

Quintix® and Secura® Standard Laboratory Balances: Reliable Weighing for Faster and Better Results



More Efficient Workflows with Quintix® and Secura®

... Based on Intuitive Operation:

The Start Assistant function and SmartTouch user interface will enable you to operate the balance straight away and feel right at ease using all the functions available. Both the convenient operating design with its easy-to-understand menu structure and the ergonomic balance design have already won several awards. **These features will significantly help accelerate your workflows right from the start for error-free results.**

... Using Integrated Applications:

Built-in standard applications let you instantly extend the options for using the balance, turning special, often complicated, lab tasks into efficient routines. These applications are set up for intuitive use, guiding you in simple steps throughout each of your workflows. **This will save you valuable time and increase the reliability of your results.**

... Made to Deliver the Best Results:

Quickly and easily transfer, evaluate, further process or save and document weighing results in perfect compliance with legal requirements – just by using Plug & Play technology, without needing any extra software or complicated settings. **In this way, you can unlock the full potential of your lab data and reliably save it. As a result, you will have more time for new, intriguing analytical procedures, while being on the safe side.**

“ We use Quintix® non-stop – it gives us the greatest support like no other – even for our most unusual applications. ”



	Quintix®	Secura®
Reliability and Data Security		
isoCAL: internal calibration and adjustment function	+	+
CAL audit trail function	+	+
Menu protection	+	+
Password protection		+
SQmin: monitoring the minimum sample weight for compliance with the USP* operating range		+
Individual security levels		+
Data Transfer and Documentation		
Plug & Play functionality via mini USB port	+	+
PC-Direct feature for data transfer to a computer	+	+
Sample and batch identification		+
Monitored data transfer		+
Leveling		
Leveling guided by the level indicator	+	
Continuous real-time level monitoring		+
Models		
Semi-micro balances	+	+
Micro balance		+
Higher weighing capacity levels**		+

“ Secura® makes it particularly easy for me to ensure I reliably comply with special requirements, such as those in regulated areas. ”



* USP - United States Pharmacopeia

** See Technical Specifications

Monitoring Functions and Security

isoCAL: Reliability Guaranteed

Temperature fluctuations can distort the accuracy of your weighing results. Quintix® and Secura® rule out this risk right from the start with their temperature- and time-controlled calibration and adjustment function, isoCAL. In addition, the internal Cal Audit Trail function documents each calibration and adjustment procedure, generating a 100% traceable audit trail for your quality assurance.

Menu Access Protection

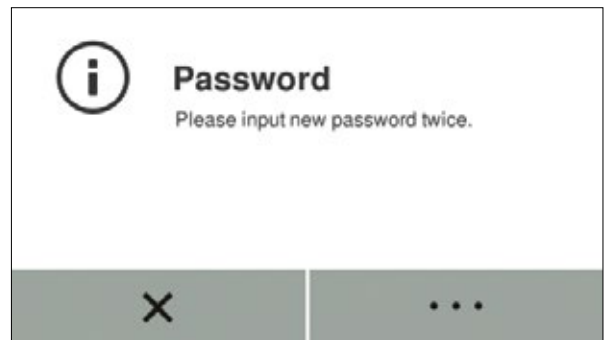
Reliably protect the system settings of your balance from any unintentional changes by activating the tamper-proof supervisor lock on Quintix®.





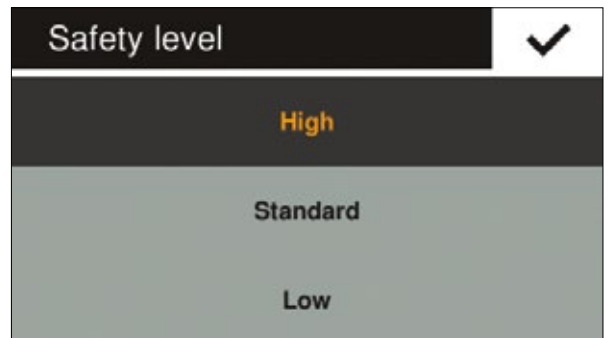
Password-Protected Menu Settings

Secura® offers you customizable security: Once entered, a password prevents any changes to the menu settings that could alter the metrological performance of the balance. In this way, Secura® restricts access to the menu to exclusively authorized persons, ensuring fully controlled protection of your weighing procedures.



