

## Entris® II Laboratory Balances



turning science **into solutions**

## 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  
For more information, please visit  
[www.sartorius.com](http://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



Find out more  
For more information, please visit  
[www.sartorius.com](http://www.sartorius.com)

# 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
- and 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



## 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	
							[mm]	[mm]	Internal	External
<b>Analytical balances</b>										
BCE64i-1x	60	0.1	0.1	≤1,5	Ø 90	240	219x317x345	•	•	
BCE64-1x	60	0.1	0.1	≤1,5	Ø 90	240	219x317x345	-	•	
BCE124i-1x	120	0.1	0.1	≤1,5	Ø 90	240	219x317x345	•	•	
BCE124-1x	120	0.1	0.1	≤1,5	Ø 90	240	219x317x345	-	•	
BCE224i-1x	220	0.1	0.1	≤1,5	Ø 90	240	219x317x345	•	•	
BCE224-1x	220	0.1	0.1	≤1,5	Ø 90	240	219x317x345	-	•	
<b>Precision balances</b>										
BCE223i-1x	220	1	1	≤1,0	Ø 120	240	219x317x345	•	•	
BCE223-1x	220	1	1	≤1,0	Ø 120	240	219x317x345	-	•	
BCE323i-1x	320	1	1	≤1,0	Ø 120	240	219x317x345	•	•	
BCE323-1x	320	1	1	≤1,0	Ø 120	240	219x317x345	-	•	
BCE423i-1x	420	1	1	≤1,0	Ø 120	240	219x317x345	•	•	
BCE423-1x	420	1	1	≤1,0	Ø 120	240	219x317x345	-	•	
BCE623i-1x	620	1	1	≤1,0	Ø 120	240	219x317x345	•	•	
BCE623-1x	620	1	1	≤1,0	Ø 120	240	219x317x345	-	•	
BCE653i-1x	650	1	1	≤1,0	Ø 120	50	219x317x145	•	•	
BCE653-1x	650	1	1	≤1,0	Ø 120	50	219x317x145	-	•	
BCE622i-1x	620	10	10	≤0,9	182x182	-	219x317x94	•	•	
BCE622-1x	620	10	10	≤0,9	182x182	-	219x317x94	-	•	
BCE822i-1x	820	10	10	≤0,9	182x182	-	219x317x94	•	•	
BCE822-1x	820	10	10	≤0,9	182x182	-	219x317x94	-	•	
BCE1202i-1x	1,200	10	10	≤0,9	182x182	-	219x317x94	•	•	
BCE1202-1x	1,200	10	10	≤0,9	182x182	-	219x317x94	-	•	
BCE2202i-1x	2,200	10	10	≤0,9	182x182	-	219x317x94	•	•	
BCE2202-1x	2,200	10	10	≤0,9	182x182	-	219x317x94	-	•	
BCE3202i-1x	3,200	10	10	≤0,9	182x182	-	219x317x94	•	•	
BCE3202-1x	3,200	10	10	≤0,9	182x182	-	219x317x94	-	•	
BCE4202i-1x	4,200	10	10	≤0,9	182x182	-	219x317x94	•	•	
BCE4202-1x	4,200	10	10	≤0,9	182x182	-	219x317x94	-	•	
BCE6202i-1x	6,200	10	10	≤0,9	182x182	-	219x317x94	•	•	
BCE6202-1x	6,200	10	10	≤0,9	182x182	-	219x317x94	-	•	
BCE2201i-1x	2,200	100	100	≤0,9	182x182	-	219x317x94	•	•	
BCE2201-1x	2,200	100	100	≤0,9	182x182	-	219x317x94	-	•	
BCE5201i-1x	5,200	100	100	≤0,9	182x182	-	219x317x94	•	•	
BCE5201-1x	5,200	100	100	≤0,9	182x182	-	219x317x94	-	•	
BCE8201i-1x	8,200	100	100	≤0,9	182x182	-	219x317x94	•	•	
BCE8201-1x	8,200	100	100	≤0,9	182x182	-	219x317x94	-	•	
BCE6200i-1x	6,200	1,000	1,000	≤0,9	182x182	-	219x317x94	•	•	
BCE6200-1x	6,200	1,000	1,000	≤0,9	182x182	-	219x317x94	-	•	
BCE8200i-1x	8,200	1,000	1,000	≤0,9	182x182	-	219x317x94	•	•	
BCE8200-1x	8,200	1,000	1,000	≤0,9	182x182	-	219x317x94	-	•	

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



Find out more  
For more information, please visit  
[www.sartorius.com](http://www.sartorius.com)

# 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.  
Kagyó 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  
565 Johnson Avenue  
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 Tepetzotlan s/n,  
Colonia Barrio Tlacateco,  
Municipio de Tepetzotlan,  
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

For more information, view our Sartorius catalog  
and price list or visit our website.



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