



Manual Filtration Skid – Essential Range

Filtration Systems for Wines,
Flavoured wines, Sparkling Wines

Product Information

The manual filtration skid of the Essential range is designed for the clarification and microorganisms retention before bottling. It uses disposable filter cartridges and can also be connected to a Clean In Place (CIP) system. Two sizes are available:

- up to 2000 Bottles/hour (BtIs/hr)
- up to 5000 BtIs/hr



Applications

The manual filtration skid Essential is designed for the clarification, filtration and microorganisms retention of wines, flavoured wines or sparkling wines.

Process Optimization

No wine oxidation with a completely closed filtration system.
Filtration for 1, 2 or 3 stages.

Repeatability

Reliable bottling and optimization of wine production with integrity checking of final membrane filter

Savings

Longer cartridge life time with individual cleaning of each filtration stage. Limited water loss during sterilization step with recovery circuit.

Efficiency and safety

- Total process control with sensors such as manometers and thermometers.
- Water leak valve and fully drainable skids by gas ensure no water comes into contact with wine bottling processes.
- Wine sampling valve to check the wine quality & filtration performances before bottling.

Filtration Skids Description

An essential filtration range consists of two or three housings dedicated to wine filtration and one housing dedicated to the filtration of cleaning water. Each housing is connected to the others by a ramp and can either be used or isolated by the manual inlet and outlet valves. The system is equipped with temperature and pressure sensors, drain and sampling valves.

Manual Filtration Skid Essential Range

	Essential 2.2S	Essential 2.2	Essential 2.3	Essential 5.2S	Essential 5.2	Essential 5.3
Filtration	1 – 2 stages	1 – 2 stages	1, 2 or 3 stages	1 – 2 stages	1 – 2 stages	1, 2 or 3 stages
Flow rate	2 000 btls/h	2 000 btls/h	2 000 btls/h	5 000 btls/h	5 000 btls/h	5 000 btls/h
Housing Cartridge	2 housings 1 × 30" cartridge per housing	2 housings 1 × 30" cartridge per housing	3 housings 1 × 30" cartridge per housing	2 housings 3 × 30" cartridges per housing	2 housings 3 × 30" cartridges per housing	3 housings 3 × 30" cartridges per housing
Pre-filters backwash	Included	Option	Option	Included	Option	Option
High pressure equipment for sparkling wine	Option	Option	Option	Option	Option	Option
Built on wheels	Option	Option	Option	Option	Option	Option
Automated sterilization processes	Option	Option	Option	Option	Option	Option
Integrity testing of final filter	Option	Option	Option	Option	Option	Option

Options

Additional options such as automated dosing of oenological adjuvants can be added when the skid is connected to a Sartorius CIP system.

CIP System

All Manual filtration skid of the Essential range can be connected to a CIP system. The CIP System produces and maintains hot water up to 90°C | 194°F. It supplies filtered water for the filling line (filtration skid, piping and bottling machine) either for cleaning purposes at 20°C | 68°F and 50°C | 122°F or for sterilization at 90°C | 194°F.

Technical Data

- Main construction materials: 304L Stainless Steel
- Housings material: 316L Stainless Steel
- Manual valves gasket material: EPDM
- Surface finish: electropolished
- Inlet and outlet connection: DIN standard
- Mobility (option): four swivelling wheels
- Food and CE compliant
- Dimensions and weight (approx.)

	W × D × H	Weight
Essential 2.2S	1705 × 465 × 1770 mm (67 1/8 × 18 5/16 × 69 11/16 in)	90 kg (198 lb)
Essential 2.2	1705 × 465 × 1770 mm (67 1/8 × 18 5/16 × 69 11/16 in)	90 kg (198 lb)
Essential 2.3	2190 × 465 × 1770 mm (86 1/4 × 18 5/16 × 69 11/16 in)	101 kg (222 lb)
Essential 5.2S	1660 × 595 × 1685 mm (65 3/8 × 18 5/16 × 69 11/16 in)	132 kg (291 lb)
Essential 5.2	1660 × 595 × 1685 mm (65 3/8 × 23 7/16 × 66 5/16 in)	132 kg (291 lb)
Essential 5.3	2050 × 595 × 1685 mm (80 11/16 × 23 7/16 × 66 5/16 in)	167 kg (368 lb)

Filter Cartridges

- Prefiltration step with Sartopure® IND: 0.45 µm; 0.65 µm; 1.2 µm; 3 µm; 5 µm; 8 µm; 20 µm and 50 µm
- Final filtration for yeast and bacterial retention with either 1 µm; 0.65 µm Sartobev®, 0.65 µm; 0.45 µm Sartocool® or 0.65 µm; 0.45 µm Vinosart® PS.