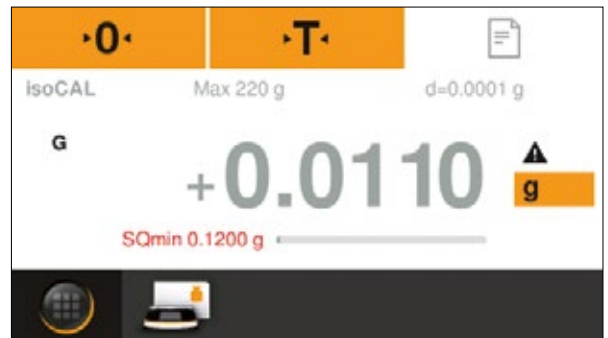
Secure Weighing Data at All Times

Benefit from the different safety level settings that you can select on Secura® for controlled data transfer. You can optimally adapt the automatic monitoring systems of Secura® to your security needs. Whichever level you opt for – high, standard or low – you can be sure that with Secura®, you will be processing only data that is correct.

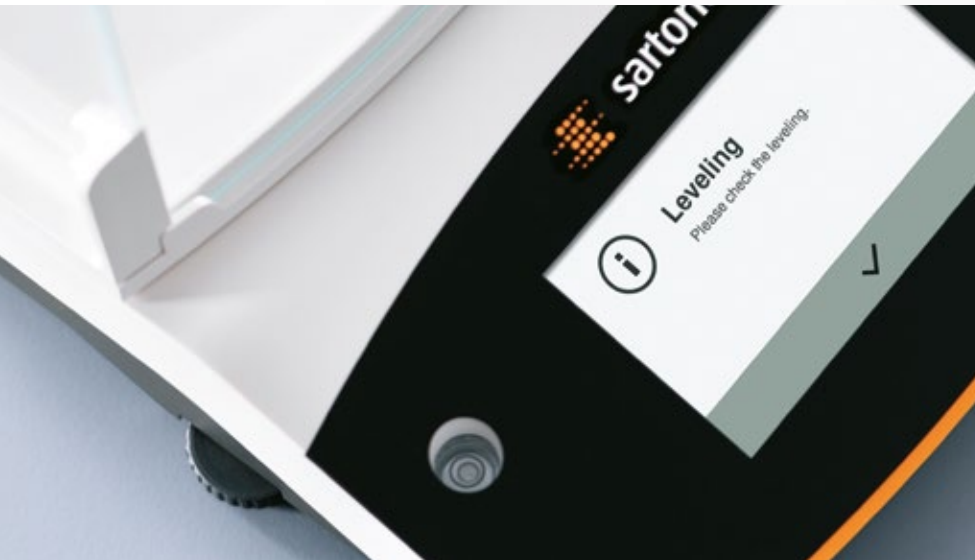


SQmin: USP-Compliant Weighing Results

Compliance with USP guidelines along with absolute process reliability? This is guaranteed by Secura®: the SQmin function checks that your sample quantity is within the operating range of your balance. Secura® will alert you whenever your sample is below the minimum net weight and will temporarily block transfer of data to a printer or a computer. Secura® ensures that only permissible weighing data will be further processed.



Correct Leveling



Level Status at a Glance

Quintix® makes it so simple for you to check on your own whether your balance is correctly level – just one quick glance at the ergonomically positioned level indicator is all it takes. Plus, the two large and easy-to-adjust feet on the front enable you to readily level Quintix®. Your weighing results will always be “rock-steady” with Quintix®.



Continuous Real-Time Level Monitoring

The Secura® LevelControl function ensures that your balance is perfectly level. It continuously monitors the level status in real time, instantly detecting any tilt. In this case, it will respond immediately to ensure data integrity: You will be alerted, and data transfer to a printer or computer will be temporarily blocked. Secura® will then reliably guide you in real time throughout the leveling procedure. With Secura® LevelControl, you will have total peace of mind knowing that your measurements are fully accurate.





Leveling

The balance needs to be leveled

Sartorius

X

START

Documentation and Data Transfer: Easy, Fast and Reliable



Data Transfer: Direct and Correct

Quickly and easily connect your Quintix® or Secura® to your computer or lab printer using the mini USB interface port and Plug & Play technology of the balance. Its PC-Direct feature will transfer your measurement results, without the need for additional software, to the desired Microsoft® Office program – in the format of your choice and ready for calculation in spreadsheets.

