

## Cubis® II: Tu balanza, tu manera



# Cubis® II: La nueva generación de balanzas modulares premium

Para los científicos en I + D o laboratorios analíticos que necesitan los resultados de pesaje de laboratorio más confiables, la plataforma Cubis® II de Sartorius es una cartera de hardware y software de pesaje de laboratorio completamente configurable y de alto rendimiento. El Cubis® II se alineará con sus demandas únicas y requisitos de cumplimiento para ayudarlo a maximizar su eficiencia operativa y resultados experimentales.

La plataforma Cubis® II es el único equilibrio de laboratorio con hardware, software y conectividad totalmente personalizables. Ofrece interfaces de usuario modernas, cumplimiento farmacéutico y GxP, que incluyen manejo de datos, integridad y conectividad de datos, manejo ergonómico de muestras, fácil integración de procesos y comunicación ilimitada con el más alto nivel de exactitud y precisión.



Conectividad



Monitoreo de clima



Integración de flujos de trabajo con QApps



Más información

[www.sartorius.com/cubis-ii](http://www.sartorius.com/cubis-ii)



### Requisitos de Compliance

### Modularidad personalizable

- Miles de configuraciones de hardware y software que admiten preferencias personales.
- Paquetes de software estructurados temáticamente para sus flujos de trabajo optimizados.
- Conectividad perfecta a diversos sistemas de información de laboratorio sin la necesidad de software adicional.
- La flexibilidad a largo plazo garantiza la adaptación del software y accesorios a las necesidades regulatorias y de laboratorio en evolución.

### Inteligencia por diseño

- Las aplicaciones integradas lo guían a través de sus flujos de trabajo.
- Nivelación rápida y fácil garantizada con la función de nivelación automática y motorizada.
- Innovadora tecnología de cabina con sensor de gestos y funciones de memoria.
- Sensor de clima para monitoreo ambiental.
- Ionizador incorporado para la eliminación electrostática automática.
- Centro de estado para notificaciones y alertas de usuarios\*.
- Ultramicro y micro balanzas con tecnología innovadora para un rendimiento de vanguardia.



### Ionizador Integrado

### Cumplimiento de Compliance\*

- Integridad completa de los datos siguiendo los principios de ALCOA+.
- Trazabilidad completa a través de un registro de auditoría avanzado y función de informes para revisiones eficientes.
- Datos listos para envío con selección remota de datos en el navegador web.
- Gestión de usuarios, incluida la gestión centralizada de contraseñas. ("Inicio de sesión único" a través de LDAP), configuración de contraseña de alta seguridad, reglas de inicio de sesión.
- Firma electrónica en cumplimiento con 21 CFR Parte 11.



### Nivelación automática motorizada

### Resultados confiables

- Rendimiento líder
- Precisión de alta velocidad
- Operaciones sin errores
- Integridad de datos completa de extremo a extremo
- Soporte confiable

\*disponible con la interfaz MCA

# Modularidad personalizable – Miles de configuraciones

En 2009, Sartorius ofreció la primera balanza modular con el Cubis®, y hoy continúa esa tradición ofreciendo la Cubis® II, la primera serie de balanzas de laboratorio con un diseño completamente modular:

## Módulos de Pesaje

Elija entre una gama de 45 módulos de pesaje diferentes que mejor se adapten a sus preferencias.

## Cabinas

Elija entre 7 opciones diferentes, dependiendo de su tipo de saldo, cada una con características únicas para admitir el máximo rendimiento del equilibrio y el diseño centrado en el usuario.

## Display

Elija entre la interfaz de usuario avanzada y la esencial, según sus requisitos y aplicaciones de pesaje.

