

Secura[®] and Quintix[®]
First-Class Performance in the 5-Digit Range



0.000001g

The New Semi-Micro Balances: Best in Class

The new Sartorius semi-micro balances feature unsurpassed weighing technology available within their 5-digit class in the standard market segment today. We offer a unique selection, giving you five different models to choose from. These semi-micro balances deliver outstanding performance at an affordable price. Plus, they are backed by the excellent quality standards of products "Made in Germany".

■ Top Performance

The latest technologies for even more performance:
Benefit from unbeatable repeatability and linearity
in the semi-micro range.

■ Easy-to-Clean Design

Removable draft shield and weighing chamber base:
All parts of the draft shield and weighing chamber can
be quickly and easily disassembled – which helps ensure
the best possible cleaning results. And if you happen to
spill any liquid or granular samples while weighing, no
problem for these five new balances. Their weighing
chamber base, featuring a revolutionary new rimmed
design, can be completely taken out of the chamber and
readily wiped cleaned.



■ Intuitive Operation

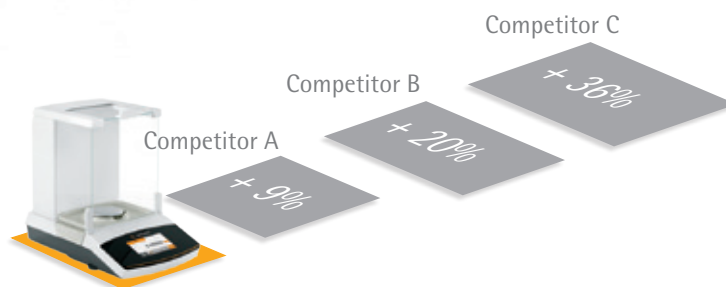
Start Assistant function and SmartTouch user interface: These easy-to-understand features make it so simple to set up and work with the new semi-micro balances. Forget having to read an instruction manual. The display prompts show you all you need to know – helping you increase weighing efficiency in your lab.

■ Ergonomic Draft Shield

One-Touch technology: The side doors of the draft shield open at a simple touch, enabling you to load a sample conveniently – without any mishaps. The draft shield is designed to compensate for ambient conditions, reducing their impact on your weighing results to a bare minimum.

■ Space-saving Footprint

Compact footprint and integrated electronics: Save up to 36% space compared with other competitor models thanks to the compact design and integrated electronics of these semi-micro balances.



Accessories

Weighing Pans

Order No.	Description
YSP01SQP	Weighing pan, 80 mm, slotted
YWP01SQP	Weighing pan, 90 mm; includes conversion kit
YFW01SQP	Filter weighing pan, 130 mm

Rechargeable Battery Pack

Order No.	Description
YRB11Z	For standard lab balances

Printer

Order No.	Description
YDP30	Premium GLP lab printer
69Y03285	Set of standard paper and ink cartridge
69Y03286	Set of self-adhesive labels and ink cartridge
69Y03287	Standard thermosensitive paper
69Y03288	Self-adhesive thermo-sensitive paper

Data Communication Cables

Order No.	Description
YCC04-D09	Data communication cable, USB USB A
YCC03-D09	Data communication cable, mini USB RS232, 9-pin
YCC03-D25	Data communication cable, mini USB RS232, 25-pin

Density Kit

Order No.	Description
YDK05	For semi-micro balances

Protective Covers

Order No.	Description
6960SE04	Dust cover
6960MA03	In-use dust cover



Weighing pan, 80 mm, for increasing the weighing performance



Weighing pan, 90 mm; includes conversion kit



Filter weighing pan, 130 mm

Unique Features for Higher Performance and Flexibility

■ All five models have the familiar Secura® and Quintix® features that will enhance the efficiency of your processes:

□ Fast Data Transfer

Just connect the balance to your computer and start weighing right away – without the need for additional software. Thanks to the balance's Plug & Play technology, your weighing results will be transferred instantly to the program you select. As a result, this makes "number-crunching" fast and easy using spreadsheet calculation programs.

□ Automatic Adjustment

The built-in adjustment function isoCAL will alert you whenever a defined interval has elapsed or a temperature limit has been exceeded. The balance will then perform fully automatic span adjustment using its internal weight. If the balance prompts you that it is ready to perform adjustment while you are weighing, just touch the button to have adjustment carried out later. Each adjustment procedure is documented by the Cal Audit Trail function and is 100% traceable for your quality assurance.

