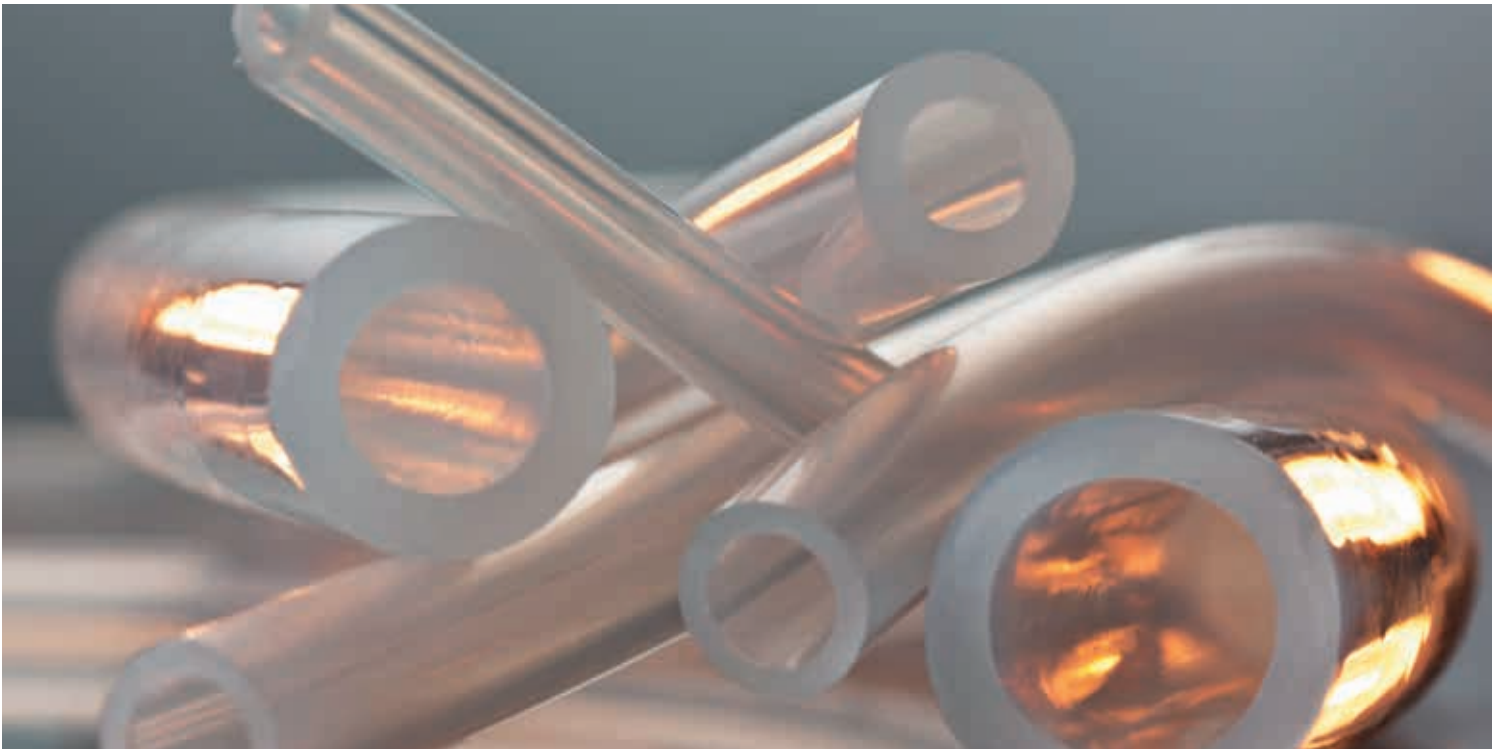




sartorius stedim
biotech

Comparability Guide for TuFlux® SIL Tubing (Sartorius Stedim Biotech) vs Pharma 50 (Dow Corning)



Applicable to:
TuFlux® SIL tubing used for the
manufacture of Flexboy®, Flexel®, and
CultiBag®, as well as standalone Coils and
all other Bag Systems manufactured by
Sartorius Stedim Biotech.

The results shown in this Comparability
Protocol and Equivalency Test Report are
indicative and do not constitute product
specifications.

This confidential document including
any attachment contains confidential
and proprietary information of Sartorius
Stedim Biotech.

Disclosure, copying or distribution to any
third party of the information included in
this document is prohibited without the
prior written consent of Sartorius Stedim
Biotech.

To obtain authorization, please contact
Sartorius Stedim Biotech through your
regular contact person.

