



# Sartocheck® mini

Filter Integrity Tester for Food & Beverage Applications



## Technical Specifications

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

## Operating Conditions

Temperature	3–30°C.
Humidity	5–95% rel.

## Measuring Ranges

Test pressure	0–3900 mbar
Max. inlet pressure	0–4500 mbar
Net volume	0.1–999 L

## Measuring Accuracy

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

## Interfaces

External printer	Centronics 25 pol
Communication port	232, 9 pole male

## Equipment Supplied

- Sartocheck® mini integrity test unit
- Low volume adapter for net volumes <5 L
- Printer paper (4 rolls)
- Pressure inlet tubing (18104)
- Pressure outlet tubing (18103)
- Carrier bag (soft case)
- Hard case

## Order Information

26292---01

## Description

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.

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
- 100 test results to 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 0.1 and 999 L
- Protection rating IP50
- Incl. bag and case

## 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 100 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.

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](http://www.sartorius-stedim.com)

USA Toll-Free +1.800.368.7178

UK +44.1372.737159

France +33.442.845600

Italy +39.055.63.40.41

Spain +34.91.3586102

Japan +81.3.3740.5407

Specifications subject to change  
without notice. Printed and copyrighted  
by Sartorius Stedim Biotech GmbH

W · G

Publication No.: SPI2015-e09061

Order No.: 85032-535-23

Ver. 06 | 2009