## Software

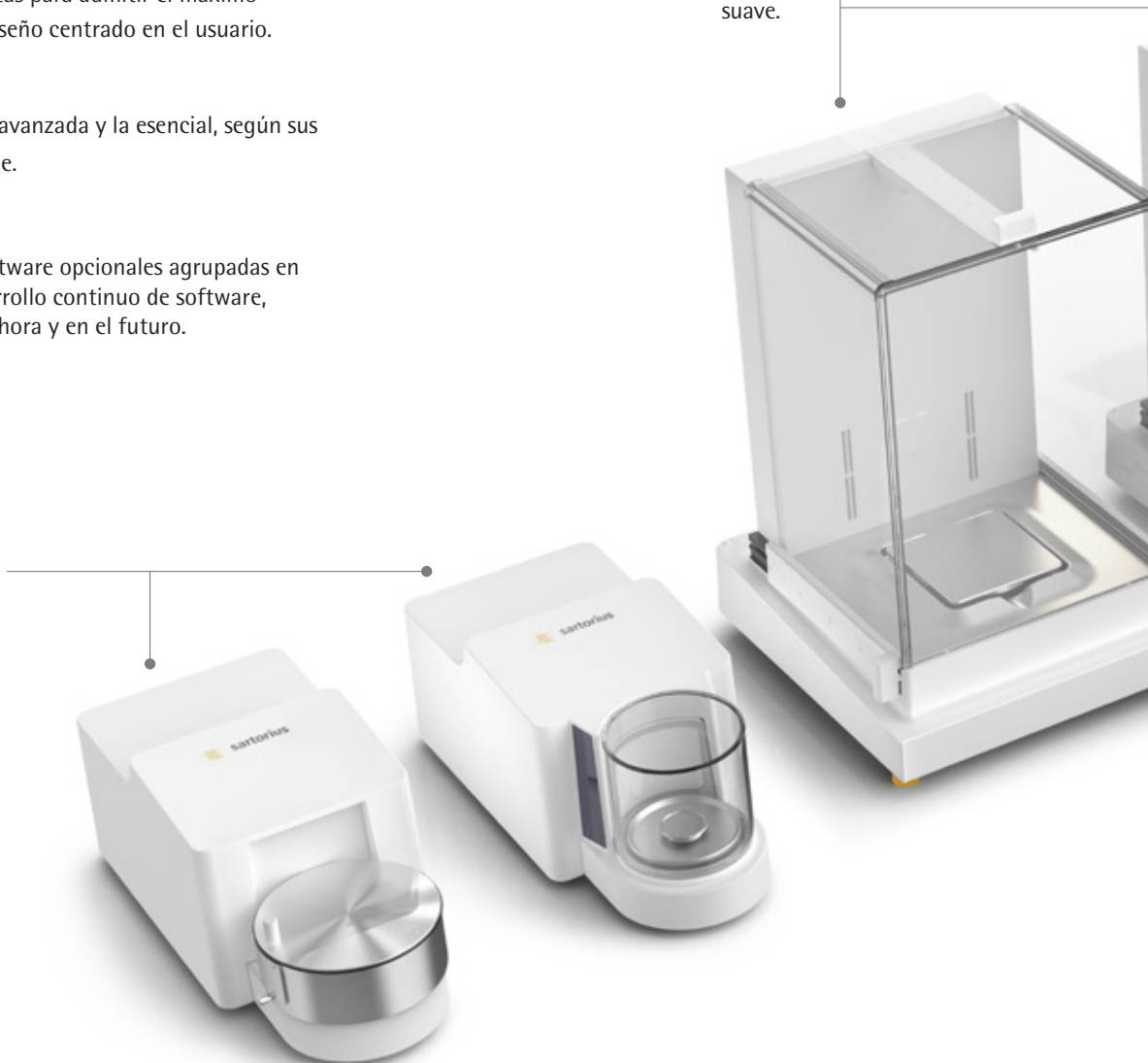
Con hasta 60 aplicaciones de software opcionales agrupadas en paquetes de QApp, y con el desarrollo continuo de software, apoyamos sus flujos de trabajo ahora y en el futuro.

