, approx.	kg	6.20	6.20	6.20	6.20	
Gross weight, approx.	kg	8.40	8.40	8.40	8.40	

 $^{^{\}star}$ upper edge of the weighing pan to the lower edge of the upper draft shield panel

¹ Levelling, x =

x = R : Real-time level support

x = M: Motorized, automatic level feet

¹ Draft shield, y = y = U : Manual, glass y = O : No draft shield

¹ Country-specific marking in model, z =

z = S: Standard balances without country-specific additions

z = SAR : Standard balances with country-specific additions for Argentina

 $z=SJP: Standard\ balances\ with\ country-specific\ additions\ for\ Japan\ z=SKR: Standard\ balances\ with\ country-specific\ additions\ for\ South\ Korea$

z = CEU : Conformity-assessed balances with EU type examination certificate without country-specific additions
z = CFR : Conformity-assessed balances with EU type examination certificate only for France

z = NUS : Balances with approval for Canada and USA

z = OBR : Balances with approval for Brazil z = OCN : Balances with approval for China

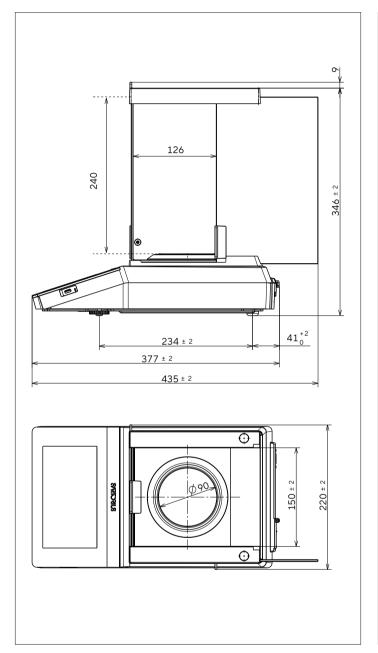
z = OIN : Balances with approval for India z = OJP : Balances with approval for Japan

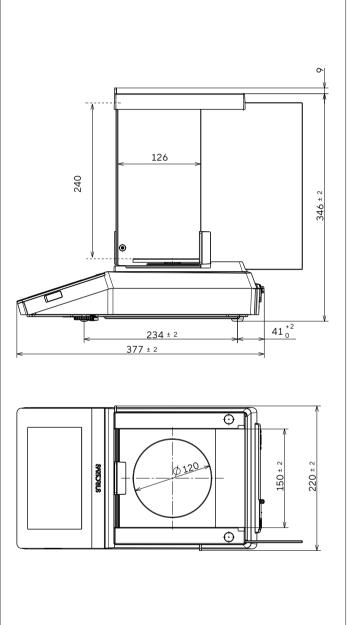
z = ORU: Balances with approval for Russia

1202lxy-1z ¹	12201lxy-1z1	10201lxy-1z ¹	8201lxy-1z1	6201lxy-1z ¹	3201lxy-1z1
10	100	100	100	100	100
1,200	12,200	10,200	8,200	6,200	3,200
EMC	EMC	EMC	EMC	EMC	EMC
5	50	50	50	50	50
10	100	100	100	50	50
20	100	100	100	100	100
6	60	60	60	60	60
500	5,000	5,000	5,000	2,000	2,000
20	200	200	200	200	200
10	100	100	100	100	100
2	4	4	4	2	2
		<100% of	f maximum capacity		
2	2	2	2	2	2
6	6	6	6	6	6
II	II	II	II	II	II
BC-QE	BC-QG	BC-QG	BC-QG	BC-QG	BC-QE
100	1,000	1,000	1,000	100	100
500	5,000	5,000	5,000	5,000	5,000
8.2	82	82	82	82	82
10	100	100	100	100	100
≤1.0	≤1.0	≤1.0	≤1.0	≤1.0	≤1.0
≤0.9	≤0.9	≤0.9	≤0.9	≤0.9	≤0.9
1,000	10,000	10,000	5,000	5,000	2,000
F1	F2	F2	F2	F2	F2
182×182	182×182	182×182	182×182	182×182	182×182
-	-	-	-	-	-
215×377×95	215×377×95	215×377×95	215×377×95	215×377×95	215×377×95
				6.20	6.20
6.20	6.20	6.20	6.20	0.20	0.20

