

The New Secura<sup>®</sup>.  
Safe & Secure Weighing.



Reduce Your Risk with Integrated  
and Intelligent



# Protection Systems

Secura® gives you the security and peace of mind of knowing that you have done everything right. Besides providing highly accurate weighing results and operating convenience, Secura® also features built-in protection systems for complete reliability and regulatory compliance – the safe and secure way.

Real-time guidance prompts for leveling, automatic internal adjustment and 100% traceable, clear documentation with sample and batch identifiers make your lab work more efficient.

First, the new operating concept of Secura® will noticeably ease your daily workload during weighing and, second, its APC function – Advanced Pharma Compliance – will relieve you from tedious and time-consuming documentation and monitoring tasks.



reddot design award  
winner 2013



product  
design award

2013



communication  
design award

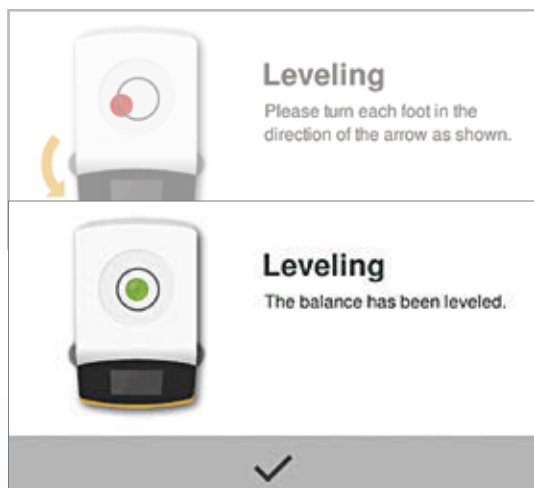
2013 GOLD

A detailed view of the Secura analytical balance scale. It has a white and grey body with a glass-enclosed weighing chamber. Inside the chamber, a small glass vial containing white powder sits on the weighing pan. The brand name 'Secura' is printed in white on a dark grey panel at the bottom of the front door.