Data Documentation: Fast and Easy

Directly connected to a Sartorius laboratory printer, Quintix® and Secura® generate fast and accurate printouts of your weighing results. You can use standard printer paper or practical, continuous-length paper label rolls or even fade-resistant paper for reliable records with unique identification of your samples.



Sample Identification: Even More Reliable and Precise with Secura®

Do you work with samples requiring unique identification in regulated environments? Then you will benefit from the built-in sample identification feature on Secura®: this function enables you to customize the printout information to your specific needs. With an interactive touch-screen display and an alphanumeric input capability, Secura® will make it easy for you to enter complicated sample IDs. Generate sample records reliably and precisely with Secura®.





Quintix® and Secura® Semi-Micro Balances

Unmatched Linearity and Repeatability

Maximize the efficiency of your lab work: the unique temperature management system ensures stable climate conditions within the weighing chamber.

Compensation of Ambient Effects

The user-friendly draft shield featuring One-Touch technology is convenient to use and compensates for even the slightest ambient effects.





Prevention of Cross-Contamination

Quintix® and Secura® semi-micro balances make it easy for you to prevent cross-contamination: For fast, thorough and effortless cleaning, just a few steps are all it takes to dismantle the draft shield and completely remove the rimmed base plate from the weighing chamber. Moreover, this base plate is specially designed to catch all spills. Quintix® and Secura® semi-micro balances enable you to successfully perform even the most sensitive analyses reliably and accurately – to five decimal places.

Applications

Filter Weighing

Quintix® and Secura® semi-micro balances are ideal for high-precision filter weighing. Their ergonomic draft shield and special weighing pan permit easy and reliable operating sequences for weighing filters to an accuracy of five decimal places.



Pipette Calibration

Calibrate your pipettes quickly and accurately at any time with Quintix® and Secura® semi-micro balances. They provide convenient access to the weighing chamber, both from the front and from the side. Used together with a specially matched pipette calibration kit, these balances make calibration of your pipettes as easy as 1-2-3.



Secura® Micro Balance

Do you place extremely high requirements on accurate results for analytical weighing? Obtain the highest quality of results in your analytical weight measurements with perfectly reliable accuracy of six decimal places. Featuring a maximum capacity of 21 g and a resolution of 2 μg over the entire weighing range, the Secura® micro balance determines the exact weight of the smallest quantities – and therefore delivers unsurpassed performance unique throughout the world.



With the Secura® micro balance, you will meet your high standards on accurate results for analytical weighing, even under difficult ambient conditions. The reason: Secura® combines a number of ergonomic features specially designed for user-friendly weighing on high-resolution balances. These include a built-in temperature management system, the proven One-Touch draft shield for convenient, yet well-protected weighing procedures and an easy-to-remove weighing chamber base plate for effortless cleaning. A Secura® micro balance will give you security, speed and six-decimal-place accuracy down to the lowest analytical weighing ranges.





Weighing Pan and Pan Draft Shield Rings Made of High-Grade Titanium

The Secura® micro balance really shines with its unique titanium weighing pan and its two titanium pan draft shield rings supplied as standard equipment. The raised rim of the pan draft shield ring enables you to perform high-precision measurements successfully, even when the balance is exposed to interfering air flow under a laboratory fume hood. Non-magnetic and absolutely inert in contact with aggressive substances, the titanium components of the Secura® micro balance ensure 100% protected weighing.



Integrated Applications

Whether you need to perform common weighing procedures or run dedicated laboratory applications, Quintix® and Secura® will give you the best of both. They provide built-in applications as standard features and offer the widest array of accessories and options to choose from.



Weighing



Perform high-precision weight measurements fast and professionally. Custom-configure the weight unit and number of decimal places. Your Quintix® or Secura® balance will do all of the above in the Weighing app.

Peak Hold



Does an oversized object on the balance hide the display or is a reading displayed only for a split second? The solution: Quintix® and Secura® lock in the reading so you can view it later, taking all the time you need.

Mixing



One formula. Several individual components. Objective: Weigh all into only one container, while keeping an eye on the total weight. Quintix® and Secura® let you weigh in each component sequentially with 100% traceability.

