

## The WZA-N series



### OEM weigh cells

WZA8202-N

#### Outstanding features:

The WZA-N series is the successor of the well established WZA-line. This series focus on weighing resolutions as low as 0.1 mg and capacities of up to 8200 g.

The weigh cell consist of two parts:

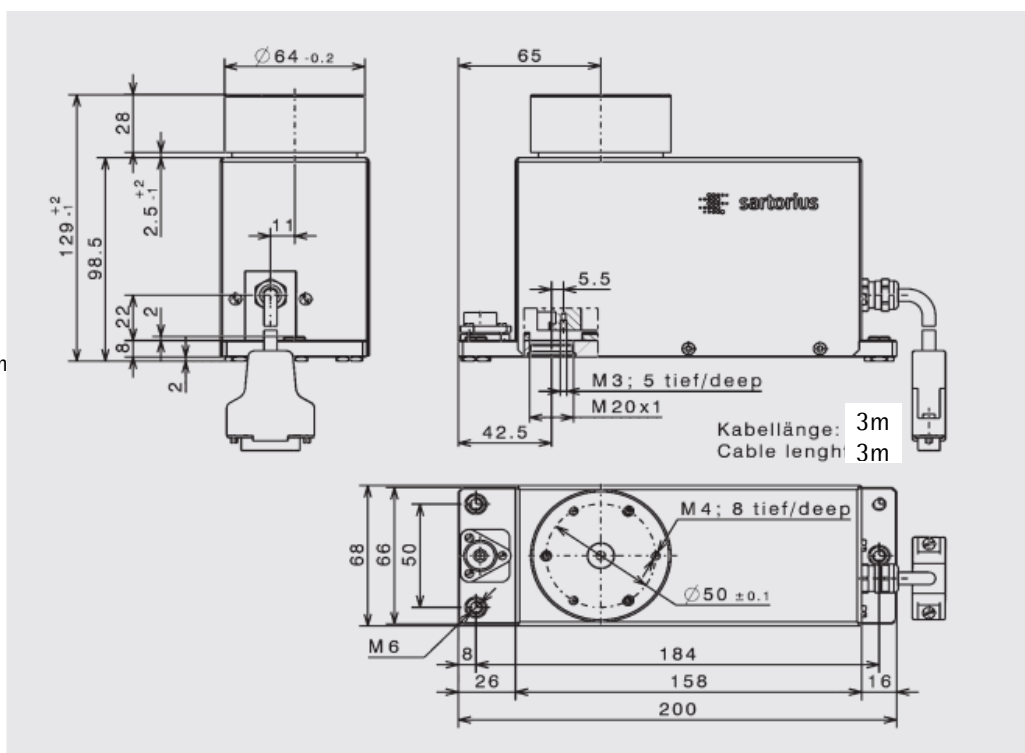
- A monolithic weighing system in stainless steel IP44 housing
- Aluminium electronic box, all connection via the front
- Force compensation technology weighing system with a resolution of 10 mg
- Excellent weighing technology bundled with easy adaptation
- Adjustable load receptor with overload protection, optional lift-off protection
- Possibility for under floor weighing
- Interconnection on electronic front panel
- LED's to control power and data transfer
- Fast warm-up by separation of weighing system and electronics
- RS232C interface for configuration and data transfer
- Optional PC-software or display to handle weighing data and configuration
- Display can be plugged in and out during operation

Model WZA...	8202-N
Capacity	8200 g
Readability	10 mg



## WZA8202-N

Dimensions in mm



### Technical Data

Model	WZA8202-N
Capacity	8200 g
Readability	10 mg
Preload <sup>1)</sup>	1300 g
Repeatability (standard deviation) <sup>2)</sup>	≤ ± 10 mg
Linearity	≤ ± 20 mg
Measuring time <sup>2)3)</sup>	0.6 s / ± 10 mg
Operating temperature range	+10 .... +30 °C
Allowable temperature range	+5 .... +40 °C
Weigh cell dimension	68 x 200 x 129 mm (base x height)
Electronic / PCB dimension	130 x 208 x 45 mm (base x height)
Load receptor (adjustable, with overload protection)	64 mm (diameter), 31 mm height above housing
Under floor load receptor	M3 thread
Cable length weigh cell to electronics	3 m
Power supply	TNG11-2 UNI / 6971993, 115/230 V <sub>AC</sub> , +15 ... - 20%, 48-60 Hz
Power supply (alternative)	12...26, optimal 15V (+15% -10%); 0.5V <sub>PP</sub> (Peak to Peak)
Power consumption	8 VA (average)
Interface	RS232C, software / hardware handshake; No system power via RS232 connector

### Options / Accessories

Windows configuration software for test and adjustment	Sartorius CAS Suite
Liquid-crystal display	YAC01ED

Sartorius Weighing Technology GmbH  
Weender Landstrasse 94-108  
37075 Goettingen, Germany

Phone +49.551.308.0  
Fax : +49.551.308.3502

info.mechatronics@sartorius.com  
www.sartorius-mechatronics.com

- 1) the preload can be configured by using the Sartorius CAS Suite software; higher preload will reduce the weighing range
- 2) depending on measurement equipment set-up and conditions;
- 3) measuring time is the time in which the measured value is in the given range around the static end value; test weight is approx. 25% of the weighing range

Specifications are subject to change without notice.

Current status: July 2011  
Version: 2.0