

# High Resolution Weighing Cells

Integrators & Original  
Equipment Manufacturers

Simplifying Progress

**SARTORIUS**

# For Fast, Robust and Reliable Integration

Sartorius offers a wide range of weigh cells that can be integrated into any machinery or instrumentation, where capturing accurate and reproduceable weighing data quickly is important.

## Wide Portfolio of Models

Sartorius manufactures each load cell using state-of-the-art techniques and precision machining from a single block of aluminum. Measurement is based on electromagnetic force compensation, which provides the highest level of weighing accuracy and reproducibility during extensive operating cycles.

All Sartorius weigh cells are designed for diverse applications in production and testing processes: dosing, dispensing, counting, statistical quality control and check weighing.

The following range is available:

- Capacities up to 8.2 kg
- Resolutions from 0.01 g to 1  $\mu$ g
- Latest electromagnetic force compensation weighing technology
- Choice of aluminum or stainless-steel housings
- Separate electronics and weigh cell for optimized performance

## Easy installation and integration

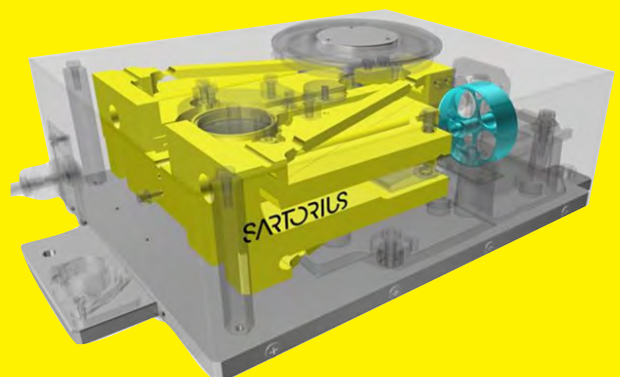
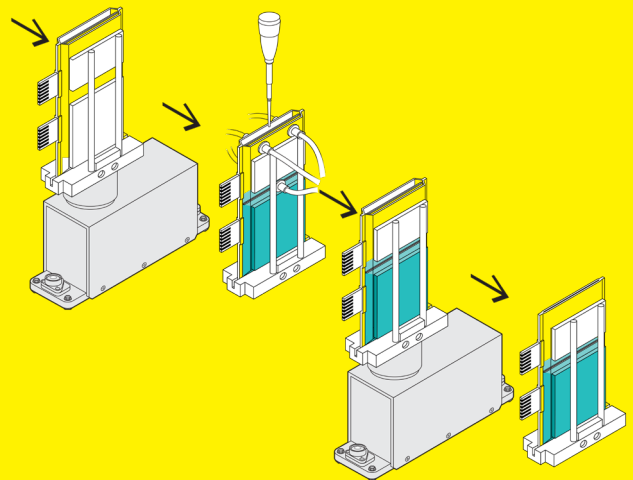
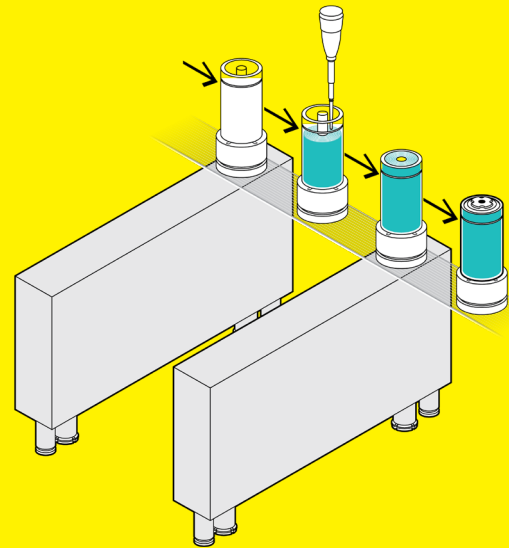
- A wide range of weigh cell constructions and designs enable integration into almost any application.
- Included PC-based software allows for easy configuration and setup of the weigh cells for optimal performance.
- Internal adjustment / calibration are available on many models.
- Additional optional display allows user to monitor weight readings and adjust setup during use.