Checkweighing



Are your samples still within the allowable tolerance range? Quintix® and Secura® will give you the answer – on a straightforward color-coded bar graph. You will know at a glance whether your samples get the “green light” so you can proceed.

Components



One formula. Several individual components. Objective: Weigh all components independently of one another. At the same time, check the entire amount weighed in. Quintix® and Secura® enable you to easily keep track of it all for perfect weighing results.

Counting



A bag with an unknown number of identical small parts: Forget counting by hand. Quintix® and Secura® will show you the exact count along with the total weight of all parts – two easy weighing steps are all it takes.

Statistics



Generate statistics on a relatively large sample size at the touch of a button? Yes, you can! Quintix® and Secura® have an internal statistics memory that saves all weighing results so you can retrieve a well-organized list of your key statistics.

Percentage



Need to compare different samples and determine their difference from a reference standard in percent? Quintix® and Secura® along with their Percentage app will take care of all this for you in no time.

Conversion



No time for complicated conversions? Quintix® and Secura® will save you from doing all the number-crunching. Just enter a conversion factor, weigh and instantly read off the weight or value along with the converted result.

Unstable Conditions



Is your work environment extremely hectic, with coworkers running in and out? Or do you have to weigh a mouse that just won't sit still? No problem with Quintix® and Secura®. They will give you absolutely rock-steady results even under unstable ambient conditions.

Density



High-precision density determination of liquids and solids – easy. Just select the Density app and use the accessory density determination kit. This winning combination permits Quintix® and Secura® to accurately determine the quality of precious metals or plastic parts, for example.

A Full Range of Services: Our Mission Is Your Productivity

From installation to a full range of services covering the entire lifecycle of our products – our global Services team focuses on maximizing the reliability and availability of your lab balances, as well as on guaranteeing confidence in your weighing results at all times. Our service specialists regularly complete in-depth training to make sure you obtain the best possible service on site and to meet your specific requirements.

Optimum Equipment Performance

- Our installation service ensures that your lab balances will be up and running perfectly right from the start and deliver full performance - without any compromises.
- Reliably trained staff is essential for accurate and efficient work. That's why our service technicians will see to it that your lab personnel is optimally trained to work with laboratory balances and related equipment.

Quality of Results

- Rely on the accuracy and precision of your weighing results by having regular professional calibration performed on your balances.
- Sartorius offers a number of ISO 17025 accredited USP and manufacturer's certificates that ensure the traceability of weighing results to your national standard and meet the requirements of the widest variety of quality management systems.

Compliance with the Standards of the Regulated Industry

- Sartorius equipment qualification services (IQ/OQ) will provide you with the quality documents you require to operate your equipment in a highly regulated environment.
- Our calibration services include issuing documents that verify the quality and traceability of measurement results in compliance with ISO 17025, GLP | GMP and FDA standards.

Trouble-Free Operation

- Regular maintenance will ensure that your lab balances will continue to operate reliably at peak performance throughout their entire life time, giving you the best cost of ownership.
- Professional repair service, including the use of original spare parts, will enable you to avoid unnecessary downtime.



Installation Service

For maximum equipment performance, rely on our regularly trained service technicians.



Extended Warranty

Reliably protect your investment and obtain full cost transparency with our extended warranty packages.



Equipment Qualification (IQ|OQ)

Our comprehensive and product-specific qualification documents provide proof of the suitability of your equipment, as required by GLP | GMP and FDA standards.



Preventive Maintenance and Service Contracts

Sartorius maintenance service ensures that your equipment will deliver dependable performance you can trust, as well as the reliable quality of your measurement results.



Calibration Service

Our calibration certificates document the quality and traceability of your measurement results in compliance with the requirements of leading quality management systems.



User Training | Academies

Sign up for professional training programs and courses so you and your staff can stay ahead of the curve to successfully meet the fast-paced challenges of evolving technologies and work efficiently.



Accessories

Extend the performance capabilities of your Quintix® and Secura® with accessories from our comprehensive range:



YDP30

Thermal transfer printer* for GxP printouts on standard paper, continuous-length label paper rolls or individual labels; for connection to a USB port or an RS232 port.



YDP40

Direct thermal printer for printouts on standard paper or continuous-length paper label rolls; for connection to a USB port or an RS232 interface port.



YRB11Z

External rechargeable battery pack for standard laboratory balances.



VF4601

Density determination kit for solids and liquids; for balances with a readability of 0.01 mg.