## Balanzas Ultra-Micro & Micro

- Max. Capacidad entre 2,1g y 10,1g | intervalo de escala con un rango de 0,1 µg a 1µg.
- Cabina de tiro de vidrio, redonda, motorizada, automática, con capacidad de aprendizaje, iluminación y sensor de movimiento integrado.
- Protector de tiro de acero inoxidable manual para aplicaciones de pesaje de filtros.

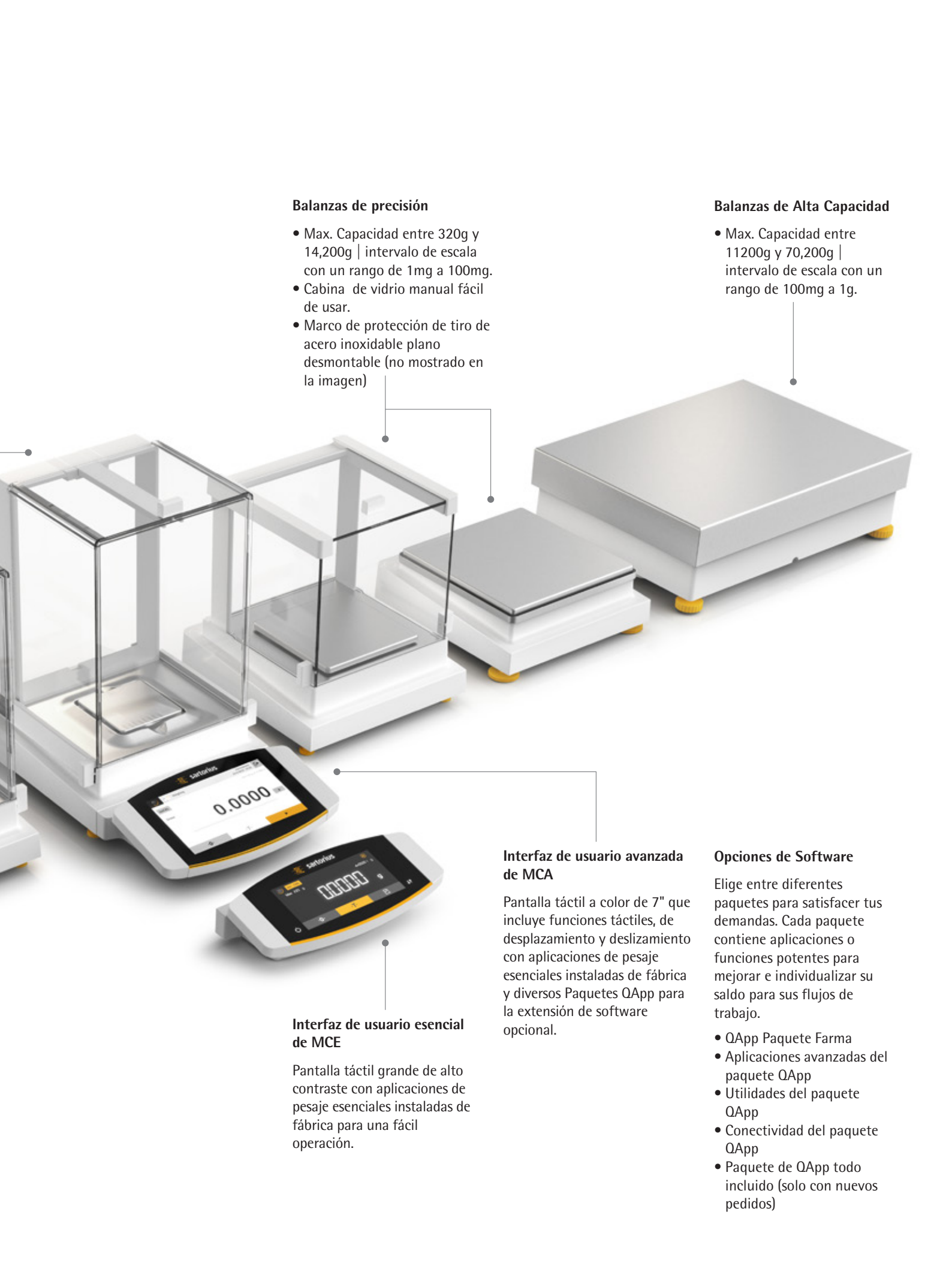
## Balanzas Analíticas & Semi-Micro

- Max. Capacidad entre 120g y 520g | intervalo de escala con un rango de 0,01mg a 0,1mg.
- Cabina de protección de vidrio automática con o sin ionizador integrada para eliminar las cargas electrostáticas que interfieren en las muestras y recipientes de muestras.
- Cabina de protección de vidrio manual con puertas de acción suave.



Más información

[www.sartorius.com/cubis-ii](http://www.sartorius.com/cubis-ii)