## Custom Solutions

If the standard weigh cell models do not meet your needs, our Special Solutions Team is on hand to work with your design engineers to ensure the perfect fit.

## Service & Support

Like all high-resolution weighing instrumentation, integrated load cells require periodic maintenance by a trained engineer to ensure continued performance to full specification. Sartorius has a global network of technicians who can support you with calibration, repair or a standard service contract when the need arises. Sartorius will cooperate closely with your engineering team to ensure seamless integration with minimal downtime.



## Precision Weigh Cells

Precision weighing cells cover a range of up to 8.2 kg with a resolution of 0.01-0.001 g. The weighing system is entirely machine-made from an aluminum block. This avoids mechanical and temperature influences created by the inclusion of other materials, ensuring optimal performance.



Model	Readability (g)	Weighing Capacity (g)	Preload (g)	Internal Calibration
WZB553-L	0.001	550	0 - 50	No
WZB653-N	0.001	650	0 - 100	No
WZB1403-N	0.001	1400	0 - 100	No
WZB1403-NC	0.001	1400	0 - 100	Yes
WZB8202-L	0.01	8200	0 - 500	No
WZB8202-N	0.01	8200	0 - 1300	No
WZB8202-NC	0.01	8200	0 - 1300	Yes

## Analytical Weighing Cells

The Analytical weighing cells cover a range of up to 600 g with a resolution of 0.1 mg. The fully automatic production of the weighing system from an aluminium block prevents assembly errors. The weighing system is also available with an internal adjustment function on request. This ensures the unit can be easily and quickly recalibrated for the environment in which it is being used.



Model	Readability (g)	Weighing Capacity (g)	Preload (g)	Internal Calibration
WZB54-L	0.0001	50	min. 5	No
WZB254-L	0.0001	250	0 - 10	No
WZB254-LC	0.0001	250	0 - 10	Yes
WZB254-ND	0.0001	250	0 - 50	No
WZA614-NC	0.0001	610	0 - 90	Yes
WZB254-N	0.0001	250	0 - 50	No
WZB254-NC	0.0001	250	0 - 50	Yes

## Semi Micro | Micro Weighing Cells

The Semi-Micro and Micro cells cover a weighing range of up to 240 g with a resolution of 0.01-0.001 mg. These high-resolution load cells use the same innovative electronics and weighing technology found in our premium laboratory balances. In addition, other weighing systems have been developed for OEM use only. These systems are characterized by increased IP protection and a slim design.



Model	Readability (g)	Weighing Capacity (g)	Preload (g)	Internal Calibration
WZB25-NC	0.00001	20	min. 5	Yes
WZA215-LC	0.00001	210	min. 13.5	Yes
WZA245-NC	0.00001	240	0	Yes
WZA16-LC	0.000001	10	12.5	Yes
WZA26-HC	0.000001	20	12.5	Yes

## Explosion Proof Weighing Cells

Explosion-Proof weighing cells are manufactured specifically for situations where potentially hazardous explosive products are present. These units are designed to meet all relevant safety standards associated with ATEX, IECEx, and related guidelines. They are also certified by FM Approvals for use in Class I, II, III, Division 1 and Class I, Zone 1. They come with specially-protected electronics and weighing modules and include all the necessary cabling and power supply for a safe installation. Refer to the relevant user manuals for details.



Model	Readability (g)	Weighing Capacity (g)	Preload (g)	Internal Calibration
WZA623-NX	0.001	620	600	No
WZA6202-NX	0.01	6200	0	No

**Germany**

Sartorius Lab Instruments  
GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

**USA**

Sartorius Corporation  
North America Inc.  
565 Johnson Avenue  
Bohemia, NY 11716  
Toll-Free +1 800 635 2906

 **For more information, please visit**

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