

SARTORIUS

Entris[®] II NTEP-Certified Balances Legal for Trade

Best Value in its Class for Basic Weighing Tasks

100% NTEP-Compliant Weighing

No matter what you're weighing, the new Entris[®] II balance is always the right choice. Offering unrivalled value and backed by more than 150 years of German engineering expertise, the Entris[®] II offers a wide range of models, so you can find the balance that meets your specific weighing needs.

Our Entris[®] II Essential line is the only balance in its class offering LED touch technology and 12 built-in applications.

Features & Benefits

- Highly accurate results—guaranteed via monolithic weigh cell technology invented by Sartorius
- Fastest stabilization time in its class—using state-of-the-art weighing sensors
- High chemical resistance—ensured using parts made from hard-wearing polybutylene terephthalate (PBT), stainless steel and glass



100% Legal for Trade

Key Features

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

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



Easy clean for operational efficiency and durability



Effective draft shield



Plug-and-play convenience

Model Information

Model	Weighing capacity	Readability	Repeatability, typical	Stabilization time, typical	Weighing pan size	Weighing chamber height*	Dimensions (W × D × H)
	[g]	[mg]	[mg]	[s]	[mm]	[mm]	[mm]
BCE124-1NUS	120	0.1	0.1	≤1.5	Ø 90	240	219×317×345
BCE224-1NUS	220	0.1	0.1	≤1.5	Ø 90	240	219×317×345
BCE323-1NUS	320	1	1	≤1.0	Ø 120	240	219×317×345
BCE623-1NUS	620	1	1	≤1.0	Ø 120	240	219×317×345
BCE653-1NUS	650	1	1	≤1.0	Ø 120	50	219×317×145
BCE622-1NUS	620	10	10	≤0.9	182×182		219×317×94
BCE1202-1NUS	1,200	10	10	≤0.9	182×182		219×317×94
BCE3202-1NUS	3,200	10	10	≤0.9	182×182		219×317×94
BCE4202-1NUS	4,200	10	10	≤0.9	182×182		219×317×94
BCE6202-1NUS	6,200	10	10	≤0.9	182×182		219×317×94
BCE5201-1NUS	5,200	100	100	≤0.9	182×182		219×317×94
BCE8201-1NUS	8,200	100	100	≤0.9	182×182		219×317×94

* upper edge of the weighing pan to the lower edge of the upper draft shield panel
 NTEP Approval no. 22-017
 Measurement Canada no. AM-6176

Please contact your Sales Representative or any of our authorized Partners for more information.

Find out more:
sartorius.com