### Balanzas de precisión

- Max. Capacidad entre 320g y 14,200g | intervalo de escala con un rango de 1mg a 100mg.
- Cabina de vidrio manual fácil de usar.
- Marco de protección de tiro de acero inoxidable plano desmontable (no mostrado en la imagen)

### Balanzas de Alta Capacidad

- Max. Capacidad entre 11200g y 70,200g | intervalo de escala con un rango de 100mg a 1g.

### Interfaz de usuario esencial de MCE

Pantalla táctil grande de alto contraste con aplicaciones de pesaje esenciales instaladas de fábrica para una fácil operación.

### Interfaz de usuario avanzada de MCA

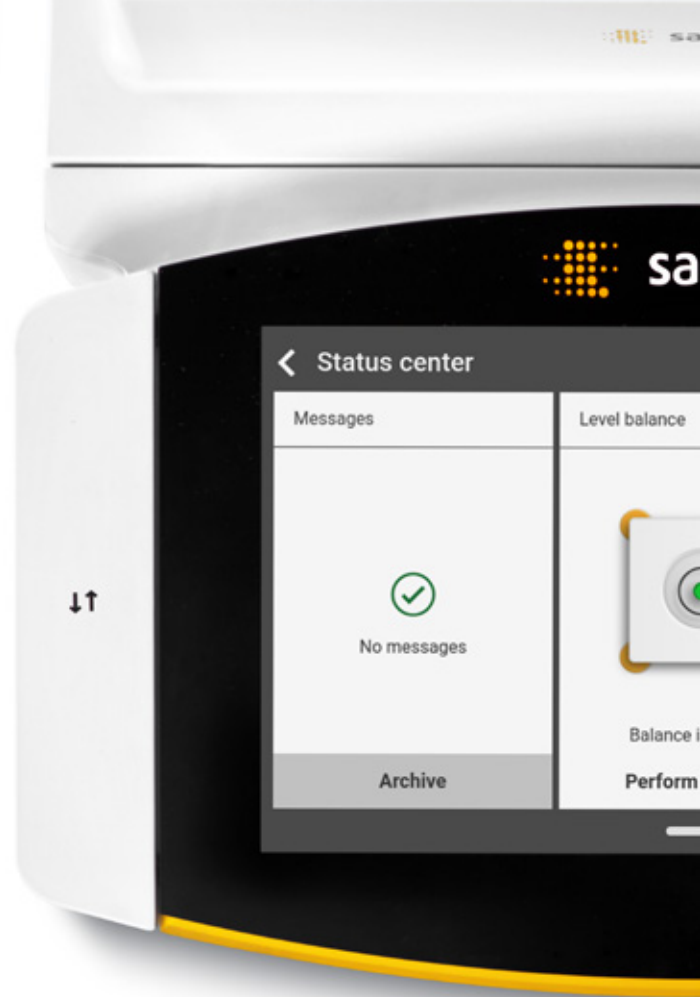
Pantalla táctil a color de 7" que incluye funciones táctiles, de desplazamiento y deslizamiento con aplicaciones de pesaje esenciales instaladas de fábrica y diversos Paquetes QApp para la extensión de software opcional.