YDK03

Density determination kit for solids and liquids; for balances with a readability of 0.1 mg or 1 mg.



YDK04

Density determination kit for solids and liquids; for balances with a readability of 10 mg (supplied without a balance).



YCP06SQ

Pipette calibration kit (hardware) for balances with a readability of 0.01 mg. The kit consists of a moisture trap and all required adapters.



YDS01SQP

Draft shield for balances with a readability of 10 mg.



YDS02SQP

Round glass draft shield for balances with a readability of 1 mg (supplied without a balance).

* White Paper entitled "GxP-compliant Printing and Archiving of Initial Weights and Related Weighing Data" available free of charge on request

**YFW01SQP**

Special pan with a 130 mm diameter for weighing filters; for balances with a readability of 0.01 mg.

**YSP01SQP**

Weighing pan with an 80 mm diameter, slotted; for increasing the weighing performance of balances with a readability of 0.01 mg.

**YWP01SQP**

Weighing pan with a 90 mm diameter; includes conversion kit; for balances with a readability of 0.01 mg.

**YHK01SQP**

Interior glass draft shield for increasing the weighing performance of balances with a readability of 0.002 mg (supplied without a balance).

**VF4589**

Stainless steel weighing pan with a 50 mm diameter for balances with a readability of 0.002 mg. The set consists of a weighing pan, a low-design and a high-design round pan draft shield (supplied without a balance).

Additional Accessories:

69Y03285	Standard paper and ink ribbon set for YDP30 laboratory printer	6960SE04	Dust cover for balances with a readability of 0.01 mg or 0.002 mg
69Y03286	Self-adhesive paper and ink ribbon set for YDP30 laboratory printer	6960SE03	Dust cover for balances with a readability of 0.1 mg or 1 mg
69Y03287	Standard paper for YDP 30 YDP40 lab printers	YIB01-OUR	Ionizing blower for eliminating electrostatic charges on sample containers and samples
69Y03288	Self-adhesive paper for direct thermal printing for YDP30 YDP40 laboratory printer	YSTP01	Stat-Pen ionizing probe for neutralizing electrostatic charges on samples and filters
YCC04-D09	Data cable, mini USB USB A	YWT09	Wooden balance table with natural stone inset for precise, reliable weighing
YCC03-D09	Data cable, mini USB RS232, 9-pin	YWT03	Balance table made of natural stone, with vibration dampeners
YCC03-D25	Data cable, mini USB RS232, 25-pin	YWT04	Wall console for micro, analytical and precision balances
YDB-SECURA-5	Demo box for 0.01 mg or 0.002 mg balances	641214	Weighing scoop made of chrome nickel steel, 90 mm × 32 mm × 8 mm
YDB-SECURA	Demo box for 0.0001 g or 0.001 g balances	6565-250	Aluminum weighing scoops, 4.5 mg (pkg. of 250) for micro balances
YDB-SECURA-P	Demo box for 0.01 g balances	6566-50	Aluminum weighing scoops, 52 mg (pkg. of 50) for micro balances
6960SE05	In-use covers for balances with a readability of 0.01 mg or 0.002 mg		
6960SE01	In-use covers for balances with a readability of 0.1 mg or 1 mg		
6960SE02	In-use covers for balances with a readability of 0.01 g 0.1 g 1 g		

