

# Sartocheck<sup>®</sup> mini

## Filter Integrity Tester for Food & Beverage Applications



### Product Information

The automatic filter integrity tester Sartocheck<sup>®</sup> 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.

## Technical Data

### 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	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

## 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.

### Measuring Accuracy

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

### 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

#### Germany

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

#### USA

Sartorius Stedim North America Inc.  
565 Johnson Avenue  
Bohemia, NY 11716  
Toll-Free +1 800 368 7178

 For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)