

The New Entris[®].
Just Weighing.



The New Entris®. Just Weighing.

Get the reliability and quality of a Sartorius balance, but only pay for the features you really need.

The new Entris® has been specially designed to deliver effective and reliable weighing results in your daily work. The easy and clearly structured user interface, the logical key assignment and excellent readability are ideal features for error-free operation.

Easy Cleaning

The top and side sliding doors and easy-to-remove stainless steel parts of the draft shield chamber make Entris® convenient for thorough cleaning.



Internal Adjustment

Each model is also available with an internal, motorized adjustment function for optimal weighing accuracy and operating convenience.





High Accuracy

The monolithic weigh cell ensures long-term high measurement accuracy and features the legendary reliability and quality of a Sartorius balance.



Basic Applications

Entris® supports you in your daily work with integrated basic applications: animal weighing, density determination, counting, conversion and weighing in percent.



User-friendly Operation

The backlit and high-contrast display ensures excellent readability. Easy navigation with function keys and the simple-to-read level indicator on the front enable user-friendly and fast operation.



Entris® Technical Specifications

Model	Weighing capacity (g)	Readability (mg)	Repeatability (mg)	Linearity (mg)	Typical stabilization time (s)	Weighing pan size (mm)	Dimensions D×W×H (mm)	Calibration and adjustment
Analytical balances								
ENTRIS224i-1S	220	0.1	0.1	0.2	2.5	Ø 90	230×303×330	Internal
ENTRIS224-1S	220	0.1	0.1	0.2	2.5	Ø 90	230×303×330	External
ENTRIS124i-1S	120	0.1	0.1	0.2	2.5	Ø 90	230×303×330	Internal
ENTRIS124-1S	120	0.1	0.1	0.2	2.5	Ø 90	230×303×330	External
ENTRIS64i-1S	60	0.1	0.1	0.2	2.5	Ø 90	230×303×330	Internal
ENTRIS64-1S	60	0.1	0.1	0.2	2.5	Ø 90	230×303×330	External
Precision balances								
ENTRIS623i-1S	620	1	1	2	1.0	Ø 115	230×303×136	Internal
ENTRIS623-1S	620	1	1	2	1.0	Ø 115	230×303×136	External
ENTRIS423i-1S	420	1	1	2	1.0	Ø 115	230×303×136	Internal
ENTRIS423-1S	420	1	1	2	1.0	Ø 115	230×303×136	External
ENTRIS323i-1S	320	1	1	2	1.1	Ø 115	230×303×136	Internal
ENTRIS323-1S	320	1	1	2	1.1	Ø 115	230×303×136	External
ENTRIS153i-1S	150	1	1	2	1.3	Ø 115	230×303×136	Internal
ENTRIS153-1S	150	1	1	2	1.3	Ø 115	230×303×136	External
ENTRIS6202i-1S	6,200	10	10	30	1.5	180×180	230×303×91	Internal
ENTRIS6202-1S	6,200	10	10	30	1.5	180×180	230×303×91	External
ENTRIS4202i-1S	4,200	10	10	30	1.5	180×180	230×303×91	Internal
ENTRIS4202-1S	4,200	10	10	30	1.5	180×180	230×303×91	External
ENTRIS3202i-1S	3,200	10	10	30	1.5	180×180	230×303×91	Internal
ENTRIS3202-1S	3,200	10	10	30	1.5	180×180	230×303×91	External
ENTRIS2202i-1S	2,200	10	10	30	1.5	180×180	230×303×91	Internal
ENTRIS2202-1S	2,200	10	10	30	1.5	180×180	230×303×91	External
ENTRIS822i-1S	820	10	10	30	1.5	Ø 150	230×303×87	Internal
ENTRIS822-1S	820	10	10	30	1.5	Ø 150	230×303×87	External
ENTRIS8201i-1S	8,200	100	100	300	1.5	180×180	230×303×91	Internal
ENTRIS8201-1S	8,200	100	100	300	1.5	180×180	230×303×91	External
ENTRIS5201i-1S	5,200	100	100	300	1.5	180×180	230×303×91	Internal
ENTRIS5201-1S	5,200	100	100	300	1.5	180×180	230×303×91	External
ENTRIS2201i-1S	2,200	100	100	300	1.5	180×180	230×303×91	Internal
ENTRIS2201-1S	2,200	100	100	300	1.5	180×180	230×303×91	External

For accessories, please see the Entris® 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