



TuFlux[®] SIL

Sartorius Stedim Biotech Silicone Tubing

Product Information

Sartorius Stedim Biotech TuFlux[®] SIL is designed to facilitate fluid transfer in the biopharmaceutical manufacturing processes. TuFlux[®] SIL is a highly resistant platinum cured silicone tubing manufactured by Raumedic. TuFlux[®] SIL is available in 7 different dimensions from 1/8" (3.2 mm) to 3/4" (19.1 mm) for the internal dimensions and a wall thickness from 1.6 to 4.8 mm.



Applications

TuFlux[®] SIL is designed to be used in many pharmaceutical and biopharmaceutical applications such as:

- Media and Buffer processing
- Filtration
- Fermentation
- Cell harvest

Flexibility

Sartorius Stedim Biotech silicone tubing TuFlux[®] SIL is available as non sterile tubing coils and can also be mounted and sterilized on any Sartorius Stedim Biotech single-use assembly. Several dimensions are available to match the process requirements in term of flow rate and to be perfectly adapted to the peristaltic pump for maximum pumping efficiency.

Features & Benefits

- Low extractable profile and unique Validation Guide for TuFlux[®] SIL available: Ease the validation of TuFlux[®] SIL in a process.
- Shore hardness A 60: Very good rupture strength in peristaltic pumps and reduced kink effect.
- Platinum Cured Silicone: Resistant against weak acids and bases and extraordinary heat and cold resistance.
- Translucent: Visual contact with the fluid.
- Printing of ID and OD dimensions on the tubing: Facilitates tubing identification.
- Coils wrapped in double PE-bags: Tubing protected and easy to introduce in clean room.
- "Low-Tack" significantly reduced surface friction: Easier handling with gloves.
- Certificate of Release: Delivered with each order for traceability

Technical Data

Specifications

Colour	Transparent
Material	Platinum cured silicone suitable for pumping application
Shore hardness A	60 ± 5
Temperature range	from -60°C to +200°C (-76°F to +392°F)
Tear strength	> 8.0 MPa
Elongation at break	> 500%
Compliant	E.P. 3.1.9, USP <88> Class VI, ADCF
Printing	Physiologically inert ink showing tubing dimensions
Sterilization	Gamma irradiation and autoclave
Tubing coils packaging	Primary and secondary PE packaging

Tubing Coil

TuFlux® SIL is available from stock as non sterile tubing coil packaged in a double PE bag. The length of a coil is depending on the tubing dimensions and varies between 100 m to 15 m.

TuFlux® Sil is compatible with Gamma Sterilization and autoclave. The Shelf life is 5 years.

Integration

- All TuFlux® SIL tubing dimensions can be integrated into any Sartorius Stedim Biotech assembly.
- TuFlux® SIL could be pre-assembled to bags, filters and then Gamma sterilized to offer a ready-to-use solution.
- TuFlux® SIL can also be assembled on tubing and filtration sets without Gamma sterilisation to offer ready-to-autoclave solutions to end users when autoclaving is required in the process.

Security of Supply

Sartorius Stedim Biotech has established multiple manufacturing sites with consistent industrial processes for its fully integrated disposable assemblies. The expertise of designing Single-Use solutions based on collaborative supplier management and customer demand planning assures a state of the art and robust supply chain that can cope with strong market growth.

Quality Assurance

Sartorius Stedim Biotech Quality Systems for Single-Use Products is compliant with ISO 9001 : 2008. Design, Manufacture and Sterilization processes are conducted under conditions that mirror biopharmaceutical operations and meet cGMP requirements.

Available from Stock

TuFlux® SIL coils are available from stock.

Validation

TuFlux® SIL has been extensively qualified in collaboration with Raumedic applying the most stringent tests. Biological, chemical and physical tests combined with extensive extractable testing provide users of TuFlux® SIL a silicone tubing with data representing a variety of processing conditions covering a lot of applications.

When tubing is mounted on a product, a full compliance with ISO11137 allows for a validated claim of sterility by Gamma irradiation on all Sartorius Stedim Biotech single-use products with a sterility assurance level of 10⁻⁶ over the shelf life.

TuFlux® SIL is tested for compliance to:

- E.P. 3.1.9
- USP <88>: Class VI Implant test, systemic toxicity tests and intracutaneous tests
- Cytotoxicity according to ISO 10993-5
- Haemolysis test according to ISO 10993-4
- USP <381>: Test for elastomeric closures for injections
- USP <661>: Tests for plastic
- FDA 21 CFR 177.2600
- Animal Derived Component Free
- REACH

Details on methodologies and equipment used as well as further tests performed are detailed in the Validation Guide.



Ordering Information

Part Number	Description	ID × OD Dimensions (mm)	Wall thickness (mm)	Tubing coil length (m)	Minimum Order Quantity
FSA117425	TuFlux [®] SIL 1/8" × 1/4"	3.2 × 6.4	1.6	100	1 coil
FSA117426	TuFlux [®] SIL 1/4" × 3/8"	6.4 × 9.5	1.6	50	1 coil
FSA117427	TuFlux [®] SIL 1/4" × 7/16"	6.4 × 11.1	2.4	50	1 coil
FSA117428	TuFlux [®] SIL 3/8" × 5/8"	9.5 × 15.9	3.2	25	1 coil
FSA117429	TuFlux [®] SIL 1/2" × 3/4"	12.7 × 19.1	3.2	20	1 coil
FSA117430	TuFlux [®] SIL 3/4" × 1"	19.1 × 25.4	3.2	15	1 coil
FSA117431	TuFlux [®] SIL 3/4" × 1 1/8"	19.1 × 28.6	4.8	15	1 coil

Part Number	Description	ID × OD Dimensions (inch)	Wall thickness (inch)	Tubing coil length (ft)	Minimum Order Quantity
FSA117425	TuFlux [®] SIL 1/8" × 1/4"	1/8" × 1/4"	0.06	328	1 coil
FSA117426	TuFlux [®] SIL 1/4" × 3/8"	1/4" × 3/8"	0.06	164	1 coil
FSA117427	TuFlux [®] SIL 1/4" × 7/16"	1/4" × 7/16"	0.09	164	1 coil
FSA117428	TuFlux [®] SIL 3/8" × 5/8"	3/8" × 5/8"	0.13	82	1 coil
FSA117429	TuFlux [®] SIL 1/2" × 3/4"	1/2" × 3/4"	0.13	66	1 coil
FSA117430	TuFlux [®] SIL 3/4" × 1"	3/4" × 1"	0.13	49	1 coil
FSA117431	TuFlux [®] SIL 3/4" × 1 1/8"	3/4" × 1 1/8"	0.19	49	1 coil

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
Modectcenterstrasse 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 A. Meucci, 4
50012 Grassano - 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

Africa

Republic of South Africa

Sartorius South Africa (Pty) Ltd
Unit 4, Alphen Square South
853 16th Road
Midrand 1685, RSA
Phone +27.11.315.5444

Americas

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
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