



Cubis[®] II Ultra-High Resolution Balances

The New Benchmark in
Performance

Simplifying Progress

SARTORIUS

Introducing the new Semi-Micro and High-Capacity Micro Balances

Uncompromized Accuracy and High Speed Under Real Laboratory Conditions

A high-resolution laboratory balance is used to weigh very small sample quantities and is broadly used across many industries, such as pharmaceutical, biopharmaceutical, or chemical industry in analytical, research, QC labs or at production sites. Most of these industries are subject to strict legal regulations or their SOPs and are looking for a laboratory balance which meets their requirements for very high accuracy and repeatability over the entire weighing range.

The new High-Capacity Microbalance and the new Semi-Micro balance are the latest addition to our Cubis® II portfolio. These instruments were designed to deliver fast results and lowest minimum sample weight at the highest level of accuracy and precision. Thanks to the novel engineering solutions, they are less prone to environmental effects, resulting in faster weighing workflows and improved performance in a daily laboratory environment. The adaptable design allows customization options for both hardware and software, also after post purchase. This level of upgradeability is unparalleled in the market and sets a new standard for investment security in premium laboratory balances.

Your Benefits

- Minimize your sample costs by lowest sample weight through the entire weighing range
- Speed up your weighing process with industry leading measurement time
- 21CFR part 11 and EU Annex 11 compliance, all controls available directly on your balance
- 100% discharged samples due to the novel ionizing technology guarantees fast stabilization
- Thanks to the intuitive disassembly and our new Cleaning App, we made cleaning easier, safer, and traceable
- Save sample costs with min weight below 10 mg and max capacity up to 220 g or below 5 mg and max capacity up to 111 g
- Choose from the following semi-micro balances:
 - 10 microgram readability up to 220g capacity
 - 5 microgram readability up to 220g capacity
- Choose from the following high-capacity microbalance:
 - 1 microgram readability up to 61g capacity
 - 2 microgram readability up to 111 g capacity

Lab-Verified Performance



Peak Performance Under Changing Laboratory Conditions

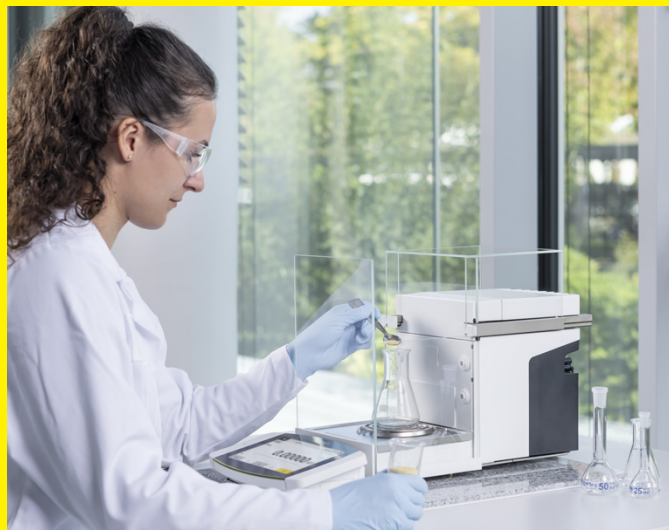
High resolutions instruments are prone to environmental influences. In a lab, changes in temperature such as air-pressure, humidity, or in air-draft can lead to unstable behavior and long measurement time. Static charges occur frequently, when working with small powder compounds. This results in drifting values. The environmental adaptability of these devices help with compensating or mitigating these effects. Built-in intelligent compensating systems manage temperature-, humidity- and air pressure changes. The powerful integrated ionizer guarantees complete electrostatic elimination. Hence, high degree of repeatability and fast measurement time is to be expected when working with these devices.

Find out more about lab-verified performance results in our App Note: Cubis® II Ultra-High Resolution Balances: From Data Sheet to Reality.

Minimize your Sample Costs with Low Minimum Sample Weight

The Cubis® II High-Capacity Microbalance and Semi-Micro Balance ensures stable and reliable performance over the entire weighing range resulting in low minimum sample weight. Minimum weight under 10 mg can be easily achieved with our 220 g capacity model 226S. For minimum weight under 5 mg we recommend our models up to 110 g capacity.

Learn More:
Real Lab Tests App Note



Ergonomic Sample Weighing Without Sample Loss

Increased weighing pan size and spacious inner weighing area supports easy handling of samples and flasks. Additionally, a great variety of sample holders are available which enable weighing in workflows where using a standard pan is difficult or even impossible. All weighing pans and sample holders are made of high-quality titanium. This ensures uncompromised accuracy due to its non-magnetic properties.