Secura





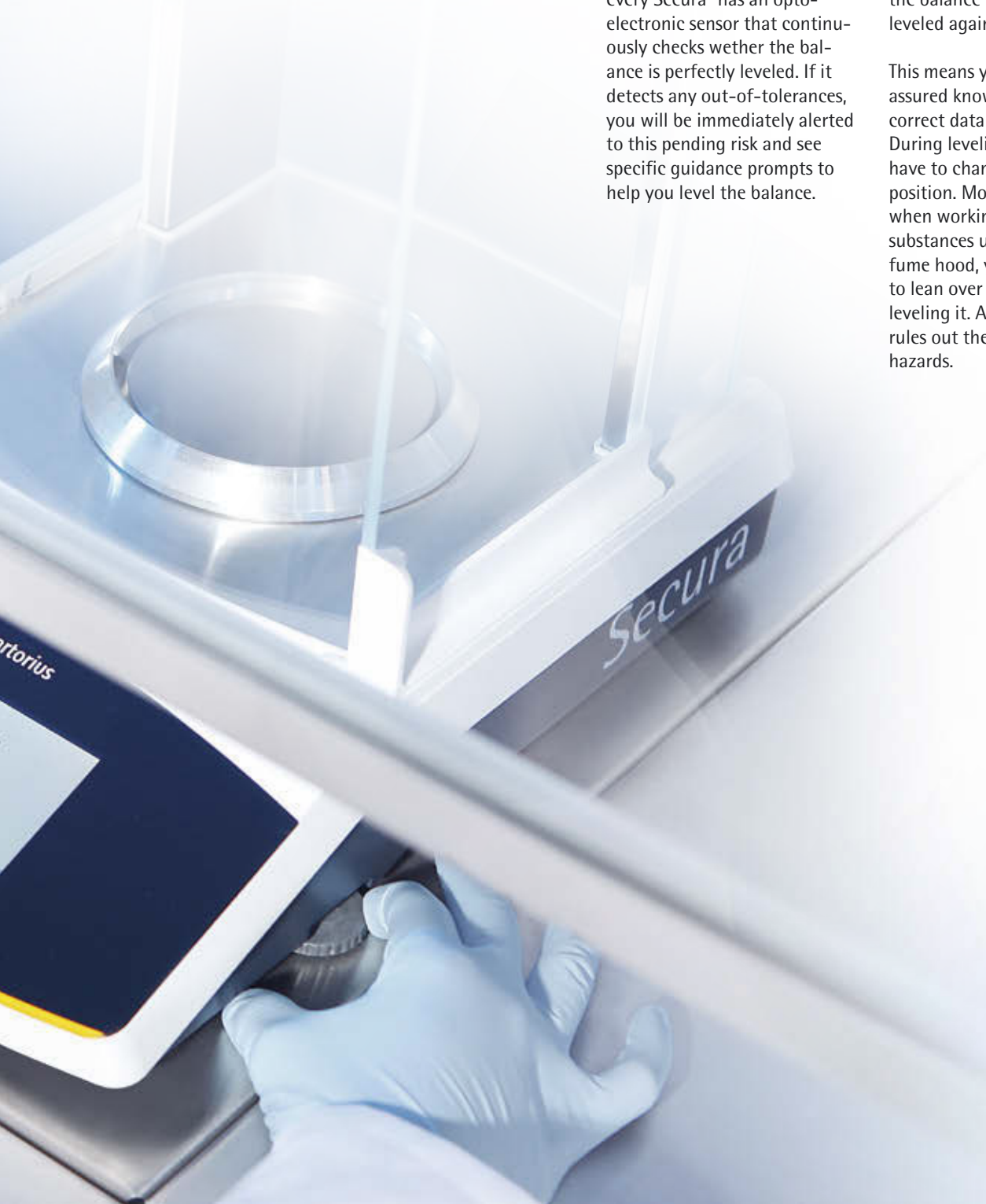


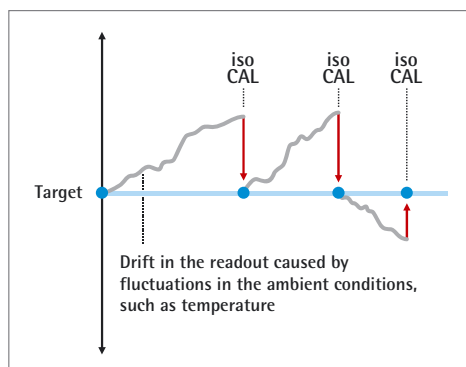
## Always Correctly Levelled with Sartorius LevelControl

Avoid errors caused by an unlevelled balance with real time leveling support - therefore every Secura® has an opto-electronic sensor that continuously checks whether the balance is perfectly leveled. If it detects any out-of-tolerances, you will be immediately alerted to this pending risk and see specific guidance prompts to help you level the balance.

Output of the weighing data over the interface port will be temporarily interrupted until the balance has been correctly leveled again.

This means you can rest assured knowing that only correct data is transmitted. During leveling, you will not have to change your seated position. More importantly, when working with toxic substances under a laboratory fume hood, you will not have to lean over the balance while leveling it. As a result, Secura® rules out the risk of health hazards.





## No Risk Due to Temperature Fluctuations

Temperature fluctuations have an influence on the accuracy of your weighing results.

This is why every Secura® comes equipped with the internal adjustment function isoCAL that ensures consistently accurate results.

Secura® can be configured to prompt you to launch this procedure or to have isoCAL take care of this as soon as an activation threshold is reached.

The Cal Audit Trail function records each isoCAL procedure performed so it can be traced for your quality assurance.



## Stress-Free USP Compliant Measurements

Compliance with the new USP, Chapter 41, requirements (valid from 1st of December, 2013) with absolute process reliability? No problem for the Secura®.

SQmin automatically detects and clearly indicates on the display if you are out of the accepted operating range. In addition, the balance will temporarily block the data transfer to a printer or a device connected to its interface port.

This gives you peace of mind knowing that inadequate data cannot be further processed by accident.

Our specially trained service staff will determine the operation range for you right at the balance's site of installation, program it into your Secura® and attach a corresponding verification seal to the balance.





## Security by Password Protection

Secura® comes with password protection to restrict adjustment of the balance settings only to authorized persons.

If you set a password, all functions that could change the metrological performance specifications are inactivated.

## Interactive User Guidance for Fast GLP- | GMP-Compliant Printouts

If you require traceable documents, Secura® will guide you through the steps for obtaining GLP- | GMP-compliant printouts with proper time and date stamps. The novel user guidance will support you especially for working with traceable sample and batch identification.

Used in conjunction with the new premium laboratory printer YDP30, which is automatically detected by Secura®, the balance generates printouts of the results on standard printer paper or on endless labels.

2013-02-21	14:46
Sartorius	
Mod.	SECURA224-1S
SerNo.	0027703716
BAC:	00-50-02
APC:	01-70-02
Comp1 +	1.7082 g
Comp2 +	1.5017 g
Comp3 +	0.0510 g
T-Comp+	3.2608 g
2013-02-21	14:47
Name:	



## Best Results – Quality Documentation Made Easy

Secura® features a Cal Audit Trail function to document the quality of each weigh-in procedure you perform.