Quintix® Technical Specifications

Model	Semi-Micro Balances			Analytical Balances	
	125D-1x ¹⁾	65-1x ¹⁾	35-1x ¹⁾	224-1x ¹⁾	
Design	1	1	1	2	
Weighing capacity Capacity levels	g	40 60 120	40 60	30	220
Readability	mg	0.01 0.01 0.1	0.01 0.01	0.01	0.1
Repeatability (standard deviation)	mg	0.03 0.04 0.07	0.03 0.04	0.03	0.1
Repeatability (standard deviation), typical	mg	0.02 0.04 0.07	0.02 0.04	0.03	0.1
Linearity	mg	0.1 0.1 0.2	0.1 0.1	0.1	0.1
Optimal starting point of the operating range*	g	0.0082**	0.0082**	0.0082**	0.082
Sensitivity drift between +10°C and +30°C	± ppm/K	1	1	1	1.5
Typical stabilization time	s	6 6 2	6 6	6	2
isoCAL:					
– Temperature change	K	1.5	1.5	1.5	1.5
– Time interval	h	4	4	4	4
Weighing pan size	mm	Ø 80	Ø 80	Ø 80	Ø 90
Weighing chamber height ²⁾	mm	218	218	218	209
Net weight, approx.	kg	7.8	7.8	7.8	4.9
Dimensions, D x W x H	mm	376 x 214 x 316			359 x 214 x 319
	inches	14.8 x 8.4 x 12.4			14.1 x 8.4 x 12.6
Version verified for legal metrology available		-	-	-	Yes
Interface port		Mini USB			
		Precision Balances			
Model		5102-1x ¹⁾	3102-1x ¹⁾	2102-1x ¹⁾	1102-1x ¹⁾
Design		4	4	4	4
Weighing capacity	g	5,100	3,100	2,100	1,100
Readability	mg	10	10	10	10
Repeatability (standard deviation)	mg	10	10	10	10
Repeatability (standard deviation), typical	mg	10	10	10	10
Linearity	mg	20	20	30	30
Optimal starting point of the operating range*	g	8.2	8.2	8.2	8.2
Sensitivity drift between +10°C and +30°C	± ppm/K	3	3	4	4
Typical stabilization time	s	1	1	1.5	1.5
isoCAL:					
– Temperature change	K	2	2	2	2
– Time interval	h	6	6	6	6
Weighing pan size	mm	Ø 180	Ø 180	Ø 180	Ø 180
Weighing chamber height ²⁾	mm	-	-	-	-
Net weight, approx.	kg	5.2	5.2	4.7	4.7
Dimensions, D x W x H	mm	359 x 214 x 94			359 x 214 x 88
	inches	14.1 x 8.4 x 3.7			14.1 x 8.4 x 3.5
Version verified for legal metrology available		Yes	Yes	Yes	Yes
Interface port		Mini USB			

* The USP (United States Pharmacopeia) Chapter 41 defines the optimal operating range as being from 820 d to the maximum capacity; This value can typically be greater, depending on the place where the balance is set up and on the ambient conditions

** Used together with the slotted 80 mm weighing pan, YSP01SQP



Milligram Balances						Precision Balances
124-1x ¹⁾	64-1x ¹⁾	613-1x ¹⁾	513-1x ¹⁾	313-1x ¹⁾	213-1x ¹⁾	6102-1x ¹⁾
2	2	3	3	3	3	4
120	60	610	510	310	210	6,100
0.1	0.1	1	1	1	1	10
0.1	0.1	1	1	1	1	10
0.1	0.1	1	1	1	1	10
0.1	0.1	2	2	2	2	20
0.082	0.082	0.82	0.82	0.82	0.82	8.2
1.5	1.5	3	3	3	3	3
2	2	1	1	1	1	1
1.5	1.5	2	2	2	2	2
4	4	6	6	6	6	6
Ø 90	Ø 90	Ø 120	Ø 120	Ø 120	Ø 120	Ø 180
209	209	209	209	209	209	-
4.9	4.9	4.9	4.9	4.9	4.9	5.2
						359 x 214 x 94 14.1 x 8.4 x 3.7
Yes	Yes	Yes	Yes	Yes	Yes	Yes
612-1x ¹⁾	412-1x ¹⁾	6101-1x ¹⁾	5101-1x ¹⁾	2101-1x ¹⁾	6100-1x ¹⁾	5100-1x ¹⁾
4	4	4	4	4	4	4
610	410	6,100	5,100	2,100	6,100	5,100
10	10	100	100	100	1,000	1,000
10	10	100	100	100	500	500
10	10	100	100	100	500	500
30	30	300	300	300	1,000	1,000
8.2	8.2	82	82	82	820	820
4	4	8	8	8	8	8
1.5	1.5	1.5	1.5	1.5	1	1
2	2	2	2	2	2	2
6	6	6	6	6	6	6
Ø 180	Ø 180	Ø 180	Ø 180	Ø 180	Ø 180	Ø 180
-	-	-	-	-	-	-
4.7	4.7	4.7	4.7	4.7	4.7	4.7
Yes	-	Yes	Yes	-	Yes	Yes

¹⁾ Placeholder for possible names of country-specific model versions

²⁾ Top edge of the weighing pan to the bottom edge of the uppermost draft shield plate

