



**Susceptometer**  
Innovation in  
mass metrology

# Innovation in mass metrology: A compact solution for testing the magnetic properties of weights

With the introduction of the new OIML Directive R111 comes the recommendation that the magnetic properties of weights be tested. This confirms that the requirements placed on weights no longer only apply to geometric quantities, the density of the material comprising the weight, or its surface qualities. On the contrary, many scientific evaluations support the idea that the susceptibility and magnetization of weights must be considered.

## Susceptometer method

There are several methods for defining the magnetic properties of weights. The so-called susceptometer method is recommended in OIML Directive R111 for testing weights of the accuracy classes E1, E2, F1 and F2. The new YSZ01C | YSZ02C susceptometer from Sartorius enables you to easily and conveniently determine the susceptibility and magnetization of weights in accordance with OIML R111.

In this procedure, the interaction between a permanent magnet and the weight being tested is determined as a force using a high-resolution balance.

Taking into account the known properties of the test magnet, the distance of the weight from the test magnet and the weight geometry, the desired magnetic quantities of the weight can be calculated.

Compared to other well-known equipment, the Sartorius susceptometer stands out on account of its highly compact design. Integration of the individual components, an innovative turning mechanism for adjusting the position of the test magnet and simple adjustment of the height of the loading platform make working with the susceptometer simple and effective. This significantly improves the repeatability of measurements. The susceptometer is available in two different resolution levels. The YSZ01C features 10-µg resolution, and the YSZ02C has 1-µg resolution.

## Application software

The software that comes standard with the susceptometer is very easy to use and enables data to be automatically transferred from the unit. It also supports the user through all required steps and provides assistance with settings before and during a measurement. Menu prompts request all relevant data and guide the user through the steps necessary for performing a measurement. The results of the test are recorded and can be exported to higher-level databases.

## A user-friendly, all-in-one solution

The YSZ01C | YSZ02C susceptometer from Sartorius rounds out the array of accessories available for high-resolution balances and mass comparators. Along with specialized software, the susceptometer provides a unique, complete solution that enables you to ensure that the limits of the magnetic properties of weights are maintained.



YSZ01C | YSZ02C susceptometer

## Technical Specifications

Base dimensions:	338 × 286 mm
Height:	249 mm
Maximum capacity:	50 kg
Dipole moment of the magnet:	m ~ 0.1 Am <sup>2</sup>
Geometry ratio of the magnet:	h/d = 0.87
Height Z <sub>0</sub> (between base of weight and mid-point of the magnet)	Adjustable in fixed steps: Z1=18 Z2=20 Z3=27 Z4=35 Z5=43 mm
Field strength due to different height Z <sub>0</sub> :	2,700; 2,000; 800; 360; 200 A/m
Readability of the susceptometer:	YSZ01C: 10 mg YSZ02C: 1 mg
Turning mechanism for magnet:	Easy to operate by external rotary knob Marking for N-S orientation of the magnet
Software:	Convenient application software, default parameters and easy, user-defined configurations possible Quick check function Printing of reports, exporting of results
Data transfer protocol:	HTML mode and data transfer to user-specific metrology software via CVS file format
Applications for the Sartorius susceptometer:	OIML R111-compliant testing of weights; please note the tables on the next page for the ranges of susceptibility and magnetization measurements

# Data sheet for YSZ01C | YSZ02C susceptometers

## Susceptibility measurements

Maximum susceptibility  $\chi$  according to OIML R111 Recommendation

Change in susceptibility  $\chi$  for  $Z_0 = 18$  mm for 1 digit  $\Delta\chi$

Measuring range of the Sartorius susceptometer\*

Nominal Mass	OIML Class				Model Order No.: YSZ01C		YSZ02C		Model Order No.: YSZ01C		YSZ02C	
	E1	E2	F1	F2	Readability:	10 $\mu\text{g}$	Readability:	1 $\mu\text{g}$	Readability:	10 $\mu\text{g}$	Readability:	1 $\mu\text{g}$
2 g	0.06	0.18			2 g	0.00084	0.000084		2 g	0.042... 1	0.0042... 1	
5 g	0.06	0.18			5 g	0.00047	0.000047		5 g	0.024... 1	0.0024... 1	
10 g	0.06	0.18			10 g	0.00031	0.000031		10 g	0.016... 1	0.0016... 1	
20 g	0.02	0.07	0.2	0.8	20 g	0.0002	0.00002		20 g	0.010... 1	0.0010... 1	
50 g	0.02	0.07	0.2	0.8	50 g	0.00012	0.000012		50 g	0.006... 1	0.0006... 1	
100 g	0.02	0.07	0.2	0.8	100 g	0.0001	0.00001		100 g	0.005... 1	0.0005... 1	
200 g	0.02	0.07	0.2	0.8	200 g	0.00008	0.000008		200 g	0.004... 1	0.0004... 1	
500 g	0.02	0.07	0.2	0.8	500 g	0.00006	0.000006		500 g	0.003... 1	0.0003... 1	
1 kg	0.02	0.07	0.2	0.8	1 kg	0.00006	0.000006		1 kg	0.003... 1	0.0003... 1	
2 kg	0.02	0.07	0.2	0.8	2 kg	0.00006	0.000006		2 kg	0.003... 1	0.0003... 1	
5 kg	0.02	0.07	0.2	0.8	5 kg	0.00006	0.000006		5 kg	0.003... 1	0.0003... 1	
10 kg	0.02	0.07	0.2	0.8	10 kg	0.00006	0.000006		10 kg	0.003... 1	0.0003... 1	
20 kg	0.02	0.07	0.2	0.8	20 kg	0.00006	0.000006		20 kg	0.003... 1	0.0003... 1	
50 kg	0.02	0.07	0.2	0.8	50 kg	0.00005	0.000005		50 kg	0.003... 1	0.0003... 1	

