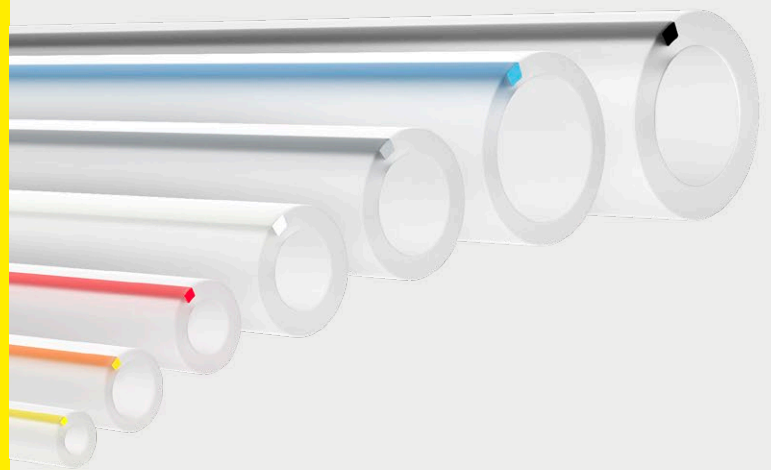


Tuflux[®] TPE

Sartorius Sealable and Weldable TPE Tubing



Description

Sartorius Tuflux[®] TPE is designed to facilitate fluid transfer in the biopharmaceutical manufacturing processes. Tuflux[®] TPE is a highly resistant sealable and weldable tubing manufactured by Raumedic AG. Tuflux[®] TPE is available in 7 different internal dimensions from 1/8" (3.2 mm) to 3/4" (19.1 mm) and from 1.6 to 4.8 mm in wall thickness.

Sartorius Tuflux[®] TPE is available as non-sterile tubing coils and can also be assembled on sterilized finished products.

Product Information

Tuflux[®] TPE is a translucent thermoplastic elastomer tubing that guarantees an easy identification due to the embedded color stripe aligned with corresponding tube holder for safe welding applications using Biowelder[®] TC devices as well as with Clipster[®] color-coding for aseptic disconnections. Tuflux[®] TPE has been designed and validated to be compatible with cell-growth and it can be co-welded to established TPE brands in the market with pre-validated welding parameters under sterile conditions.

Applications

Tuflux® TPE is designed to be used in pharmaceutical and biopharmaceutical applications such as:

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

Features & Benefits



- **Unique color coding**, aligned with Biowelder® TC tube holders, and printing of ID and OD dimensions on the tubing facilitates tubing **identification and ease of use** to mitigate operator errors.



- The antioxidant package in the tubing formulation is **optimized to obtain the best cell growth results**. No inhibition of cell growth even under worst case surface to volume ratios in small scale applications



- Validated Biowelder® TC parameters for **co-welding of Tuflux® TPE** to other TPE tubes. Welding parameters have been **validated for all combinations of gamma-irradiated & autoclaved TPE tubes**.

- **Validated for sealing** with the Biosealer® TC for sterile disconnections in selected sizes.
- **Validation & Extractable Guides** for an easy integration of Tuflux® TPE products in your processes.
- **Tuflux® TPE is available from stock** as non-sterile tubing coil double packaged in PE bags.
- **Tuflux® TPE is suitable to be used with peristaltic pumps** for liquid transfer applications.

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® TPE coils are available from stock.

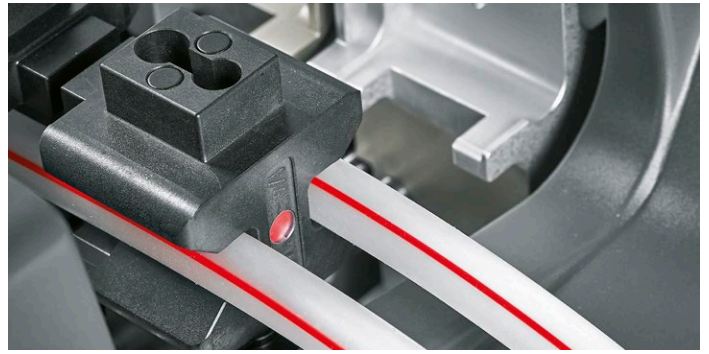
Validation

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

Technical Data

Specifications

Color	Translucent
Material	Synthetic thermoplastic elastomer
Hardness	60 ± 5 Shore A
Tensile strength at break	≥ 5.0 MPa
Elongation at break	≥ 450%
Temperature range	-55°C to 135°C
Shelf Life	5 years non sterile 3 years sterile
Compliant to	<ul style="list-style-type: none"> ▪ USP<87> & USP<88> Class VI ▪ ISO 10993-5 ▪ ADCF ▪ USP<661> ▪ REACH ▪ 2011/65/EC for heavy metals
Sterilization	Gamma irradiation (max 50 kGy) and autoclave (max 123 °C, 30 min)



Ordering Information

Part Number	Product Description	Coil [m ft]	Diameter [ID × OD]	ID [mm inch]	WT [mm inch]	Color stripe
FSA121869	Tuflux® TPE 1/8"ID × 1/4"OD yellow	200 m 656 ft	1/8" × 1/4"	3.2 0.125	1.6 0.06	Yellow
FSA121870	Tuflux® TPE 1/4"ID × 3/8"OD orange	125 m 410 ft	1/4" × 3/8"	6.4 0.25	1.6 0.06	Orange
FSA121871	Tuflux® TPE 1/4"ID × 7/16"OD red	100 m 328 ft	1/4" × 7/16"	6.4 0.25	2.4 0.09	Red
FSA121872	Tuflux® TPE 3/8"ID × 5/8"OD white	50 m 164 ft	3/8" × 5/8"	9.5 0.375	3.2 0.125	White
FSA121873	Tuflux® TPE 1/2"ID × 3/4"OD grey	30 m 98 ft	1/2" × 3/4"	12.6 0.5	3.2 0.125	Grey
FSA121874	Tuflux® TPE 3/4"ID × 1"OD blue	20 m 66 ft	3/4" × 1"	19.1 0.75	3.2 0.125	Blue
FSA121875	Tuflux® TPE 3/4"ID × 1 1/8"OD black	20 m 66 ft	3/4" × 1 1/8"	19.1 0.75	4.8 0.188	Black




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

Specifications subject to change without notice.

*2023 Sartorius Stedim FMT S.A.S., Avenue de Jouques - Zone Industrielle Les Paluds - C.S. 91051, 13781 Aubagne Cedex, France
Order No.: 1000049950 | Status: 05 | 2023