

Sartocheck® mini

Filter Integrity Tester for Food & Beverage Applications

Product Information

The automatic filter integrity tester Sartocheck® mini can be used to verify the integrity of membrane filters which are used in the food & beverage environment.



Description

Taken those specific needs into account, this unit offers the following main features:

- Automatic filter integrity tester
- Pressure Drop Test &
- Diffusion Test
- Small, portable unit
- 19 different test programs
- 200 test results can be stored
- LCD display
- Automatic venting after the test
- Thermo-printer (57 mm paper)
- Easy and reliable data transfer to PC
- High capacity batteries for up to 4 hrs work
- For filter housings with a net volume between 10 mL and 999 L
- Ingress protection IP54
- Incl. bag and case

Performance

Test Result Documentation

Test results are automatically printed using the built-in thermo printer. An additional port allows the connection of an external printer.

Data Storage

The unit stores up to 200 test results in the internal memory. To avoid the oldest data to be overwritten, electronic data can be stored on a connected PC with user-friendly software. The same software can be used for programming the device.

Technical Data

Specifications

Power requirements	100 – 240 V AC, 50 60 Hz
Max. Power Input	20 W
Max. inlet pressure	4500 mbar
Dimensions	$315 \times 150 \times 280 \text{ mm}$
Weight	ca. 3.900 g
Languages	German, English, French, Italian, Spanish, Portuguese

Operating Conditions

Temperature	0 – 40 °C
Humidity	5 – 95% rel.

Measuring Ranges

Test pressure	50 – 4000 mbar
Max inlet pressure	4500 mbar
Over pressure protection	8000 mbar
Net volume	10 mL-999 L

Measuring Accuracy

Rel. deviation pressure	< 0.2%
measurement	
Abs. deviation pressure	max. ±4 mbar (@20°C)
measurement	

Interfaces

Communication port	Mini USB	
--------------------	----------	--

Equipment Supplied

- Sartocheck® mini integrity test unit
- Printer paper (4 rolls)
- Pressure inlet tubing (18104)
- Pressure outlet tubing (18103)
- Carrier bag (soft case)
- Hard case

Order Information

26292---02

Sartorius Stedim Biotech GmbH August-Spindler-Strasse 11 37079 Goettingen, Germany Phone +49.551.308.0 Fax +49.551.308.3289 www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178 Argentina +54.11.4721.0505 Brazil +55.11.4451.6226 Mexico +52.5555.62.1102 UK +44.1372.737159 France +33.442.845600 Italy +39.055.63.40.41 Spain +34.913.586.098 Russian Federation +7.812.327.5.327 Japan +81.3.4331.4300 China +86.21.68782300

Specifications subject to change without notice. Copyright Sartorius Stedim Biotech GmbH. Printed in the EU on paper bleached without chlorine. Publication No.: SPI2015-e160702 Order No.: 85032-535-23 Ver. 07 | 2016