**YSH02-3 Adjustable
Flask Holder**



**YCP07MC Pipette
Calibration Kit**



YSH22-3 Vial Holder



**YSH30-3 Filter Holder
(150 mm)**

Compliant Cleanability

Cleaning of the analytical balances is often neglected, despite of the chance of cross-contamination when uncleaned regularly and properly. The higher the resolution, the greater the risk, because even a small particle can negatively impact the weighing accuracy.

Tool-Free Assembly and High Chemical Resistance

All draft-shields, including the front-shield, as well as the weighing pans and the base plate, can be easily removed allowing a tool-free, intuitive assembly. Both the weighing floor and the weighing pan show high chemical resistance against the most frequently used cleaning agents in the laboratory. The inner weighing system is protected thanks to the design of the weighing floor.

Software Guidance and Process Documentation

Follow the steps in the new cleaning QApp for error free cleaning execution. It supports your daily and advanced cleaning routines, shows the cleaning status and allows tracing the process in the audit trail. The Cleaning QApp is configurable with user management allowing the documentation, including electronic signature. The cleaning QApp comes with every MCA Ultra-High Resolution Balance. Find out more cleaning in Best Cleaning Practices.

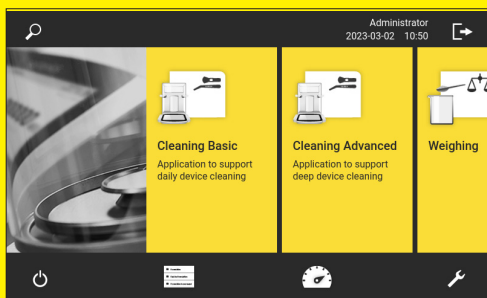


**Learn More:
Cleaning Practices for Cubis® II
Ultra-High Resolution Balances**

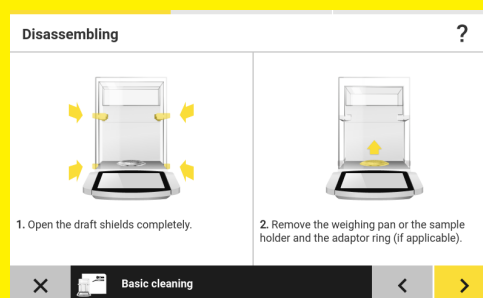




Integrated Cleaning QApp Makes Cleaning Easy and Traceable



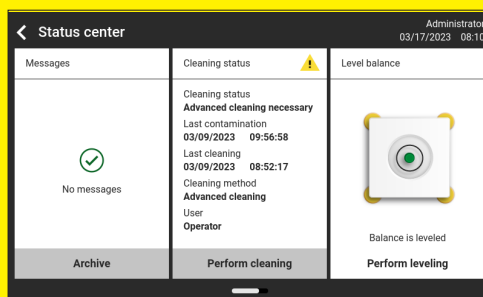
Daily and Advanced Cleaning



Guided Cleaning Process adapted to your balance configuration



Traceable documentation in audit trail



Cleaning status shown in the status center

When Pharma Compliance Matters

Industries depend on strict legal regulations and follow the US or European Pharmacopoeias. The Pharma QApp Package (QP1) of the Cubis® II MCA offers applications which are needed to comply with pharmaceutical-relevant guidelines such as 21 CFR Part 11 and USP 39, Chapter 41. It includes: user management, electronic signature, audit trail, minimum weight, USP advanced, measurement uncertainty, and user calibration.

Pharma (QP1)		License
Q	User management Version: 1.0.1	✓
Q	Electronic signature Version: 1.0.1	✓
Q	Audit trail Version: 1.0.1	✓
Q	Minimum weight Version: 1.0.2	✓
min	USP Advanced Version: 2.4.3	✓
u,w	Measurement uncertainty	✓