\*) The quantity contributed by the display resolution to the total uncertainty of measurement is  $\leq 2\%$  within the range indicated

## Magnetization measurements

Maximum magnetization  $\mu_0 M_z$  according to the OIML R111 Recommendation

Change in magnetization  $\Delta\chi$  for  $Z_0 = 18$  mm for 1 digit  $\Delta\mu_0 M_z$

Measuring range of the Sartorius susceptometer\*\*

Nominal Mass	OIML Class				Model Order No.: YSZ01C		YSZ02C		Model Order No.: YSZ01C		YSZ02C	
	E1	E2	F1	F2	Readability:	10 $\mu\text{g}$	Readability:	1 $\mu\text{g}$	Readability:	10 $\mu\text{g}$	Readability:	1 $\mu\text{g}$
2 g	2.5 $\mu\text{T}$	8 $\mu\text{T}$			2 g	1.8 $\mu\text{T}$	0.18 $\mu\text{T}$		2 g	18 $\mu\text{T}... 1$ mT	1.8 $\mu\text{T}... 1$ mT	
5 g	2.5 $\mu\text{T}$	8 $\mu\text{T}$			5 g	0.9 $\mu\text{T}$	0.09 $\mu\text{T}$		5 g	9 $\mu\text{T}... 1$ mT	0.9 $\mu\text{T}... 1$ mT	
10 g	2.5 $\mu\text{T}$	8 $\mu\text{T}$			10 g	0.6 $\mu\text{T}$	0.06 $\mu\text{T}$		10 g	6 $\mu\text{T}... 1$ mT	0.6 $\mu\text{T}... 1$ mT	
20 g	2.5 $\mu\text{T}$	8 $\mu\text{T}$			20 g	0.4 $\mu\text{T}$	0.04 $\mu\text{T}$		20 g	4 $\mu\text{T}... 1$ mT	0.4 $\mu\text{T}... 1$ mT	
50 g	2.5 $\mu\text{T}$	8 $\mu\text{T}$			50 g	0.2 $\mu\text{T}$	0.02 $\mu\text{T}$		50 g	2 $\mu\text{T}... 1$ mT	0.2 $\mu\text{T}... 1$ mT	
100 g	2.5 $\mu\text{T}$	8 $\mu\text{T}$	25 $\mu\text{T}$	80 $\mu\text{T}$	100 g	0.2 $\mu\text{T}$	0.01 $\mu\text{T}$		100 g	2 $\mu\text{T}... 1$ mT	0.2 $\mu\text{T}... 1$ mT	
200 g	2.5 $\mu\text{T}$	8 $\mu\text{T}$	25 $\mu\text{T}$	80 $\mu\text{T}$	200 g	0.1 $\mu\text{T}$	0.01 $\mu\text{T}$		200 g	1 $\mu\text{T}... 1$ mT	0.1 $\mu\text{T}... 1$ mT	
500 g	2.5 $\mu\text{T}$	8 $\mu\text{T}$	25 $\mu\text{T}$	80 $\mu\text{T}$	500 g	0.1 $\mu\text{T}$	0.01 $\mu\text{T}$		500 g	1 $\mu\text{T}... 1$ mT	0.1 $\mu\text{T}... 1$ mT	
1 kg	2.5 $\mu\text{T}$	8 $\mu\text{T}$	25 $\mu\text{T}$	80 $\mu\text{T}$	1 kg	0.1 $\mu\text{T}$	0.01 $\mu\text{T}$		1 kg	1 $\mu\text{T}... 1$ mT	0.1 $\mu\text{T}... 1$ mT	
2 kg	2.5 $\mu\text{T}$	8 $\mu\text{T}$	25 $\mu\text{T}$	80 $\mu\text{T}$	2 kg	0.1 $\mu\text{T}$	0.01 $\mu\text{T}$		2 kg	1 $\mu\text{T}... 1$ mT	0.1 $\mu\text{T}... 1$ mT	
5 kg	2.5 $\mu\text{T}$	8 $\mu\text{T}$	25 $\mu\text{T}$	80 $\mu\text{T}$	5 kg	0.1 $\mu\text{T}$	0.01 $\mu\text{T}$		5 kg	1 $\mu\text{T}... 1$ mT	0.1 $\mu\text{T}... 1$ mT	
10 kg	2.5 $\mu\text{T}$	8 $\mu\text{T}$	25 $\mu\text{T}$	80 $\mu\text{T}$	10 kg	0.1 $\mu\text{T}$	0.01 $\mu\text{T}$		10 kg	1 $\mu\text{T}... 1$ mT	0.1 $\mu\text{T}... 1$ mT	
20 kg	2.5 $\mu\text{T}$	8 $\mu\text{T}$	25 $\mu\text{T}$	80 $\mu\text{T}$	20 kg	0.1 $\mu\text{T}$	0.01 $\mu\text{T}$		20 kg	1 $\mu\text{T}... 1$ mT	0.1 $\mu\text{T}... 1$ mT	
50 kg	2.5 $\mu\text{T}$	8 $\mu\text{T}$	25 $\mu\text{T}$	80 $\mu\text{T}$	50 kg	0.2 $\mu\text{T}$	0.02 $\mu\text{T}$		50 kg	2 $\mu\text{T}... 1$ mT	0.2 $\mu\text{T}... 1$ mT	

\*\*) The quantity contributed by the display resolution to the total uncertainty of measurement is  $\leq 10\%$  within the range indicated



YSZ01C | YSZ02C software interface

Sartorius AG  
Weender Landstrasse 94–108  
37075 Goettingen, Germany

Phone +49.551.308.0  
Fax +49.551.308.3289

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