

## The New Quintix. Redefining Standard.



"We found it really simple to use  
the new Quintix right away.  
I don't know of any other balance  
that is this easy to  
operate intuitively."



The new Sartorius Quintix sets new benchmarks in every aspect for standard lab balances: features, technology, ergonomic style, and, above all, the entirely new, user-friendly operating design of the Quintix that helps make workflows much more efficient. What it boils down to is the Quintix is fail-safe: it rules out weighing errors due to incorrect operation and eliminates the time-consuming chore of fine-tuning the settings. Fractions of a second are all it takes for any user to adapt the Quintix to specific requirements and obtain traceable results much faster than ever.

Sartorius teamed up with experienced lab users to develop the user interface of the new Quintix. Lab users' No. 1 priority was to ensure easy, intuitive operation without having read the manual first. They did it! The self-explanatory icons and plain-text prompts on the large touch screen show you all the information you need to know for the procedure – no more, no less. For example, the touch-activated print key will appear only when Quintix has detected that a Sartorius printer is connected. The Quintix also excels in compliance with workplace safety requirements: Even if you are wearing several pairs of lab gloves, you can still correctly use all the functions of the balance.





Quintix comes standard with convenient, built-in applications. These, along with the extensive accessories Sartorius offers, enable you to customize Quintix to a variety of weight measurement tasks. Take specific gravity. With the built-in Density application and the Sartorius Density Kit, you can effortlessly determine the density of solids and liquids.



#### Mixing

Enables you to weigh in various components of a formula quickly with 100% traceability.



#### Components

Conveniently lets you weigh components of a formula into separate containers and retrieve the total amount weighed-in at any time.



#### Statistics

Saves you work when you need to know the standard deviation and other statistics on an entire group of different samples, but don't have the time to do the number-crunching yourself.



#### Checkweighing

Helps you zip through series of samples when you have to quickly determine whether each one is within a specific tolerance range.



#### Density

Takes the effort out of determining the density of a solid, irregularly-shaped object.



#### Weighing

In addition to this app, all Quintix balances feature isoCAL, the fully automatic internal calibration and adjustment function engineered by Sartorius to ensure total accuracy.



#### Conversion

Takes care of the math when you need to convert a weight using a factor, say, to calculate the weight per unit area.



#### Peak Hold

Locks in the reading so you can later view the maximum force released in milliseconds when a switch is activated during an experiment, or the weight of a bulky sample that hides the display while weighing.



#### Unstable Conditions

Delivers rock-steady results when you have to weigh in an extremely unstable environment or measure the weight of a hyper little mouse that just won't sit still.



#### Percentage

Makes comparisons of samples easy when you need to determine their difference in percent from a reference standard.



#### Counting

Gives you the exact count when you want to know how many identical parts, such as tablets, are in a bag.



“ With the apps in Quintix, we can now take care of exotic applications much more easily and efficiently. ”



“ I just transfer weights and other app data directly into an Excel file at a mouse click, without any additional software or bothering with tweaking this really couldn't be more convenient.”



needing  
the settings –

Every advanced analytical laboratory requires GLP-compliant records, as well as fast and secure processing of the data generated. The Plug and Work technology in Quintix saves you considerable time and effort in generating fast and reliable documents.

Quintix substantially simplifies processing data in spreadsheets. How? After connecting the USB port on Quintix to a computer, just open the Microsoft® Office program you need. Quintix delivers data in your choice of format, either text or numbers, enabling effortless and accurate spreadsheet calculations. No additional software or special configuration settings are needed.

The Sartorius GLP-compliant premium printer, YDP30, is automatically detected by Quintix. Thanks to interactive user guidance, Quintix provides help in generating error-free printouts, for instance, by inputting the correct sample ID numbers.



“ We use Quintix non-stop. It's worth  
its weight in gold because it's a  
balance you can always depend on. ”







Repeatable results and linearity are the key prerequisites to obtaining accurate and traceable weighing results. That's why Quintix has been subjected to extensive automated repeatability, linearity and temperature drift tests to guarantee that it meets our high metrological specifications. Inside a virtual "rigorous testing chamber", Quintix test samples undergo more than 1,000,000 load stress tests to ensure that after years of standing up to the rigours of daily use with different weight loads, it will still deliver accurate results just like it did on day one. Quintix is crafted in Germany with

high-grade components and materials. For example, the weigh cells are mounted on a solid, special aluminum alloy base plate, which is extremely impervious to ambient conditions, such as temperature fluctuations and vibration.

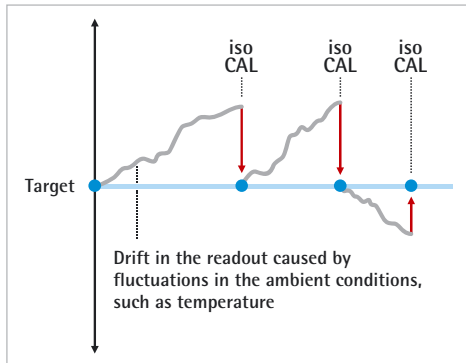
All these tests and features add up to user benefits like the legendary reliability of the Sartorius Quintix, a major advantage it shares with every other Sartorius balance.



Is the linearity still correct even after 1 million cycles? Absolutely! Shown here is the Quintix undergoing an endurance test.



# Peace of Mind with isoCAL



## No Risk Due to Temperature Fluctuations

The slightest fluctuations in temperature have a major impact on the repeatability and the quality of weighing results.

This is why every Quintix comes equipped with the isoCAL internal calibration and adjustment function that ensures consistently accurate results.

