



Entris® II

Laboratory Balances

Simplifying Progress

SARTORIUS



Entris® II – Best Value in Its Class for Basic Weighing Tasks

Every Sartorius Balance Offers Quality, Value and Consistency

No matter what you're weighing, the new Entris® II is always the right choice. Offering unrivalled value at a budget price, this is the only balance in its class featuring isoCAL, LED touch technology and 12 built-in applications. Backed by almost 150 years of German engineering expertise, and available in 40 different models, you'll easily find an Entris® II balance that exactly meets your specific weighing needs.

 Find out more
www.sartorius.com



Quality, Value and Consistency

For Accurate Weighing Results You Can Trust Time After Time

Highest Quality Standards Engineered in Germany

- Highly accurate results year in and year out—guaranteed via monolithic weigh cell technology invented by Sartorius*
- Fastest stabilization time in its class—using state-of-the-art weighing sensors
- Best repeatability
 - with rectangular weighing pan
- Overload protection
 - rugged design weighs up to a pre-set amount
- Guaranteed reliability—with the self-test “@start”
- Made in Germany

Easy Clean for Operational Efficiency and Durability

- High chemical resistance—ensured using parts made from hard-wearing polybutylene terephthalate (PBT), stainless steel and glass
- Prevents cross contamination—with wipe clean design and easy-to-remove parts

Effective Draft Shield

- Minimizes weighing errors caused by electrostatically charged samples—using specially coated glass parts

Innovations in This Weighing Class

For Efficient, User-Friendly Operation

isoCAL (Internal Calibration and Adjustment) for Total Assurance of Accurate Weighing Results

- Optimal accuracy and operating convenience—using fully automated internal temperature and time-controlled calibration and adjustment feature, unique in this weighing class
- Assures (SOP) compliant operation—with self-notification if calibration is outside the normal range

Hybrid Screen for Excellent Readability and Use

- Combines intuitive, wear-resistant LED and touch technology—provides an easy and clear structured user interface

Plug-And-Play Convenience

- Automatically detects Sartorius accessories (e.g., printer, second display)
- Real “PC-direct feature” for easy connection to a PC to transfer weighing data directly into spreadsheets or documents such as Microsoft® Excel or Word

Data Output for Dynamic Weighing Applications

- Configurable time interval for data output

Easily Adapts to Your Ambient Conditions

- With just one click on the home screen



* may differ for a few models

Universal Application

For Future Proof Assurance

The Right Model for Each and Every Task

- Offering a range of 40 models (the largest portfolio in its class), starting from 60 g up to 8,200 g and a readability from 0.1 mg to 1 g—finding the right model to suit any weighing application is simple

12 Built-In Applications with GLP | GMP Compliant Printout | Data Output

- Weighing | Dosing
- Counting
- Percentage weighing
- Mixing | Net Total
- Components | Totalizing
- Animal weighing
- Calculation | Free Factor
- Density determination
- Statistics
- Peakhold
- Checkweighing
- Mass unit conversion
- Underfloor weighing feature for bigger samples

Weighing Chamber (Where Applicable)

- Highest usable weighing chamber height for easy handling of even large containers
- Convenient access to weighing pan
- Easy-to-(re)move top and side sliding doors
- Draft shield can be removed completely

8 Built-In Languages for International Operation

- English, German, French, Italian, Spanish, Portuguese, Russian, Polish

Modern, Up-To-Date Connectivity Methods

- Future-proof USB Type C interface Industry-proof RS232 9-pin interface
- Backwards compatibility (by using RS232 port)

Password Protection for Secure Operation

- Protects the balance against unintentional changes

Find out more
www.sartorius.com



Entris® II Technical Specifications



Model	Weighing Capacity	Readability	Repeatability, Typical	Stabilization Time, Typical	Weighing Pan Size	Weighing Chamber Height*	Dimensions (W x D x H)	Calibration and Adjustment	
	[g]	[mg]	[mg]	[s]	[mm]	[mm]	[mm]	Internal	External
Analytical Balances									
BCE64i-1x	60	0.1	0.1	≤1.5	Ø 90	240	219×317×345	<input type="checkbox"/>	<input type="checkbox"/>
BCE64-1x	60	0.1	0.1	≤1.5	Ø 90	240	219×317×345		<input type="checkbox"/>
BCE124i-1x	120	0.1	0.1	≤1.5	Ø 90	240	219×317×345	<input type="checkbox"/>	<input type="checkbox"/>
BCE124-1x	120	0.1	0.1	≤1.5	Ø 90	240	219×317×345		<input type="checkbox"/>
BCE224i-1x	220	0.1	0.1	≤1.5	Ø 90	240	219×317×345	<input type="checkbox"/>	<input type="checkbox"/>
BCE224-1x	220	0.1	0.1	≤1.5	Ø 90	240	219×317×345		<input type="checkbox"/>
Precision Balances									
BCE223i-1x	220	1	1	≤1.0	Ø 120	240	219 x 317 x 345	<input type="checkbox"/>	<input type="checkbox"/>
BCE223-1x	220	1	1	≤1.0	Ø 120	240	219 x 317 x 345		<input type="checkbox"/>
BCE323i-1x	320	1	1	≤1.0	Ø 120	240	219 x 317 x 345	<input type="checkbox"/>	<input type="checkbox"/>
BCE323-1x	320	1	1	≤1.0	Ø 120	240	219 x 317 x 345		<input type="checkbox"/>
BCE423i-1x	420	1	1	≤1.0	Ø 120	240	219 x 317 x 345	<input type="checkbox"/>	<input type="checkbox"/>
BCE423-1x	420	1	1	≤1.0	Ø 120	240	219 x 317 x 345		<input type="checkbox"/>
BCE623i-1x	620	1	1	≤1.0	Ø 120	240	219 x 317 x 345	<input type="checkbox"/>	<input type="checkbox"/>
BCE623-1x	620	1	1	≤1.0	Ø 120	240	219 x 317 x 345		<input type="checkbox"/>
BCE653i-1x	650	1	1	≤1.0	Ø 120	50	219 x 317 x 145	<input type="checkbox"/>	<input type="checkbox"/>
BCE653-1x	650	1	1	≤1.0	Ø 120	50	219 x 317 x 145		<input type="checkbox"/>
BCE622i-1x	620	10	10	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE622-1x	620	10	10	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE822i-1x	820	10	10	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE822-1x	820	10	10	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE1202i-1x	1,200	10	10	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE1202-1x	1,200	10	10	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE2202i-1x	2,200	10	10	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE2202-1x	2,200	10	10	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE3202i-1x	3,200	10	10	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE3202-1x	3,200	10	10	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE4202i-1x	4,200	10	10	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE4202-1x	4,200	10	10	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE6202i-1x	6,200	10	10	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE6202-1x	6,200	10	10	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE2201i-1x	2,200	100	100	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE2201-1x	2,200	100	100	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE5201i-1x	5,200	100	100	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE5201-1x	5,200	100	100	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE8201i-1x	8,200	100	100	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE8201-1x	8,200	100	100	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE6200i-1x	6,200	1,000	1,000	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE6200-1x	6,200	1,000	1,000	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>
BCE8200i-1x	8,200	1,000	1,000	≤0.9	182 x 182	-	219 x 317 x 94	<input type="checkbox"/>	<input type="checkbox"/>
BCE8200-1x	8,200	1,000	1,000	≤0.9	182 x 182	-	219 x 317 x 94		<input type="checkbox"/>

* upper edge of the weighing pan to the lower edge of the upper draft shield panel
x = term for country-specific models

Sales and Service Contacts

For further contacts, visit
www.sartorius.com

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