

Tuflux® SIL Braided

Sartorius Pressure Resistant Silicone Tubing



Product Information

Sartorius Tuflux® SIL Braided is designed to facilitate fluid transfer in the biopharmaceutical manufacturing processes with high pressure requirements. Tuflux® SIL reinforced tubing is a highly resistant platinum cured silicone tubing with a medical grade PET mesh material manufactured by Raumedic. It is available in six different internal dimensions from 1/8" (3.2 mm) to 1" (25.4 mm) and from 0.11" (2.9 mm) to 0.19" (4.8 mm) in wall thickness.

Applications

Tuflux® SIL Braided is designed to be used in many pharmaceutical and biopharmaceutical applications such as:

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

Flexibility

Sartorius Tuflux® SIL Braided is available as non-sterile tubing coils and can also be assembled on sterilized finished products. Several dimensions are available to match process requirements in terms of flow rate.

Features and Benefits

- Pressure resistance up to 60 bar depending on tube diameter
- Medical grade mesh reinforcement enabling high pressure resistance while ensuring fluids are kept safe
- Fully Validated Tubing:
 - Same platinum cured silicone tubing as Tuflux® SIL reducing validation effort
 - Available document package to ease integration into processes:
 - Validation guide
 - Extractable guide
 - BPOG reports
 - Comparability guide
 - Certificate of release delivered with each order
- Includes all the benefits of Tuflux® SIL: purity and cleanliness due to its low extractable profile, robustness due to its hardness, easy to handle due to its unique low-tack treatment



Technical Data

Specifications

Colour	Translucent
Material	Platinum cured silicone braided with medical grade PET mesh material
Hardness	60 ± 5 Shore A
Temperature range	-20 °C to + 135 °C (-4 °F to + 275 °F)
Tensile strength	> 8.0 MPa
Elongation at break	> 500%
Compliant	E.P. 3.1.9, USP <88> Class VI, ADCF
Printing	Biocompatible and inert ink showing tubing dimensions
Sterilization	Gamma irradiation and autoclaving
Tubing coils packaging	Primary and secondary PE packaging
Surface treatment	Low tack treatment to reduce tube stickiness and surface friction

Tubing Coil

Tuflux® SIL Braided is available from stock as non-sterile tubing coil double-packaged in PE bags. The length of a coil depends on the tubing dimensions and varies between 10 m to 25 m.

Tuflux® SIL Braided is compatible with gamma sterilization and autoclaving. The shelf life of the non-sterile coils is five years.

Quality Assurance

Sartorius Quality Systems for Single-Use Products is compliant with ISO 9001. Design, manufacture and sterilization processes are conducted under conditions that mirror biopharmaceutical operations and meet cGMP requirements.

Available From Stock

Tuflux® SIL Braided coils are available from stock.

Validation

Braided Tuflux® SIL has been extensively qualified in collaboration with Raumedic. Biological, chemical and physical tests combined with extensive extractable testing provide users of Tuflux® SIL Braided with data representing a large variety of processing conditions.

Tuflux® SIL Braided is compliant to:

- E.P. 3.1.9
- USP<88> Class VI: Biological reactivity tests, in-vivo
- USP<85>: Bacterial endotoxins testes
- USP<87>: Biological reactivity tests, in-vitro (equivalent to ISO10993-5)
- Haemolysis test according to ISO 10993-4
- USP<661>: Containers-Plastics tests
- Animal Derived Component Free (ADCF)
- REACH

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

Pressure Resistance Data

Part Number	ID × OD Dimensions [inch]	Water Burst Pressure [bar]	Maximum Recommended Working Pressure For Single Tube [bar]
FSA309865	⅛" × 0.355"	≥ 60	24
FSA309864	¼" × ½"	≥ 43	17
FSA309863	⅜" × ⅝"	≥ 39	15
FSA309862	½" × ⅞"	≥ 29	11
FSA309861	¾" × 1 - ⅛"	≥ 18	7
FSA309860	1" × 1 - ⅜"	≥ 11	4

Order Information


Part Number	Description	ID × OD Dimensions [mm]	Wall Thickness [mm]	Tubing Coil Length [m]	Minimum Order Quantity
FSA309865	Tuflux® SIL Braided ⅛" × 0.355"	3.2 × 9.0	2.95	25	1 coil
FSA309864	Tuflux® SIL Braided ¼" × ½"	6.4 × 12.7	3.20	15	1 coil
FSA309863	Tuflux® SIL Braided ⅜" × ⅝"	9.5 × 15.9	3.20	15	1 coil
FSA309862	Tuflux® SIL Braided ½" × ⅞"	12.7 × 19.1	4.70	15	1 coil
FSA309861	Tuflux® SIL Braided ¾" × 1 - ⅛"	19.1 × 28.6	4.80	10	1 coil
FSA309860	Tuflux® SIL Braided 1" × 1 - ⅜"	25.4 × 34.9	4.80	10	1 coil

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