Each time isoCAL is performed, the Cal Audit Trail function in Quintix records all data on this procedure so it can be traced for your quality assurance.

## Designs



Design 1



Design 2



Design 3

# Technical Specifications

[illegible]

# Sales and Service Contacts for Industrial Weighing

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

## Europe

### Germany

Sartorius Weighing Technology GmbH  
Weender Landstrasse 94–108  
37075 Goettingen

Phone +49.551.308.0  
Fax +49.551.308.3289

[info.mechatronics@sartorius.com](mailto:info.mechatronics@sartorius.com)  
[www.sartorius.com](http://www.sartorius.com)

Sartorius Mechatronics  
C&D GmbH & Co. KG.  
Am Gut Wolf 11  
52070 Aachen

Phone +49.241.1827.0  
Fax +49.241.1827.213

Sartorius Mechatronics T&H GmbH  
Meiendorfer Strasse 205  
22145 Hamburg

Phone +49.40.67960.303  
Fax +49.40.67960.383

### Austria

Sartorius Mechatronics Austria GmbH  
Franzosengraben 12  
1030 Vienna

Phone +43.1.7965760.0  
Fax +43.1.7965760.24

[info.austria@sartorius.com](mailto:info.austria@sartorius.com)

### Belgium

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

Phone +32.2.756.06.71  
Fax +32.2.253.45.95

[info.belgium@sartorius.com](mailto:info.belgium@sartorius.com)

### France & Suisse Romande

Sartorius Mechatronics France SAS  
Parc d'affaires Technopolis –  
L.P.: 853 Bâtiment SIGMA –  
RDC 3 Avenue du Canada  
91940 Les Ulis, France

Phone +33.1.6919.2100  
Fax +33.1.6920.0922

[internet.france@sartorius.com](mailto:internet.france@sartorius.com)

### Hungary

Sartorius Mechatronics Hungária Kft.  
Kagyló u. 5.  
2092 Budakeszi

Phone +3623.457.227  
Fax +3623.457.147

[mechatronika@sartorius.hu](mailto:mechatronika@sartorius.hu)

### Ireland

Sartorius Mechatronics Ireland Limited  
Unit 41, The Business Centre  
Stadium Business Park  
Ballycoolin Road  
Dublin 11

Phone +353-(0)1-8089050  
Fax +353-(0)1-8089388

[info.ireland@sartorius.com](mailto:info.ireland@sartorius.com)

### Italy

Sartorius Mechatronics Italy S.r.l.  
Viale A. Casati, 4  
20853 Muggiò (MB)

Phone +39.039.46591  
Fax +39.039.465988

[info.italy@sartorius.com](mailto:info.italy@sartorius.com)

### Netherlands

Sartorius Mechatronics  
Netherlands B.V.  
Edisonbaan 24  
3439 MN Nieuwegein

Phone +31.30.6053001  
Fax +31.30.6052917

[weegtechniek.nl@sartorius.com](mailto:weegtechniek.nl@sartorius.com)

### Poland

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

Phone +48.61.6473830  
Fax +48.61.6473839

[info.pl@sartorius.com](mailto:info.pl@sartorius.com)

### Spain

Sartorius Mechatronics Spain S.A.U.  
Offices in Madrid:  
c/ Isabel Colbrand, 10–12, of. 70  
28050 Madrid

Phone +34.91.212.33.67  
Fax +34.91.358.84.85

[spain.weighing@sartorius.com](mailto:spain.weighing@sartorius.com)

### Switzerland

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

Phone +41.44.746.50.00  
Fax +41.44.746.50.50

[mechatronics.switzerland@sartorius.com](mailto:mechatronics.switzerland@sartorius.com)

### U.K.

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

Phone +44.1372.737102  
Fax +44.1372.729927

[uk.customerservice@sartorius.com](mailto:uk.customerservice@sartorius.com)

## America

### USA

Sartorius Mechatronics Corporation  
5 Orville Drive, Suite 200  
Bohemia, New York 11716

Phone +1.631.254.4249  
Toll-free +1.800.635.2906  
Fax +1.631.254.4253

[wt.sales@sartorius.com](mailto:wt.sales@sartorius.com)

### Argentina

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

Phone +54.11.4721.0505  
Fax +54.11.4762.2333

[sartorius@sartorius.com.ar](mailto:sartorius@sartorius.com.ar)

### Brazil

Sartorius do Brasil Ltda  
Av. Dom Pedro I, 241  
Bairro Vila Pires  
Santo André  
São Paulo  
Cep 09110-001

Phone +55.11.4451.6226  
Fax +55.11.4451.4369

[sartorius@sartorius.com.br](mailto:sartorius@sartorius.com.br)

### Canada

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

Phone +1.905.569.7977  
Toll-Free +1.800.668.4234  
Fax +1.905.569.7021

[sales.canada@sartorius.com](mailto:sales.canada@sartorius.com)

### 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

[sartorius@sartomex.com.mx](mailto:sartorius@sartomex.com.mx)

## Asia | Pacific

### 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

[ssil@sartorius.com](mailto:ssil@sartorius.com)

### India

Sartorius Weighing India Pvt Ltd.  
No. 69/2 & 69/3, Jakkasandra,  
NH48, Nelamangala, Kunigal Road,  
Bangalore-562 123

Phone +91.80.4350.5250

[mechatronics.india@sartorius.com](mailto:mechatronics.india@sartorius.com)

### Japan

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

Phone +81.3.3740.5408  
Fax +81.3.3740.5406

[info@sartorius.co.jp](mailto:info@sartorius.co.jp)



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