□ GLP | GMP Documentation

Print your weighing results within seconds in compliance with GLP | GMP. In addition, the date and time stamp enables you to generate traceable records. On Secura® balances, you can even customize the printout according to your specific requirements.



- [illegible]

The integrated LevelControl function checks whether your Secura® is level. It detects any tilt immediately, showing you on the display the steps you need to take to level the balance correctly.

Use the built-in identification function to assign unique alpha-numeric IDs to your samples and batches. With the YDP30 laboratory printer connected to the semi-microbalance, you can print your data easily and individually on standard paper or continuous labels.

Activate the password function to protect all your settings from any unintentional changes that could alter the metrological performance of your balance.

In this way, you can be sure that your Secura® will deliver secure weighing results.



Model		225D	125
Weighing capacity ranges	g	60 120 220	60 120
Readability ranges	mg	0.01 0.01 0.1	0.01 0.01
Repeatability (typical standard deviation)	mg	0.02 0.04 0.07	0.02 0.04
Linearity	mg	0.1 0.1 0.2	0.1 0.1
Starting point of the operating range according to USP (typical)*	mg	25	25
Sensitivity drift between +10°C...+30°C (50°F...86°C) ± ppm/K		1	1
Typical stabilization time	s	6 6 2	6 6
isoCAL:			
– Temperature change	K	1.5	1.5
– Time interval	h	4	4
Display update rate (depends on filter level selected)	mm	0.2 0.4	0.2 0.4
Weighing pan size		Ø 80 (Ø 90 optional)	Ø 80 (Ø 90 optional)
Weighing chamber height	mm	218	218
Net weight, approx.	kg	7.8	7.8
Dimensions, D × W × H	mm inches	376 × 214 × 316 14.8 × 8.4 × 12.4	

Three identical analytical balances are shown side-by-side. Each balance has a white body with a transparent draft shield and a black base. The digital display on the front of each balance shows the number '0.0000'.

Model		125D	65	35
Weighing capacity ranges	g	40 60 120	40 60	30
Readability ranges	mg	0.01 0.01 0.1	0.01 0.01	0.01
Repeatability (typical standard deviation)	mg	0.02 0.04 0.07	0.02 0.04	0.03
Linearity	mg	0.1 0.1 0.2	0.1 0.1	0.1
Sensitivity drift between +10°C...+30°C (50°F...86°F)	± ppm/K	1	1	1
Typical stabilization time	s	6 6 2	6 6	6
isoCAL:				
– Temperature change	K	1.5	1.5	1.5
– Time interval	h	4	4	4
Display update rate (depends on filter level selected)	s	0.2 0.4	0.2 0.4	0.2 0.4
Weighing pan size	mm	Ø 80 (Ø 90 optional)	Ø 80 (Ø 90 optional)	Ø 80 (Ø 90 optional)
Weighing chamber height	mm	218	218	218
Net weight, approx.	kg	7.8	7.8	7.8
Dimensions, D × W × H	mm inches	376 × 214 × 316 14.8 × 8.4 × 12.4		

* in combination with weighing pan, 80 mm, slotted YSP01SQP

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Weender Landstrasse 94–108
37075 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
Modectcenterstrasse 22
1030 Vienna

Phone +43.1.7965760.0
Fax +43.1.7965760.24

Belgium

Sartorius Belgium N.V.
Leuvensesteenweg, 248/B
1800 Vilvoorde

Phone +32.2.756.06.71
Fax +32.2.253.45.95

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 ICR" and LLC "Biohit"
Uralskaya str. 4, Lit. B
199155, Saint-Petersburg

Phone +7.812.327.5.327
Fax +7.812.327.5.323

Spain & Portugal

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

Phone Spain +34.902.123.367
Phone Portugal +351.800.855.800
Fax Spain +34.91.358.96.23
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

LLC "Biohit"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918
Fax +380.50.623.3162

America

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.
Circuito Circunvalación Poniente
No. 149
Ciudad Satélite
53100, Estado de México
México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

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 Scientific
Instruments (Beijing) Co., Ltd.
33 Yu An Road, Airport Industrial Park
Zone B,
Shunyi District, Beijing 101300,
P.R.China

Phone +86.10.8042.6300
Fax +86.10.8042.6486

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