### Opciones de Software

Elige entre diferentes paquetes para satisfacer tus demandas. Cada paquete contiene aplicaciones o funciones potentes para mejorar e individualizar su saldo para sus flujos de trabajo.

- QApp Paquete Farma
- Aplicaciones avanzadas del paquete QApp
- Utilidades del paquete QApp
- Conectividad del paquete QApp
- Paquete de QApp todo incluido (solo con nuevos pedidos)

# Intelligence by Design – Adaptable Lab Balance to Your Specific Needs



## Onboard QApp Center

Choose your applications or individual balance functions from QApp packages clustered by different topics. License your selected package and integrate them into the tasks you can run on the balance.

## Automatic, Motorized Leveling

Avoid the burdensome process of preparing your balance for use with the Cubis® II balance's quick and easy, motorized, auto-levelling function.

## Gesture Control

The gesture sensor recognizes four gestures and can be positioned near the balance, directly in the flow of motion at your work space. It allows you to open and close the automatic draft shield, start the ionizer, or send a document to a device.

## Ultra-Micro and Micro Balances

The world's first ultra-micro and micro balances with automatic motorized levelling, motion controlled and illuminated draft shield, and titanium inner parts. Weighing ranges from 0.1 µg to 10.1 g.



### Find out more

For more information, please visit  
[www.sartorius.com/cubis-ii](http://www.sartorius.com/cubis-ii)



### Built-in Ionizer for Electrostatic Elimination

Remove balance stabilization frustrations caused by electrostatic effects with the built-in ionizing technology of the Cubis® II balance.

Electrostatic charges on samples and vessels are removed effectively, just by pressing one key or automatically when the draft shield door closes.



### Climate Monitoring\*,\*\*

The climate sensor monitors influencing variables of the environment and ensures that temperature, pressure, and humidity do not negatively affect your results.



### Status Center Notifications\*

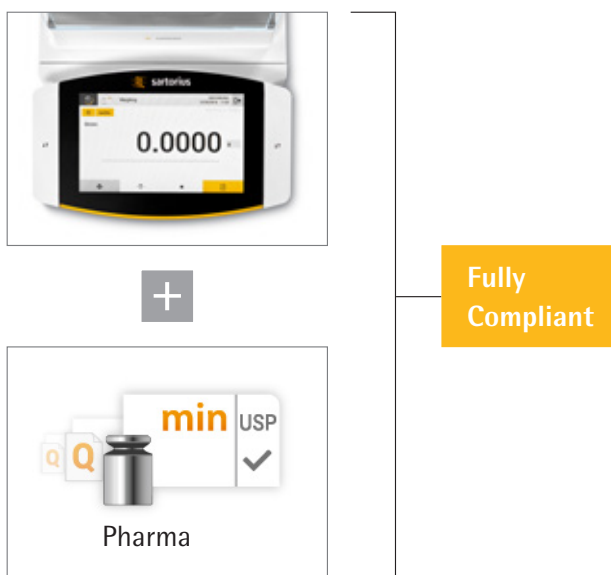
The Status Center shows all information about your balance and environmental conditions, e.g. calibration, leveling, temperature, humidity, air pressure and service, centralized in a dashboard. In case of warnings or errors, you get detailed help and support.



\* available with the MCA interface  
 \*\*available from July 2019

# Confident Compliance – Complete Traceability

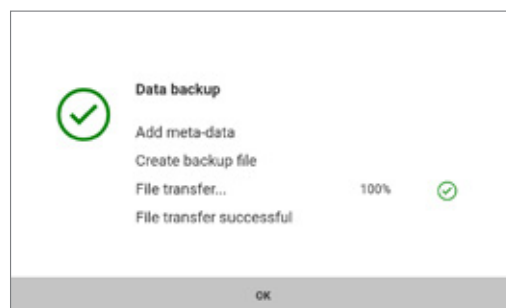
All data generated by the Cubis® II balance follow the key principles ALCOA (+), which defines data integrity standards and relates to GxPs. Therefore, the Cubis® II balance ensures safe and accurate data storage and supports all requirements for FDA directive 21 CFR part 11 and EU Annex 11 compliance.



