

Sartorius Susceptometer Model YSZ01C | YSZ02C

Susceptometer



Special performance features:

The susceptometer measures the susceptibility and magnetization of weights of accuracy classes E1, E2, F1 and F2.

With a convenient turning mechanism for the internal neodymium-boron-iron magnet with a dipole moment of approx. 0.1 Am²; distances Z0 between the center points of the weight and magnet):

Z0 = distance,	Class,	Field strength
Z1 = 18 mm,	E1,	2700 A/m
Z2 = 20 mm,	E1,	2000 A/m
Z3 = 27 mm,	E2,	800 A/m
Z4 = 35 mm,	E2,	360 A/m
Z5 = 43 mm,	F1,F2,	200 A/m

Model	YSZ01C	YSZ02C
Maximum capacity	50 kg	50 kg
Application range	2 g - 50 kg	2 g - 50 kg
OIML R111	E2 F1 F2 M1	E1 E2 F1 F2 M1
Readability	10 µg	1 µg

*) The typical standard deviation "s" is the repeatability, calculated from 6 ABBA-cycles, after elimination of the drift. Prerequisites for the standard deviation are good climatic conditions in accordance with OIML R111 for an E1 mass laboratory.

Technical Specifications

Model	YSZ01C	YSZ02C
Maximum capacity	50 kg	50 kg
Application range	2 g – 50 kg	2 g – 50 kg
Readability	10 µg	1 µg
Repeatability, s*	20 µg	8 µg
Repeatability (typical), s*	10 µg	5 µg
Electronic weighing range & taring range	10 g	10 g
Linearity	0.1 mg	0.05 mg
Range sensitivity	20 µg/1 g	10 µg/1 g
Stabilization time	15 s	15 s

Basic Equipment

Interfaces	RS232	RS232
Draft shield	x	x
PC software	x	x
Test certificate	Sartorius	Sartorius
PC connecting cable	x	x

Ambient Conditions

Permissible operating temperature range	17-27 °C	17-27 °C
Recommended operating temperature	22 °C	22 °C
Temperature fluctuation	0.3°C/h 0.5°C/12h	0.3°C/h 0.5°C/12h
Max. air movement	< 0.2 m/s	< 0.2 m/s
Humidity range	40-60 %	40-60 %
Humidity fluctuation	5%/4h	5%/4h
Power supply	100-240V AC/50-60Hz	100-240V AC/50-60Hz
Power consumption	< 35 VA	< 35 VA

*] The typical standard deviation "s" is the repeatability, calculated from 6 ABBA-cycles, after elimination of the drift. Prerequisites for the standard deviation are good climatic conditions in accordance with OIML R111 for an E1 mass laboratory.

Dimensions

Sample size(DxH)	280x300 mm	280x300 mm
External dimensions (WxDxH)	338x286x249 mm	338x286x249 mm
Net weight	20 kg	20 kg
Gross weight	26.4 kg	26.4 kg
Number of packages	1	1
Packing data	59x47x75 cm	59x47x75 cm

Applications

OIML R111	E2 F1 F2 M1	E1 E2 F1 F2 M1
-----------	-------------	----------------

Optional Accessories

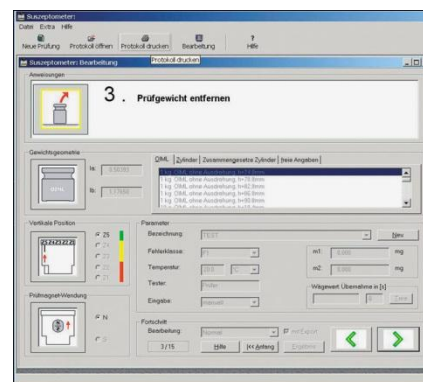
External calibration weight	10 g E2 YCW412-02	10 g E2 YCW412-02
Weighing table	YWT03	YWT03
Crane with chain hoist	YLD01C	YLD01C
Gripper for weights with handle	YLD02C	YLD02C
Lifting device for 10 kg	YAW51	YAW51
Lifting device for 20 kg	YAW52	YAW52
Lifting device for 50 kg	YAW53	YAW53
1 kg PTB susceptibility reference	YSZ01RSC	YSZ01RSC
Magnet calibration set for susceptometer	YSZ01RMC	YSZ01RMC



YSZ01RMC



YSZ01RSC



YSZ01C | YSZ02C

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

Phone +49.551.308.0
Fax +49.551.308.3289

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



Specifications subject to change without notice.
Version: 01|2013