



The New Gold and Carat Balance Series II

The Highest Precision
"Made in Germany"

Simplifying Progress

SARTORIUS

Secure operating procedures and precise weighing results are indispensable for determining the weight of precious metals and gemstones. Fast work processes and harsh environmental conditions require robust balances that provide consistently reliable and accurate results.

Sartorius offers you a wide array of gold and carat balances. This selection of analytical and precision weighing instruments includes verified models that meet the metrological requirements for legal-for-trade weight measurements.



Sartorius GCL Carat Balances

- Draft shield with a weighing chamber height of 160 mm
- Automatic, built-in calibration and adjustment function
- Monolithic weigh cell
- Hybrid display (LED and touch technology)

Model	Weighing Capacity	Readability	Weighing Pan Size
1103i*	1,100 ct	0.001 ct	Ø 90 mm
603i*	600 ct	0.001 ct	Ø 90 mm



Sartorius GL Gold Balances

- Milligram balances feature a draft shield with a weighing chamber height of 160 mm
- Automatic, built-in calibration and adjustment function available as an option
- Hybrid display (LED and touch technology)

Model	Weighing Capacity	Readability	Weighing Pan Size
523*	520 g	0.001 g	Ø 120 mm
323*	320 g	0.001 g	Ø 120 mm
5202*	5,200 g	0.01 g	Ø 180 mm
4202*	4,200 g	0.01 g	Ø 180 mm
2202*	2,200 g	0.01 g	Ø 180 mm
822*	820 g	0.01 g	Ø 180 mm
8201*	8,200 g	0.1 g	Ø 180 mm



* Model optionally available as verified for legal for trade

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	YDC10
Dust cover for balances with an analytical draft shield	1	6960BC02
Density determination set for solids and liquids for balances with a readability of 0,001 ct 1 mg	1	YDK06
"Kensington Lock" anti-theft device	1	YKL01
Pedal button foot switch	1	YFS03
Second display remote display	1	YSD01
Below-balance weighing (not for models in legal metrology)		
Hook M5	1	69EA0039
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

Item	Quantity	Order number
Thermal printer (USB-B)	1	YDP40
Thermal transfer printer (USB-B, RS232)	1	YDP30
Dot matrix printer (RS232)**	1	YDP20-OCE
Data cable USB-C > USB-B (>YDP30, YDP40)	1.5 m	YCC-USB-C-B
Data cable USB-C > USB-A (> PC)	1.5 m	YCC-USB-C-A
Data cable RS232 (9-pin) > USB-A (> PC)	1.5 m	YCC-D09M-USB-A
Data cable RS232 (9-pin) male > RS232 (9-pin) male (>YDP30)	1.5 m	YCC-D09MM
Data cable RS232 (9-pin) male > RS232 (9-pin) female (>YDP20-OCE, YSD01)	1.5 m	YCC-D09MF
Y-adaptor RS232 (9-pin) male > 2x RS232 (9-pin) female	0.5 m	YCC-D09M-2D09F

** additional power supply e.g. YEPS01-PS4 or YEPS01-PS5 is required

Weighing bowls for gold balances

1000 ml, wt. ~240 g, Ø 186 mm, ht. 77 mm, stainless steel	641211
500 ml, wt. ~113 g, Ø 151 mm, ht. 60 mm, stainless steel	641212
350 ml, wt. ~75 g, Ø 180 mm, ht. 22 mm, stainless steel	YWPO4G

Weighing bowls for carat balances

85 ml, wt. ~11 g, Ø 83 mm, ht. = 23 mm, aluminum	YWPO6G
--	--------


For more accessories and options, please see the Sartorius catalog and price list or visit our website.

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