Each combined calibration and adjustment procedure and each leveling process, including any

non-conformities, are recorded by the balance, can be viewed on the touch screen and, if necessary, can be transferred over the balance's interface port to a computer or to a Sartorius laboratory printer.

2012-05-08 [1/5]



Date and Time:	08-May-2012 13:44
Calibration function:	Internal calibration
Start calibration:	Start: isoCAL (Tim)
Calibration deviation:	Dev 0.0002 g
Adjustment function:	Internal adjustment
Adjustment deviation:	Dev 0.0000 g



PREV

NEXT



## Cleanability to Secure the Best Results

Cleanliness is one of the main prerequisites in the laboratory to prevent contamination and ensure work area safety.

That's why Sartorius placed a premium on designing Secura® to enable easy and thorough cleaning.

Sleek, resistant surfaces plus clear contours and few edges ensure that Secura® will help you stay compliant with all hygiene requirements. Even the chemically resistant housing makes cleaning the Secura® a breeze.



## Regulatory Compliance with Expert Support

Do you use your Secura® or other instruments in a regulated area and have to comply with a multitude of requirements, regulations and directives? Relax. Sartorius Lab Products & Services will provide you with the appropriate expert services specifically tailored to ensure compliance with the strict requirements of e.g. the pharmaceutical industry – and not just for balances.



## Qualification – IQ | OQ Documentation

Do you need qualification documents as proof of suitability to be able to use your Secura® in a regulated area? Sartorius Lab Products & Services will support you by providing trained specialists on site. We will

generate the documents you need according to risk-based test plans and test reports for installation and operational qualification.





## Secura® Technical Specifications

Model		224	124	613	513	313	213	6102	5102	3102	2102	1102	612
Weighing capacity	g	220	120	610	510	310	210	6,100	5,100	3,100	2,100	1,100	610
Readability	mg	0.1	0.1	1	1	1	1	10	10	10	10	10	10
Repeatability	mg	0.1	0.1	1	1	1	1	10	10	10	10	10	10
Linearity	mg	0.2	0.2	2	2	2	2	20	20	20	20	20	20
Min. sample weight according to USP, typical	g	0.12	0.12	1.5	1.5	1.5	1.5	12	12	12	12	12	12
Temperature limits for verified models with activated isoCAL		+10°C   +30°C											
Typical stabilization time	s	2	2	1	1	1	1	1	1	1	1	1	1
Weighing pan size	mm	Ø 90	Ø 90	Ø 120	Ø 120	Ø 120	Ø 120	Ø 180	Ø 180	Ø 180	Ø 180	Ø 180	Ø 180
Weighing chamber height	mm	209	209	209	209	209	209	–	–	–	–	–	–
Net weight, approx.	kg (lb)	5.1 (11.2)	5.1 (11.2)	5.1 (11.2)	5.1 (11.2)	5.1 (11.2)	5.1 (11.2)	5.2 (11.5)	5.2 (11.5)	5.2 (11.5)	5.2 (11.5)	5.2 (11.5)	5.2 (11.5)
Dimensions, D × W × H mm (Inches)		360 × 216 × 320 (14.1 × 8.5 × 12.6)						360 × 216 × 95 (14.1 × 8.5 × 3.75)					
Interface		mini USB – Automatic recognition of Sartorius printer models YDP30 and YDP40 – Direct data transfer to Microsoft® application programs – Programmable interval for data output – Choice of SBI, xBPI, table format or text format data transfer protocols											
APC functions (Advanced Pharma Compliance)		– Monitoring if you are within the accepted operating range according to the requirements of the new USP, Chapter 41 – SQmin – Password protection of setup settings – Fully automatic temperature- and time-controlled internal adjustment – isoCAL – Temporary blockage of data transfer to printer or computer when uncertain weighing results are detected, such as when you are out of the accepted operating range according to USP, the balance is not levelled or isoCAL adjustment needs to be performed – Storage of all calibration, adjustment and levelling procedures data – Cal Audit Trail											
Leveling		Intelligent, optoelectronic level sensor with alarm function and interactive user guidance for reliable and safe leveling											
Display		Touchscreen with Sartorius graphical user interface											
Built-in application programs		Special lab applications such as Mixing, Components, Statistics and Conversion; plus standard lab applications such as Weighing, Density, Percentage, Checkweighing, Peak Hold, Counting and Unstable Conditions											
Accessories		See Secura® list of accessories											



◀ www.sartorius.com

**Germany**  
 Sartorius Lab Instruments GmbH & Co. KG  
 Weender Landstrasse 94–108  
 37075 Goettingen  
 Phone +49.551.308.0  
 Fax +49.551.308.3289