The Cubis® II balance, with pharma package, contains all the technical controls to support compliance with common regulations. Full compliance can be achieved with additional procedural controls and systems for long-term data storage.

## Backup

Cubis® II can automatically execute time-controlled actions for backup. The system is able to upload data on a file share or export the data to other systems. Archiving is easily possible and fully compliant.



## Safe Data Transfer

An electronic record should be protected against any manipulation. The Cubis® II balance saves all files, together with a calculated MD5 checksum which is stored in the audit trail files and additionally in a separate MD5 file.



### Find out more

For more information, please visit

[www.sartorius.com/cubis-ii](http://www.sartorius.com/cubis-ii)

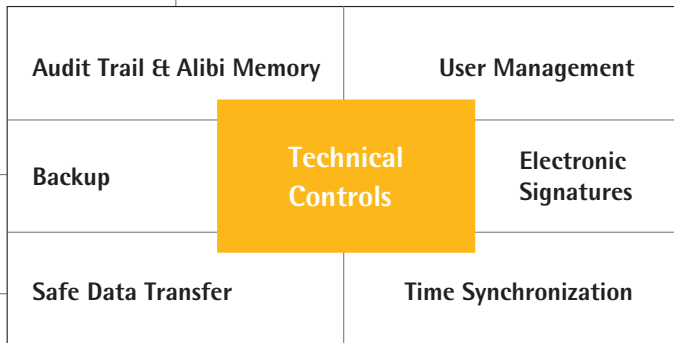


## Audit Trail & Alibi Memory

The Cubis® II balance's audit trail is a tamper-protected time-stamped electronic logfile, that allows reconstruction of events related to the creation, modification, and |or deletion of records. The data can be displayed, filtered, and exported in an easy way.

Additionally, the Cubis® II balance contains an "Alibi Memory". The system automatically stores weighing data in a ring buffer.

Audit trail				
2019-02-27	11:58:12 (UTC)	Administrator	isoCAL	CALIBRATIONADJUSTMENT
2019-02-27	11:55:47 (UTC)	Administrator	Created	TASKMANAGEMENT
2019-02-27	11:54:50 (UTC)	Administrator	Created	QAPPMANAGEMENT
2019-02-27	10:20:41 (UTC)	Administrator	Operation End	CONNECTORMANAGEMENT
2019-02-27	10:05:44 (UTC)	Administrator	Electronic Signature	ACCESSCONTROL
2019-02-27	10:03:52 (UTC)	Administrator	Operation End	CONNECTORMANAGEMENT



## Time Synchronization

The Cubis® II balance supports automatic time synchronization via Network Time Protocol (NTP). An accurate time stamp is elemental for trustable records.

NTP configuration	
NTP operating status	Enabled
Server IP	172.125.14.39

## User Management

The Cubis® II balance provides two options for complete user management with access control.

The local user management can be configured in accordance with your password policy.

The password rules for 21 CFR Part 11 compliance are easily implemented. Moreover, Cubis® II can be integrated in the company's domain to allow single sign-on. In this case the company defined password rules are implemented automatically.

