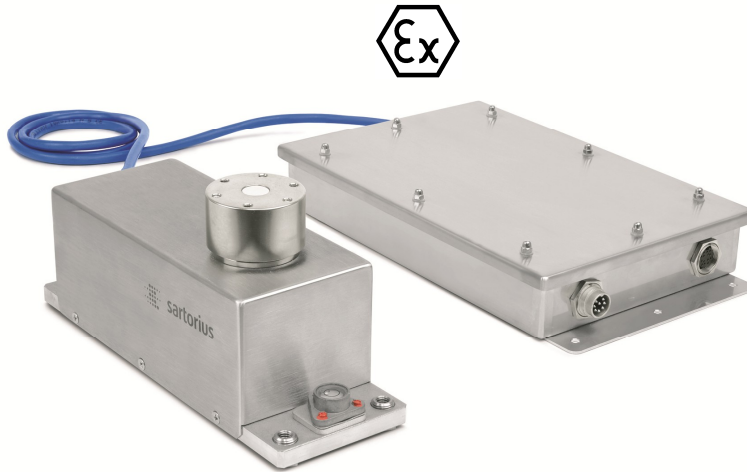




The WZA-EX series



OEM Weigh Cell

Explosion protected

WZA6202-NX
WZA623-NX

ATEX approved

OEM weigh cells

Outstanding features:

Explosion protected weigh cell for use in hazardous areas

- Zone 1 and 2 with gasses of group IIC and temperature class T4
- Zone 21 and 22 for dust atmospheres (Tmax =80°C)



II 2 G Ex ib IIC T4
II 2 D Ex ib IIIB T80 °C
-10°C ≤ Ta ≤ +40°C

FM15ATEX0013X

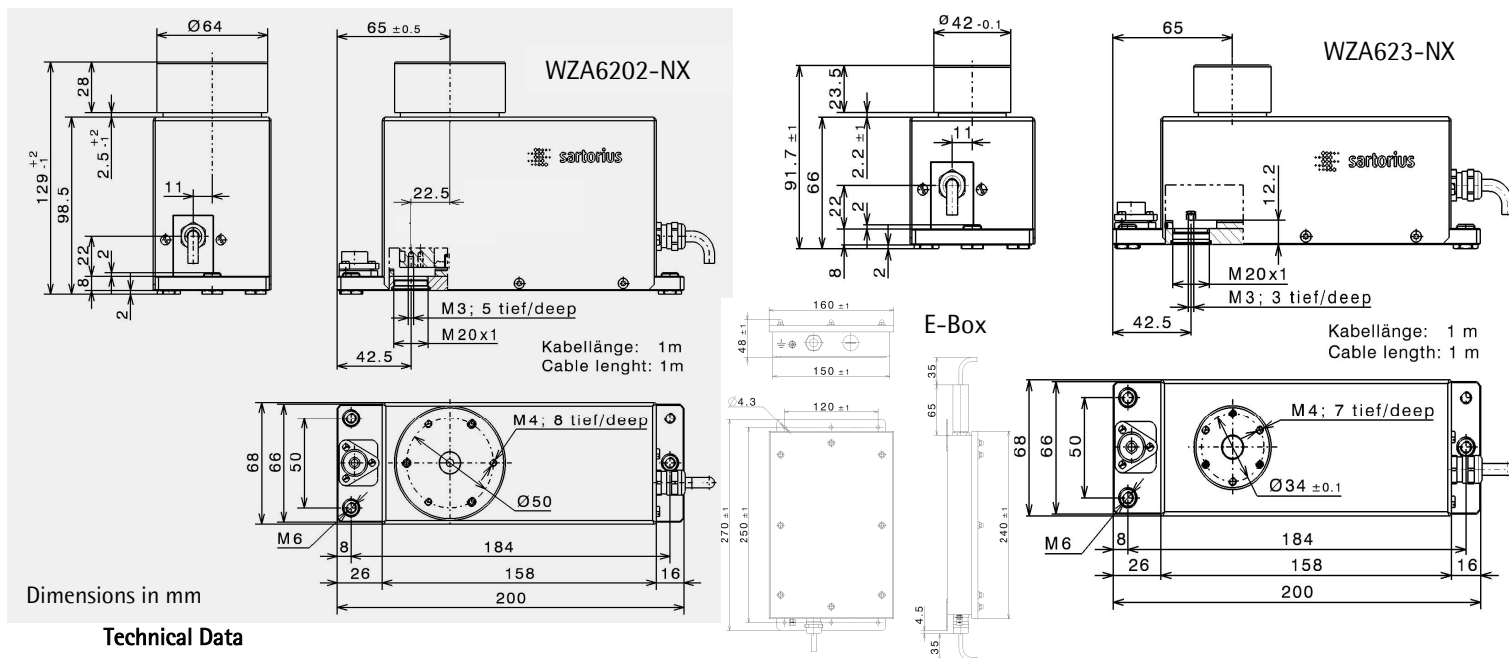
The weigh cell consist of two parts:

- Weigh cell mechanical system in stainless steel IP44 housing
 - Stainless steel electronic box
- Important accessories
- The interface converter YDI05-Z installed in the non hazardous area
 - The power supply
 - Monolithic weighing system with a resolution of 1 mg and 10 mg
 - Excellent weighing technology bundled with easy adaptation
 - Adjustable load receptor with overload protection, for WZA623-NX with optional lift- off protection

- Possibility for under floor weighing
- 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

Model WZA...	6202-NX	623-NX
Capacity	6200 g	620 g
Readability	10 mg	1 mg





Technical Data

Model	WZA6202-NX	WZA623-NX
Capacity	6200 g	620 g
Readability	10 mg	1 mg
Preload, minimal ¹⁾	0 g	600 ± 8 g
Repeatability (standard deviation) ²⁾	≤ ± 10 mg	≤ ± 1 mg
Linearity	≤ ± 20 mg	≤ ± 2 mg
Measuring time ^{2) 3)}	1.5 s / ± 10mg	1.5 s / ± 1 mg
Operating temperature range	+10 +30 °C	
Allowable temperature range	+5 +40 °C	
Weigh cell dimension	68x 200x 129 mm	68 x 200 x 92 mm (base x height)
Electronic / PCB dimension	270 x 160 x 48 mm (base x height)	
Cable length weigh cell to electronics	1 m	
Load receptor (adjustable, with overload protection)	64 mm x 30.5 mm	42 mm x 26 mm (diameter x height)
Under floor load receptor	M3 thread	

Options / Accessories

Power supply unit YPSC01-Z for non-hazardous area	Option N10 and N5; Configurable power supply and communication system for use in non-hazardous areas (USA) 100 – 240 VAC, 40VA, Cable length: AC 3m / DC 20m Option N9 and N5; Configurable power supply and communication system for use in non-hazardous areas (EU) 100 – 240 VAC, 40VA, Cable length: AC 3m / DC 20m
Interface converter in the non-hazardous area YDI05-Z Option M52 and M56	RS232C cable for PC with 9-PIN COM Port, length 3m cable from electronic box to interface converter 20m RS-485 interface available on request
Windows configuration software for test and adjustment	SartoCasSuite public
Display for the hazardous area with 2.5m cable	YAC01NX
Weigh pan to realize lift off protection	Not available
	YWPO1WZA

Sartorius Lab Instruments GmbH&Co.KG
Weender Landstrasse 94-108
37075 Goettingen, Germany

Phone +49.551.308.0
Fax : +49.551.308.3502

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

- 1) the preload can be configured by using the SartoCasSuite 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: May 2015
Version: 1.0