Ordering Information

Product code	Designation
Manual Filtration skid Essential range	
FSKID-ESS-2.2S	Manual filtration skid essential up to 2000 btls/h 2 stages of filtration – Basic version
FSKID-ESS-2.2	Manual filtration skid essential up to 2000 btls/h 2 stages of filtration
FSKID-ESS-2.3	Manual filtration skid essential up to 2000 btls/h 3 stages of filtration
FSKID-ESS-5.2S	Manual filtration skid essential up to 5000 btls/h 2 stages of filtration – Basic version
FSKID-ESS-5.2	Manual filtration skid essential up to 5000 btls/h 2 stages of filtration
FSKID-ESS-5.3	Manual filtration skid up to 5000 btls/h 3 stages of filtration
Manual Filtration skid Essential range for sparkling wine	
FSKID-ESS-2.2SSW	Manual filtration skid essential up to 2000 btls/h 2 stages of filtration for sparkling wine – Basic version
FSKID-ESS-2.2-SW	Manual filtration skid essential up to 2000 btls/h 2 stages of filtration for sparkling wine
FSKID-ESS-2.3-SW	Manual filtration skid essential up to 2000 btls/h 3 stages of filtration for sparkling wine
FSKID-ESS-5.2S-SW	Manual filtration skid essential up to 5000 btls/h 2 stages of filtration for sparkling wine – Basic version
FSKID-ESS-5.2-SW	Manual filtration skid essential up to 5000 btls/h 2 stages of filtration for sparkling wine
FSKID-ESS-5.3-SW	Manual filtration skid essential up to 5000 btls/h 3 stages of filtration for sparkling wine
Manual Filtration skid Essential range - options list	
FSKID-ESS-OPT1	Pre-filters backwash
FSKID-ESS-OPT2	Built on wheels
FSKID-ESS-OPT3	Automated sterilization program
FSKID-ESS-OPT4	Automated sterilization program + automated integrity test
Clean In Place system Essential range	
CIP-ESS-1000	Manual clean in place system 1000L
CIP-ESS-1500	Manual clean in place system 1500L
CIP-ESS-2000	Manual clean in place system 2000L
Clean In Place system Essential range - options list	
FSKID-ESS-CIP-OPT1	Automated pump start
FSKID-ESS-CIP-OPT2	Automated sterilization program
FSKID-ESS-CIP-OPT3	Automated sterilization program + water recovery
FSKID-ESS-CIP-OPT4	Automated sterilization program + automated integrity test
FSKID-ESS-CIP-OPT5	Automated sterilization program+ oenologic product dosage

Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49.551.308.0

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5-7
34302 Guxhagen
Phone +49.5665.407.0

France

Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques - CS 91051
13781 Aubagne Cedex
Phone +33.442.845600

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques - CS 71058
13781 Aubagne Cedex
Phone +33.442.845600

Austria

Sartorius Stedim Austria GmbH
Modecenterstrasse 22
1030 Vienna
Phone +43.1.7965763.18

Belgium

Sartorius Stedim Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles
Phone +32.2.756.06.80

Hungary

Sartorius Stedim Hungária Kft.
Kagyló u. 5
2092 Budakeszi
Phone +36.23.457.227

Italy

Sartorius Stedim Italy S.r.l.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)
Phone +39.055.63.40.41

Netherlands

Sartorius Stedim Netherlands B.V.
Phone +31.30.60.25.080
filtratie.nederland@sartorius-stedim.com

Poland

Sartorius Stedim Poland Sp. z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.647.38.40

Russian Federation

LLC "Sartorius Stedim RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg
Phone +7.812.327.53.27

Spain

Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)
Phone +34.913.586.098

Switzerland

Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen
Phone +41.52.354.36.36

U.K.

Sartorius Stedim UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9 QQ
Phone +44.1372.737159

Ukraine

LLC "Sartorius Stedim RUS"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918

Americas

USA

Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716
Toll-Free +1.800.368.7178

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil
Phone +55.11.4362.8900

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepotzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepotzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
leadsmex@sartorius.com

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800

China

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China
Phone +86.21.6878.2300

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Beijing Branch Office
No. 33 Yu'an Road
Airport Industrial Park Zone B
Shunyi District, Beijing 101300
Phone +86.10.8042.6501

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
Guangzhou Branch Office
Room 1105
Xing Guang Ying Jing Building
No. 119, Shui Yin Road
Yue Xiu District, Guangzhou 510075
Phone +86.20.3836.4193

India

Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250

Japan

Sartorius Stedim Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan
Phone +81.3.4331.4300

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia
Phone +60.3.8996.0622

Singapore

Sartorius Stedim Singapore Pte. Ltd.
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684
Phone +65.6872.3966

South Korea

Sartorius Korea Biotech Co., Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, BunDang-Gu
SeongNam-Si, GyeongGi-Do, 463-400
Phone +82.31.622.5700



▶ www.sartorius-stedim.com