Table of Contents

1.	Introduction	4
2.	General Features	5
3.	Biocompatibility	7
4.	Physico-chemical Properties	9
5.	Barrier Properties to Water	11
6.	Pressure Resistance	12
7.	Pumping Life Time	13
8.	Flow Rates	14
9.	Tubing Benchmarking after Recirculation	15
9.1	Pumping Time vs pH	15
9.2	Pumping Time vs Conductivity	15
9.3	Pumping Time vs TOC Content	16
9.4	Pumping Time vs Particle Content	16
10.	Extractable Studies	17
10.1	Results from Exposure of Silicone Tubing to WFI Solution	18
10.1.1	Volatiles by Headspace-GC MS	18
10.1.2	Semi Volatile Compounds (GC MS)	18
10.1.3	Metal Analysis (ICP-MS OES)	18
10.2	Results from Exposure of Silicone tubing to NaOH 1M Solution	19
10.2.1	Semi Volatile Compounds (GC MS)	19
10.2.2	Metal Analysis (ICP-MS OES)	19
10.3	Results from Exposure of Silicone Tubing to HCl 1M Solution	20
10.3.1	Semi Volatile Compounds (GC MS)	20
10.3.2	Metal Analysis (ICP-MS OES)	20
10.4	Results from Exposure of Silicone Tubing to EtOH Solution	21
10.4.1	Semi Volatile Compounds (GC MS)	21
10.4.2	Metal Analysis (ICP-MS OES)	21

1. Introduction

Sartorius Stedim Biotech bags and systems are widely used in biopharmaceutical processes for a variety of unit operations of the commercial production of drug products such as vaccines, recombinant proteins and monoclonal antibodies and for the development of future biomolecules in clinical phases.

Buffers and media are increasingly formulated, sterile filtered and stored in single-use Fluid Management Systems (FMS) that involve Flexel® and Flexboy® Bags integrated with filters, impeller mixers, tubing, connectors and monitoring tools. Product intermediates are also filtered and stored between UF|DF and chromatography purification steps in Gamma sterile Fluid Management Systems.

Fluid Management Systems are also adopted for the formulation, filtration and aseptic processing of final drug products. From buffer media preparation, cell culture operations, purification operations up to final formulation, filtration and transfer, the silicone tubing TuFlux® SIL is a key element for the successful implementation of disposable manufacturing processes.

The TuFlux® SIL, Sartorius Stedim Biotech silicone tubing, is qualified, manufactured and released according to stringent product validation protocols and Quality Control testing to offer safe and robust single-use processes to the end users of the biopharmaceutical industry.

This Guide is to show comparability and variances between Platinum cured Silicone tubing from different suppliers (e.g. Raumedic and Dow Corning).

TuFlux® SIL will be the Sartorius Stedim Biotech standard tubing and replace the Dow Corning Pharma 50 in all new Fluid management Systems

2. General Features

Physical Properties

The following values are determined on standard test specimens punched from a press plate.

	Colour	Hardness Shore A	Operating Temperature Range	Ultimate Tensile Strength (according to ISO 527)	Elongation at Break (according to ISO 527)
TuFlux® SIL (Sartorius Stedim Biotech)	Translucent	60 ± 5	-60 °C to +200 °C (-76 °F to +392 °F)	> 8 MPa	> 600%
Pharma 50 (Dow Corning)	Translucent	50 ± 5	-51 to 232 °C (-60 °F to 460 °F)	> 6.9 MPa	> 650%

Material Hardness:

Purpose and Test Method

A measure of the indentation resistance of elastomeric or soft plastic materials based on the depth of penetration of a conical indenter. Hardness values range from 0 (for full penetration) to 100 (for no penetration).

Tensile Properties:

Purpose and Test Method

A tensile test consists in applying an elongation to a tubing specimen and measuring the resulting strength. Mechanical properties can then be defined from the stress-strain curve.

Ultimate Tensile Strength (UTS):

The maximum stress, a material can withstand, is calculated by dividing the maximum load by the original cross sectional area of the specimen. The tensile strength test is performed with a tensile machine in stretching or elongation mode.

Elongation at Break:

The elongation is recorded at the moment of specimen rupture and often expressed as a percentage of the original length. Materials with high elongation at break withstand a high deformation before rupture. A high elongation at break often means high flexibility.

	Printing of ID and OD Dimensions on the Tubing	Coils Wrapped in Double Packaging	Low-Tack	Post Curing	Sterilization Resistance
TuFlux® SIL (Sartorius Stedim Biotech)	Yes	Yes	Yes	Yes	γ-irradiation Autoclave ETO
Pharma 50 (Dow Corning)	Not for Sartorius Stedim Biotech	Yes	No	Yes	γ-irradiation Autoclave ETO

Tubing Printing

"a|b × c|d" where a|b is the Internal Dimension and c|d the Outer Dimension of the tubing in inch.

The innocuity of the ink is proven on printed tubing by the tests performed according to ISO 10993-5.

Low Tack

The surface is coated in a plasma process. This coating provides a less sticky surface of these silicone tubing (Low Tack) in comparison with common non-coated silicone products.

3. Biocompatibility

Silicone tubing meets the requirements of following biocompatibility tests:

	E.P. 3.1.9	USP <88> Class VI	USP <87>	USP <85>	Other ISO
TuFlux® SIL (Sartorius Stedim Biotech)	Yes	Yes	Yes	Yes	ISO10993-4 ISO10993-5
Pharma 50 (Dow Corning)	Yes	Yes	No info	No info	No info

Purpose and Test Method

Biocompatibility tests are performed to demonstrate that the Tubing is biocompatible and meets or exceeds the current USP and ISO requirements.

