

## The WZA-SE series



### OEM weigh cells

WZA26-HC

#### Outstanding features:

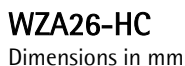
The WZA-HC series is the successor of the well established WZA...CW-line. The products provide a readability of 1 µg for the weighing capacity of 20 g. The weighing system as well the electronic are designed for space sensitive applications in stainless steel housings.

The weigh cell consist of two parts:

- A monolithic weighing system in stainless steel IP44 housing
- Stainless steel electronic box, all connection via the front
- Force compensation technology system with a resolution of 1 µg
- Superb electronics for 24 Mio. step resolution
- Internal calibration weight for easy checks in automated environments
- Load receptor with overload protection
- Possibility for under floor weighing
- Fast warm-up by separation of weighing system and electronics
- USB/ RS232C interface for configuration and data transfer
- Optional PC-software with menu setting, weigh data display and internal calibration functions
- Applications can be found in weighing of samples in automated system, manual pipette calibration and in automated liquid control, e.g. LCD productions.

Model WZA...	26-HC
Capacity	20 g
Readability	1 µg





- 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