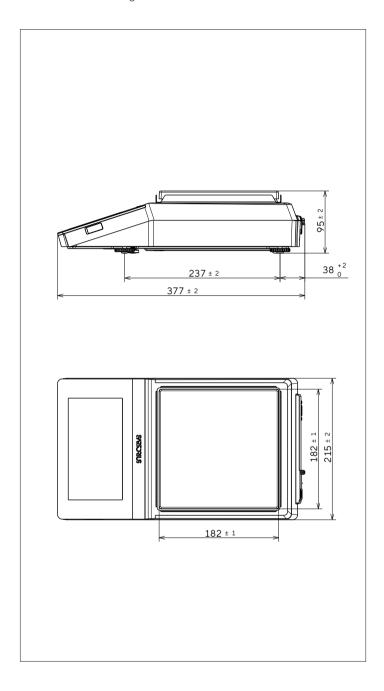
Technical Drawings

Models with a readability of 0.1 mg All dimensions are given in millimeters Models with a readability of 1 mg All dimensions are given in millimeters





Models with a readability of ≥ 10 mg All dimensions are given in millimeters



Accessories

These tables contain an excerpt of the accessories that can be ordered. For information on other products, contact Sartorius.

Balance Accessories

Item	Quantity	Order number
Display protection film (set of 5)	1	YDC40
Dust cover for balances with an analytical draft shield	1	6960SE06
Dust cover for balances with rectangular weighing pan	1	6960SE07
In-use cover for balances with an analytical draft shield (1 piece)	1	6960SE08
In-use cover for balances with an analytical draft shield (set of 5)	1	6960SE08-5
In-use cover for balances with rectangular weighing pan (1 piece)	1	6960SE09
In-use cover for balances with rectangular weighing pan (set of 5)	1	6960SE09-5
Set protection cover for 90 mm weighing pan and draft shield base plate (set of 5)	1	YIC02-A
Set protection cover for 120 mm weighing pan and draft shield base plate (set of 5)	1	YIC02-M
Set protection cover for 182 × 182 mm weighing pan (set of 5)	1	YIC02-P
Round glass draft shield for balances with a readability of 1 mg	1	YDS02QTX
Density determination set		
for solids and liquids for balances with a readability of 0.1 mg \mid 1 mg	1	YDK03
for solids and liquids for balances with a readability of ≥10 mg	1	YDK04QTX
Demo transportation case	1	YDB01QTX
"Kensington Lock" anti-theft device	1	YKL01
Pedal button foot switch	1	YFS04
Second display remote display	1	YSD01
Converter cable 5 VDC > 15 VDC	1	YCC-5V-15V
Weighing table		
Made from wood with natural stone	1	YWT09
Made from natural stone, with vibration dampening	1	YWT03
Wall console made from natural stone	1	YWT04

Printer and Accessories for Data Communication

Quantity	Order number
1	YDP30
1	YDP40
1	YDP50
1	YDP20-0CE
1	YCC-USB-C-B
1	YCC-USB-C-A
1	YCC-D09M-USB-A
1	YCC-D09MM
1	YCC-D09MF
1	YCC-D09M-2D09F
	Quantity 1 1 1 1 1 1 1 1 1 1 1 1 1

 $^{^{\}star} additional\ power\ supply\ e.g.\ YEPS01-PS4\ or\ YEPS01-PS5\ is\ required$

External Calibration and Adjustment Weights

QTX model	Weight	Accuracy class	Order number
324 224	200 g	E2	YCW522-AC-02
124	100 g	E2	YCW512-AC-02
1503 1203	1,000 g	E2	YCW612-AC-02
623	500 g	F1	YCW553-AC-02
423 323	200 g	F1	YCW523-AC-02
6202	5,000 g	F1	YCW653-AC-02
4202 3202 2202	2,000 g	F1	YCW623-AC-02
1202	1,000 g	F1	YCW613-AC-02
12201 10201	10,000 g	F2	YCW714-AC-02
8201 6201	5,000 g	F2	YCW654-AC-02
3201	2,000 g	F2	YCW624-AC-02
2201	2,000 g	F2	YCW624-AC-02

Germany

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



⊕ For further information, visit

www.sartorius.com

USA

Sartorius Corporation 565 Johnson Avenue Bohemia, NY 11716 Phone +1 631 254 4249 Toll-free +1 800 368 7178