Secura® Technical Specifications

Model		Micro Balance	Semi-Micro Balances	
		26-1x ¹⁾	225D-1x ¹⁾	125-1x ¹⁾
Design		1	2	2
Weighing capacity Capacity levels	g	21	60 120 220	60 120
Readability	mg	0.002	0.01 0.01 0.1	0.01 0.01
Repeatability (standard deviation)	mg	0.004	0.03 0.04 0.07	0.03 0.07
Repeatability (standard deviation), typical	mg	0.003	0.02 0.04 0.07	0.02 0.07
Linearity	mg	0.01	0.1 0.1 0.2	0.1 0.1
Optimal starting point of the operating range*	g	0.00164**	0.0082***	0.0082***
Sensitivity drift between +10°C and +30°C	± ppm/K	1	1	1
Typical stabilization time	s	8	6 6 2	6 6
isoCAL:				
– Temperature change	K	1.5	1.5	1.5
– Time interval	h	4	4	4
Weighing pan size	mm	Ø 50	Ø 80	Ø 80
Weighing chamber height ²⁾	mm	218	218	218
Net weight, approx.	kg	8.0	7.8	7.8
Dimensions, D x W x H	mm inches	376 x 214 x 316 14.8 x 8.4 x 12.4		
Version verified for legal metrology available		Yes	Yes	Yes
Interface port		Mini USB		
Model		Milligram Balances	Precision Balances	
		213-1x ¹⁾	6102-1x ¹⁾	5102-1x ¹⁾
Design		4	5	5
Weighing capacity	g	210	6,100	5,100
Readability	mg	1	10	10
Repeatability (standard deviation)	mg	1	10	10
Repeatability (standard deviation), typical	mg	1	10	10
Linearity	mg	2	20	20
Optimal starting point of the operating range*	g	0.82	8.2	8.2
Sensitivity drift between +10°C and +30°C	± ppm/K	2	2	2
Typical stabilization time	s	1	1	1
isoCAL:				
– Temperature change	K	2	2	2
– Time interval	h	6	6	6
Weighing pan size	mm	Ø 120	Ø 180	Ø 180
Weighing chamber height ²⁾	mm	209	-	-
Net weight, approx.	kg	5.1	5.2	5.2
Dimensions, D x W x H	mm inches	359 x 214 x 319 14.1 x 8.4 x 12.6	359 x 214 x 94 14.1 x 8.4 x 3.7	
Version verified for legal metrology available		Yes	Yes	Yes
Interface port		Mini USB		

* The USP (United States Pharmacopeia) Chapter 41 defines the optimal operating range as being from 820 d to the maximum capacity; This value can typically be greater, depending on the place where the balance is set up and the ambient conditions

** Used together with the interior glass draft shield, YHK01QP

*** Used together with the slotted 80 mm weighing pan, YSP01SQP



Analytical Balances

Milligram Balances

324-1x¹⁾	224-1x¹⁾	124-1x¹⁾	1103-1x¹⁾	613-1x¹⁾	513-1x¹⁾	313-1x¹⁾
2	3	3	4	4	4	4
320	220	120	1,100	610	510	310
0.1	0.1	0.1	1	1	1	1
0.1	0.1	0.1	1	1	1	1
0.1	0.1	0.1	1	1	1	1
0.3	0.2	0.2	2	2	2	2
0.082	0.082	0.082	0.82	0.82	0.82	0.82
1	1.5	1.5	1,5	2	2	2
2	2	2	1,5	1	1	1
1.5	1.5	1.5	1.5	2	2	2
4	4	4	4	6	6	6
Ø 90	Ø 90	Ø 90	Ø 120	Ø 120	Ø 120	Ø 120
218	209	209	209	209	209	209
7.9	5.1	5.1	5.9	5.1	5.1	5.1
	359 x 214 x 319 14.1 x 8.4 x 12.6					
Yes	Yes	Yes	Yes	Yes	Yes	Yes

3102-1x¹⁾	2102-1x¹⁾	1102-1x¹⁾	612-1x¹⁾	6101-1x¹⁾	3101-1x¹⁾
5	5	5	5	5	5
3,100	2,100	1,100	610	6,100	3,100
10	10	10	10	100	100
10	10	10	10	50	50
10	10	10	10	50	50
20	20	20	20	100	100
8.2	8.2	8.2	8.2	82	82
2	2	2	2	2	2
1	1	1	1	1	1
2	2	2	2	2	2
6	6	6	6	6	6
Ø 180	Ø 180	Ø 180	Ø 180	Ø 180	Ø 180
-	-	-	-	-	-
5.2	5.2	5.2	4.7	5.2	5.2
Yes	Yes	Yes	Yes	Yes	Yes