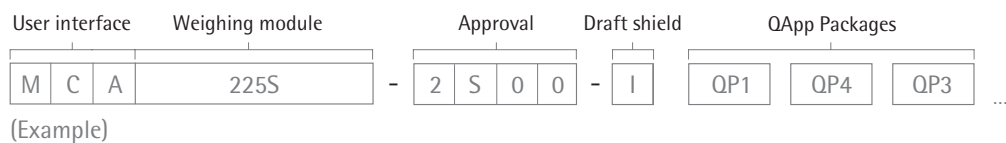
User management	
<b>Administrator</b>	User account for configuration and management
<b>Guest</b>	User account with restricted access
<b>Operator</b>	User account for everyday work
<b>Service</b>	User account for service technicians

## Electronic Signatures

In the Cubis® II balance, an electronic signature can be used to sign the final report for weighing processes. Based on a secure password, the combination of username and password is fully compliant with 21CFR.

Electronic signature	
Sign report	On

# Configure Your Cubis® II Balance



Balance type Weighing module*	Leveling		Approval			Draft shield				Software packages***		
	auto- matic	manual	S00, S01, CEU, CFR, CCN, OBR, OIN, OJP, ORU	A	E	F	I	M	R	U	O	QP1, QP2, QP3, QP4 or QP99
<b>Ultramicro</b> 2.7S	•		•			•		•				•
<b>Micro</b> 10.6S, 6.6S, 3.6P	•		•			•		•				•
<b>Semimicro</b> 225S, 225P, 125S, 125P	•		•	•			•			•		•
<b>Analytical</b> 524S, 524P, 324S, 324P, 224S, 124S	•		•	•			•			•		•
<b>Precision 1 mg</b> 5203S, 5203P, 3203S, 2203S, 2203P, 1203S, 623S, 623P, 323S	•		•	•	•		•		•	•		•
<b>Precision 10 mg</b> 5202S	•		•	•	•		•		•	•		•
<b>Precision 10 mg</b> 8202S, 6202S, 6202P, 4202S, 2202S, 1202S	•		•								•	•
<b>Precision 10 mg</b> 32202P**, 14202S, 14202P, 10202S		•	•								•	•
<b>Precision 100 mg</b> 5201S, 8201S	•		•								•	•
<b>Precision 100 mg</b> 12201S		•	•								•	•
<b>High Capacity</b> 70201S, 50201S, 36201S, 36201P, 20201S, 11201S, 70200S, 36200S		•	•								•	•



## Find out more

For more information, please visit

[www.sartorius.com/cubis-ii](http://www.sartorius.com/cubis-ii)

\*All weighing modules available with MCA or MCE display

\*\*Weighing module 32202P available with approvals S00 or S01 only

\*\*\*Software packages available for balances with MCA display only

User Interface	Description
MCA	Advanced user interface
MCE	Essential user interface

Draft Shield	Description
O	Balance without draft shield
E	Small glass draft shield
R	Stainless steel draft shield frame
U	Manual glass draft shield
A	Motorized draft shield
I	Motorized draft shield + Ionizer
M	Motorized draft shield micro balances
F	Filter draft shield micro balances

Approval	Description
S00	Standard version worldwide
S01	Standard only metric units
CEU	Verified version Europe (except FR)
CFR	Verified Version France
CCN	Verified Version China
OBR	Verifiable Version Brazil
OIN	Verifiable Version India
OJP	Verifiable Version Japan
ORU	Verifiable Version Russia

## QApp Packages



Pharma

**QP1** 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.



Advanced

**QP2** 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.



Utilities

**QP3** 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.



Connectivity

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

**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)).

# Sales and Service Contacts

For further contacts, visit [www.sartorius.com](http://www.sartorius.com)

## Europe

### Germany

Sartorius Lab Instruments  
GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49.551.308.0

### France & Suisse Romande

Sartorius France  
2, rue Antoine Laurent de Lavoisier  
ZA de la Gaudrée  
91410 Dourdan  
Phone +33.1.70.62.50.00

### Austria

Sartorius Austria GmbH  
Modectcenterstrasse 22  
1030 Vienna  
Phone +43.1.7965760.0

### Belgium

Sartorius Belgium N.V.  
Rue Colonel Bourg 105  
1030 Bruxelles  
Phone +32.2.756.06.90

