

Entris[®]



turning science **into solutions**

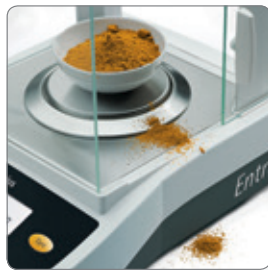
Entris[®]. Just Weighing.

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

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.





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	
								Internal	External
Analytical balances									
ENTRIS224i-1S	220	0.1	0.1	0.2	2.5	∅ 90	303×230×330	•	•
ENTRIS224-1S	220	0.1	0.1	0.2	2.5	∅ 90	303×230×330	–	•
ENTRIS124i-1S	120	0.1	0.1	0.2	2.5	∅ 90	303×230×330	•	•
ENTRIS124-1S	120	0.1	0.1	0.2	2.5	∅ 90	303×230×330	–	•
ENTRIS64i-1S	60	0.1	0.1	0.2	2.5	∅ 90	303×230×330	•	•
ENTRIS64-1S	60	0.1	0.1	0.2	2.5	∅ 90	303×230×330	–	•
Precision balances									
ENTRIS623i-1S	620	1	1	2	1.0	∅ 115	303×230×136	•	•
ENTRIS623-1S	620	1	1	2	1.0	∅ 115	303×230×136	–	•
ENTRIS423i-1S	420	1	1	2	1.0	∅ 115	303×230×136	•	•
ENTRIS423-1S	420	1	1	2	1.0	∅ 115	303×230×136	–	•
ENTRIS323i-1S	320	1	1	2	1.1	∅ 115	303×230×136	•	•
ENTRIS323-1S	320	1	1	2	1.1	∅ 115	303×230×136	–	•
ENTRIS153i-1S	150	1	1	2	1.3	∅ 115	303×230×136	•	•
ENTRIS153-1S	150	1	1	2	1.3	∅ 115	303×230×136	–	•
ENTRIS6202i-1S	6,200	10	10	30	1.5	180×180	303×230×91	•	•
ENTRIS6202-1S	6,200	10	10	30	1.5	180×180	303×230×91	–	•
ENTRIS4202i-1S	4,200	10	10	30	1.5	180×180	303×230×91	•	•
ENTRIS4202-1S	4,200	10	10	30	1.5	180×180	303×230×91	–	•
ENTRIS3202i-1S	3,200	10	10	30	1.5	180×180	303×230×91	•	•
ENTRIS3202-1S	3,200	10	10	30	1.5	180×180	303×230×91	–	•
ENTRIS2202i-1S	2,200	10	10	30	1.5	180×180	303×230×91	•	•
ENTRIS2202-1S	2,200	10	10	30	1.5	180×180	303×230×91	–	•
ENTRIS822i-1S	820	10	10	30	1.5	∅ 150	303×230×87	•	•
ENTRIS822-1S	820	10	10	30	1.5	∅ 150	303×230×87	–	•
ENTRIS8201i-1S	8,200	100	100	300	1.5	180×180	303×230×91	•	•
ENTRIS8201-1S	8,200	100	100	300	1.5	180×180	303×230×91	–	•
ENTRIS5201i-1S	5,200	100	100	300	1.5	180×180	303×230×91	•	•
ENTRIS5201-1S	5,200	100	100	300	1.5	180×180	303×230×91	–	•
ENTRIS2201i-1S	2,200	100	100	300	1.5	180×180	303×230×91	•	•
ENTRIS2201-1S	2,200	100	100	300	1.5	180×180	303×230×91	–	•

For accessories, please see the Entris® list of accessories.



◀ www.sartorius.com

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716

Phone +1.631.254.4249
Toll-free +1.800.635.2906
Fax +1.631.254.4253

Canada

Sartorius Canada Inc.
2179 Dunwin Drive #4
Mississauga, ON L5L 1X2

Phone +1.905.569.7977
Toll-Free +1.800.668.4234
Fax +1.905.569.7021

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289