¹⁾ Placeholder for possible names of country-specific model versions

²⁾ Top edge of the weighing pan to the bottom edge of the uppermost draft shield plate

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

France & Suisse Romande

Sartorius France
2, rue Antoine Laurent de Lavoisier
ZA de la Gaudrée
91410 Dourdan

Phone +33.1.70.62.50.00
Fax +33.1.64.59.76.39

Austria

Sartorius Austria GmbH
Modecenterstrasse 22
1030 Vienna

Phone +43.1.7965760.0
Fax +43.1.7965760.24

Belgium

Sartorius Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles

Phone +32.2.756.06.90
Fax +32.2.481.84.11

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki

Phone +358.9.755.951
Fax +358.9.755.95.200

Hungary

Sartorius Hungária Kft.
Kagyló u. 5.
2092 Budakeszi

Phone +3623.457.227
Fax +3623.457.147

Ireland

Sartorius Ireland Ltd.
Unit 41, The Business Centre
Stadium Business Park
Ballycoolin Road
Dublin 11

Phone +353.1.8089050
Fax +353.1.8089388

Italy

Sartorius Italy S.r.l.
Viale A. Casati, 4
20835 Muggiò (MB)

Phone +39.039.4659.1
Fax +39.039.4659.88

Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001
Fax +31.30.60.52.917

info.netherlands@sartorius.com

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn

Phone +48.61.6473830
Fax +48.61.6473839

Russian Federation

LLC "Sartorius RUS"
Uralskaya str. 4, Lit. B
199155 St. Petersburg

Phone +7.812.327.53.27
Fax +7.812.327.53.23

Spain & Portugal

Sartorius Spain, S.A.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)

Phone Spain +34.913.586.095
Phone Portugal +351.800.855.800
Fax Spain +34.913.589.623
Fax Portugal +351.800.855.799

Switzerland

Sartorius Mechatronics Switzerland AG
Ringstrasse 24a
8317 Tagelswangen (ZH)

Phone +41.44.746.50.00
Fax +41.44.746.50.50

U.K.

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9QQ

Phone +44.1372.737159
Fax +44.1372.726171

Ukraine

LLS "Sartorius RUS"
Post Box 440 "B"
01001 Kiev, Ukraine

Phone +380.44.411.4918
Fax +380.50.623.3162

Americas

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716

Phone +1.631.254.4249
Toll-free +1.800.635.2906
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900
Fax +55.11.4362.8901

Canada

Sartorius Canada Inc.
2179 Dunwin Drive #4
Mississauga, ON L5L 1X2

Phone +1.905.569.7977
Toll-Free +1.800.668.4234
Fax +1.905.569.7021

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepetzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepetzotlan,
Estado de México,
C.P. 54605

Phone +52.55.5562.1102
Fax +52.55.5562.2942

leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C.
Av. Emilio Cavenecia 264 San Isidro
15073 Lima, Perú

Phone +51.1.441 0158
Fax +51.1.422 6100

Asia | Pacific

Australia

Sartorius Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius (Shanghai) Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China

Phone +86.21.6878.2300
Fax +86.21.6878.2882

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwun Tong
Kowloon, Hong Kong

Phone +852.2774.2678
Fax +852.2766.3526

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
562 123 Bangalore, India

Phone +91.80.4350.5250
Fax +91.80.4350.5253

Japan

Sartorius Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.3740.5408
Fax +81.3.3740.5406

Malaysia

Sartorius Malaysia Sdn. Bhd
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil

57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Singapore Pte. Ltd
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park II
Singapore 117528

Phone +65.6872.3966
Fax +65.6778.2494

South Korea

Sartorius Korea Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, Bundang-Gu
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700
Fax +82.31.622.5799

Thailand

Sartorius (Thailand) Co. Ltd.
129 Rama 9 Road,
Huaykwang
Bangkok 10310

Phone +66.2643.8361-6
Fax +66.2643.8367



◀ www.sartorius.com