### Finland & Baltics

Sartorius Biohit Liquid Handling Oy  
Laippatie 1  
00880 Helsinki  
Phone +358.9.755.951

### Hungary

Sartorius Hungária Kft.  
Kagyló u. 5.  
2092 Budakeszi  
Phone +3623.457.227

### Ireland

Sartorius Ireland Ltd.  
Unit 41, The Business Centre  
Stadium Business Park  
Ballycoolin Road  
Dublin 11  
Phone +353.1.8089050

### Italy

Sartorius Italy S.r.l.  
Via Torino 3/5  
20814 Varedo (MB)  
Phone +39.0362.5557.11

### Netherlands

Sartorius Netherlands B.V.  
Phone +31.30.60.53.001  
[info.netherlands@sartorius.com](mailto:info.netherlands@sartorius.com)

### Poland

Sartorius Poland sp.z o.o.  
ul. Wrzesinska 70  
62-025 Kostrzyn  
Phone +48.61.6473830

### Russian Federation

LLC "Sartorius RUS"  
Vasilyevsky Island  
5<sup>th</sup> line 70, Lit. A  
199178 St. Petersburg  
Phone +7.812.327.53.27

### 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

### Switzerland

Sartorius Mechatronics Switzerland AG  
Ringstrasse 24a  
8317 Tagelswangen (ZH)  
Phone +41.44.746.50.00

### U.K.

Sartorius UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9QQ  
Phone +44.1372.737159

### Ukraine

LLS "Sartorius RUS"  
Post Box 440 "B"  
01001 Kiev, Ukraine  
Phone +380.44.411.4918

## Americas

### USA

Sartorius Corporation  
5 Orville Drive, Suite 200  
Bohemia, NY 11716  
Phone +1.631.254.4249  
Toll-free +1.800.635.2906

### Argentina

Sartorius Argentina S.A.  
Int. A. Ávalos 4251  
B1605ECS Munro  
Buenos Aires  
Phone +54.11.4721.0505

### Brazil

Sartorius do Brasil Ltda  
Avenida Senador Vergueiro 2962  
São Bernardo do Campo  
CEP 09600-000 - SP- Brasil  
Phone +55.11.4362.8900

### Canada

Sartorius Canada Inc  
1173 North Service Road West, D4  
Oakville, ON L6M 2V9  
Phone +1.905.569.7977  
Toll-Free +1.800.668.4234

### Mexico

Sartorius de México, S.A. de C.V.  
Libramiento Norte de Tepotzotlan s/n,  
Colonia Barrio Tlacateco,  
Municipio de Tepotzotlan,  
Estado de México,  
C.P. 54605  
Phone +52.55.5562.1102  
[leadsmex@sartorius.com](mailto:leadsmex@sartorius.com)

## Asia | Pacific

### Australia

Sartorius Australia Pty. Ltd.  
Unit 5, 7-11 Rodeo Drive  
Dandenong South Vic 3175  
Phone +61.3.8762.1800

### 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

### Hong Kong

Sartorius Hong Kong Ltd.  
Unit 1012, Lu Plaza  
2 Wing Yip Street  
Kwun Tong  
Kowloon, Hong Kong  
Phone +852.2774.2678

### India

Sartorius Weighing India Pvt. Ltd.  
#69/2-69/3, NH 48, Jakkasandra,  
Nelamangala Tq  
562 123 Bangalore, India  
Phone +91.80.4350.5250

### 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

### 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

### Singapore

Sartorius Singapore Pte. Ltd  
10 Science Park Rd  
The Alpha #02-13/14  
Singapore Science Park II  
Singapore 117684  
Phone +65.6872.3966

### 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

### Thailand

Sartorius (Thailand) Co. Ltd.  
129 Rama 9 Road,  
Huaykwang  
Bangkok 10310  
Phone +66.2643.8361-6



◀ [www.sartorius.com](http://www.sartorius.com)