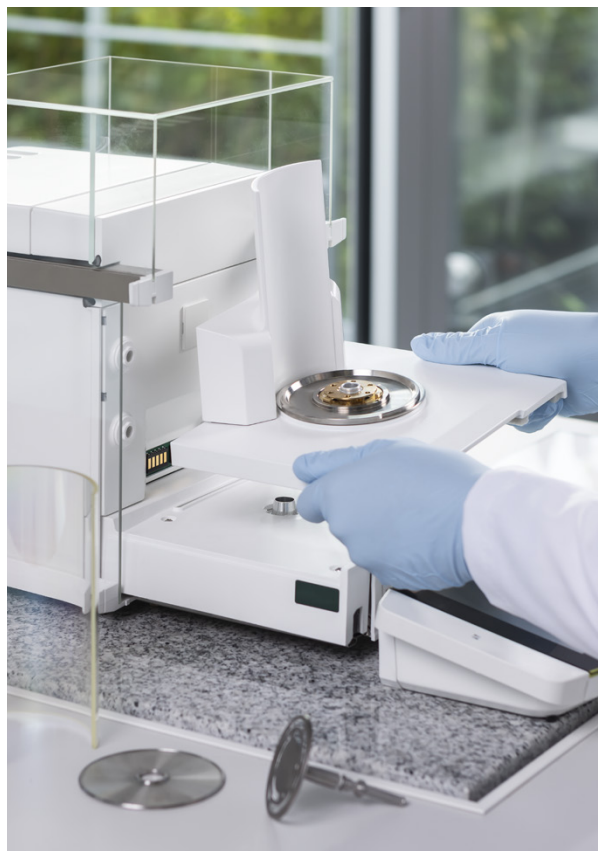
Secure Your Investment with Next Level Upgradeability

Sometimes it is difficult to predict what kind of challenge you will be facing in the future. Is the humidity in your facility too low, or your samples are prone to static charges, but you don't have an ionizer, or wish you had an auto-draftshield but you already purchased your balance?

Wish you had an inner draftshield to protect from drafty conditions? Upgrade these features, anytime

**Upgrade ionizer (QAPP1001),
Motorized draftshield (QAPP1002)**

**Motorized Inner Draftshield (YDS125A)
Click-in Module**



Configure Your Cubis® II Ultra-High Resolution Balance

User interface Weighing module Approval Draft shield QApp Packages Hardware Options

M	C	A	116S	-	3	S	0	0	-	D	QP1	QP2	QP3	HWL
---	---	---	------	---	---	---	---	---	---	---	-----	-----	-----	-----

	Leveling	Approval	Draft shield	Software packages*	Hardware options
Balance type Weighing module**	automatic	S00, S01, CEU, CFR, CCN, OBR, OIN, or OJP	D	QP1, QP2, QP3, QP4, QP10 or QP99	HWL, ION or MDS
High-capacity micro 36S, 36P, 66S, 66P, 116S	■	■	■	■	■
Semi-micro 226S, 125S, 225S, 225P	■	■	■	■	■

* MCA models only

** Weighing modules 116S and 226S available with MCA display only

User Interface	Description	Draft Shield	Description	Approval	Description	Hardware Options	Description
MCA	Advanced user interface	D	Draft shield for High-capacity micro balances with upgradable hardware option	S00	Standard version worldwide	HWL	All hardware licenses
MCE	Essential user interface			S01	Standard only metric units	ION	Ionizer
				CEU	Verified version Europe (except FR)	MDS	Motorized draft shield
				CFR	Verified Version France		
				CCN	Verified Version China		
				OBR	Verifiable Version Brazil		
				OIN	Verifiable Version India		
				OJP	Verifiable Version Japan		



QP1 Pharma The Pharma software application package contains applications concerning the topic compliance with pharmaceutical-relevant guidelines as 21 CFR Part 11 and USP 39, Chapter 41. The Pharma package includes applications such as user management, digital signatures, audit trail, USP minimum weight.



QP2 Advanced Applications The Advanced software application package includes various complex weighing applications incl. evaluation. This includes applications used for density determination, percentage weighing, counting, backweighing, residual dirt analysis, residue on ignition, sieve analysis, filter weighing, checkweighing, formulation, averaging, etc.



QP3 Utilities The Utility software application package contains weighing applications and function extensions such as bootscreen, color scheme, free formula, fiber coarseness, diameter determination, air buoyancy correction, paper weight, statistics and printing of QR | bar codes.



QP4 Connectivity The Connectivity software application package includes applications for data exchange, for example to Windows file server, FTPS, StarLims, etc.



QP10 Hardware The Hardware package offers licensable hardware functions such as motorized draft shield and ionizer at high-capacity micro & semi-micro balances.


QP99 The software application package All inclusive includes 4 different sub-packages for compliance (Pharma (QP1)), complex weighing applications (Advanced (QP2)), weighing applications and helpful tools (Utilities (QP3)) and connectors for data exchange (Connectivity (QP4)).

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
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 contacts, visit
www.sartorius.com