



Sartocheck® Cleaning Kit

Reliable cleaning of the internal pneumatics



The Cleaning Kit

The Cleaning Kit is a set of components that are necessary to flush the unit with suitable cleaning liquid and to dry the system afterwards. The following parts are included:

- 50 L modified plastic bottle (drain)
- T-piece connector with shut-off valve
- Pressure reducer
- Set of 5 tubings (PU) with quick connectors

The ends of the tubings are clearly labeled to ensure proper hookup. Together with the on-screen instructions, the installation and the process of cleaning & drying is easy to follow.

In addition to the Cleaning Kit, a stainless-steel pressure vessel is needed for the filling procedure.

Recommended Cleaning Fluids

We recommend the usage of one of the following fluids for cleaning:

- Water (cold, warm or hot up to 80°C)
- NaOH solutions up to 1 M (0.1 M recommended)

Duration of Cleaning & Drying

The entire cleaning program (that includes steps for cleaning, rinsing and drying) only takes approximately 1.5 hours in total. This is a fraction of the down time that would otherwise be incurred for factory cleaning. The Sartocheck's® automated cleaning program provides a huge advantage from a cost of ownership perspective over alternative integrity testers.

Efficacy of Cleaning & Drying

The efficacy of the cleaning procedure and the validity of the timing sequence was proven in an elaborated test series.

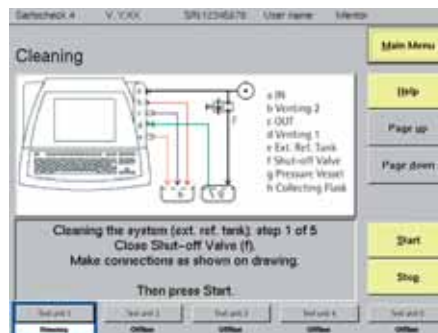
Use this unique feature of Sartocheck® to make your process of filter integrity testing as safe as possible.

Description

Sartocheck® offers the outstanding feature of an embedded automated cleaning cycle program which allows the reliable cleaning of the complete internal pneumatics even with aggressive cleaning liquids.

During the routine process of filter integrity testing, the possibility of cross contamination exists. Therefore, it is critical to have the ability to reliably clean the internal pneumatics of such an integrity tester. Having this function automated avoids the high cost and downtime associated with factory service calls.

In order to allow reliable cleaning, only materials of highest quality and broadest chemical resistance (stainless steel, PTFE) are used for those components which have direct contact with the air used for pressurization. This unique feature allows the use of even aggressive liquids (up to 1 M NaOH) for the internal cleaning of Sartocheck® integrity testers.



Patented Cleaning & Drying Procedure

The cleaning process of the Sartocheck® is fully automated and software guided. The cleaning procedure is divided into multiple simple interactive steps. Intuitive on-screen instructions clearly indicate the required user actions for each step until the entire cycle is complete.

After the successive cleaning and rinsing steps are finished, the Sartocheck® automatically engages the drying program to remove any existing water residue.

The Drying program can be started separately in order to allow a potential particle and moisture removal by using compressed air.



Order Information

Cleaning Kit complete	26288--CK
Pressure Vessel (if required)	26288--PV

Individual Components of 26288--CK

Pressure manifold	26288--PD
Tubing "Out"	18103
Tubing "Venting 1"	18116
Tubing Venting 2	18117
Tubing "Ext. Ref. Tank"	18118
Plastic bottle (50 L) (modified)	26288--CF

Which Units can be cleaned?

The Cleaning Kit can be used for the following integrity testers:

- Sartocheck® 3 plus
- Sartocheck® 4 (Version 2.03)
- Sartocheck® 4 plus (Version 1.0)
- Multiunit

The equipment of the Cleaning Kit was adapted to the latest version of the Sartocheck® software. In order to use the kit for older versions, a separate instruction has to be used. Therefore, we strongly recommend to upgrade the units before using the new Cleaning Kit.

Service

As an alternative to the end user performing the automated cleaning procedure, Sartorius Stedim Biotech has a world-class Service Team of highly trained professionals that are available to perform the cleaning of your Sartocheck®, perhaps during the annual maintenance & calibration work. Please contact our Service Team for further information.

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
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.90.2110935
Japan +81.3.3740.5407

Specifications subject to change without notice. Printed and copyrighted by Sartorius Stedim Biotech GmbH
W · G
Publication No.: SPI2017-e10031
Order No.: 85032-536-83
Ver. 03 | 2010