Tests are carried out on Tubing samples before and after Gamma irradiation (50 kGy). The tubing samples were supplied to an independent testing facility for evaluation under the current USP <88> and ISO 10993-5 biocompatibility standards.

USP <88> Class VI

Tubing material is implant tested, meets the requirements of implant test to USP <88> Class VI as well as the intracutaneous test and the acute toxicity test to USP <88> Class VI, meaning that biological neutrality has been proven via these animal experiment tests on sterile and Gamma or autoclave sterilized samples. The following tests were performed on samples with and without ink: Cytotoxicity test intracutaneous test Systemic injection test Implantation test (7 days).

Test Results of USP <88> Class VI

All material used in the construction of the silicone tubing meet or exceed the requirements of the USP <88> Class VI- 121 °C Plastics tests and are considered as non cytotoxic and non haemolytic.

USP <87> – in vitro compared to USP <88> in vivo

Physico Chemical Test

All testing presented in this comparability guide have been performed on gamma irradiated tubing at 50 kGy that represents the maximum doses. If different the tubing conditions will be precised.

European Pharmacopoeia: 3.1.9

Purpose and Test Method

With regard to their basic material, additives and properties, tubing in silicone rubber has been tested in compliance with the recommendations and guidelines on health assessment of plastics within the scope of Food and Drug Act, sections A XV and B II XV, European Pharmacopoeia (E.P.) 3.1.9. as well as FDA regulation 21 CFR, § 177.2600.

Test Results

Tubing in silicone rubber meets the requirements of the European Pharmacopoeia 3.1.9. and regulation 21 CFR, § 177.2600.

The test methods, limits and results are those described by the E.P. monograph and listed in the table below.

Test Description	E.P. 3.1.9 Limits
Appearance of solution	Colourless
Acidity	≤ 2.5 mL NaOH 0.01M
Alkalinity	≤ 1.0 mL HCl 0.01M
Reducing substances	< 1 mL
Substance soluble in hexane	< 3%
Volatile matter	< 2%
Mineral oils	< 1 ppm
Platinum	< 30 ppm

4. Physico-chemical Properties

	USP <381> compliant	USP <661> compliant	ADCF	REACH compliant	FDA 21 CFR 177.2600 compliant
TuFlux® SIL (Sartorius Stedim Biotech)	Yes	Yes	Yes	Yes	Yes
Pharma 50 (Dow Corning)	Yes	Yes	No info	Yes	Yes

USP <381>

Elastomeric Closures for Injections

Purpose and Test Method

Elastomeric closures for containers are made of materials obtained by vulcanization (cross-linking) polymerization, polyaddition, or polycondensation of macromolecular organic substances (elastomers). Elastomeric closures shall conform to biological, physicochemical, and functionality requirements. The tests are performed according to USP <381> recommendations.

Test Description	USP <381> Limits
Chearlessness	
Reducing substances	0
Lead	0 ppm
pH value change	5-7
Dry residue	< 2 mg

USP <661>

**Containers, Physicochemical Tests
– Plastic Purpose**

Physicochemical tests are designed to determine physical and chemical properties of TuFlux® SIL tubing and their extracts. They are performed on TuFlux® SIL samples before and after irradiation and accelerated ageing conditions.

Test Method

Tubing samples were cut in small portions previously Gamma irradiated (50 kGy) and extracted (ratio of 30 g per 150 mL) in UltraPure water at 70 °C (158 °F) for 24 hours. The tests are conducted in order to determine physical and chemical properties of the testarticle and its extracts. The same test have been performed on tubing Gamma irradiated (50 kGy) stored during a period corresponding to a shelf life of 3 years (accelerated conditions).

Test Results

The silicone tubing TuFlux® SIL meets the USP <661> requirements when sterilized at 50 kGy with and without aging conditions corresponding to a shelf life of 3 years.

Test Description	USP <661> Limits
Non-volatile residue	< 15 mg
Residue in ignition	< 5 mg
Heavy metals	< 1 ppm
Buffering capacity	< 10 mL

ADCF Certified

Tubing Material does not contain any animal derived components.

REACH

Tubing Material is free from any substances defines as SVHC – Substances of Very High Concern – by the European REACH regulation, Annex XIV.

21 CFR, § 177.2600

With regard to their basic material, additives and properties, tubing in silicone rubber is in compliance with the recommendations and guidelines on health assessment of plastics within the scope of Food and Drug Act, sections A XV and B II XV, as well as is FDA regulation 21 CFR, § 177.2600 compliant.

5. Barrier Properties to Water

The aim of this test is to evaluate the permeability of silicone tubes irradiated at 50 kGy to WiFi at 60 °C (140 °F) for 2, 7 and 14 days representing 1, 3 and 6 months respectively in normal room temperature conditions.

Tube	Mass loss at 60 °C [%]			pH (calculated*)		Conductivity [μ S/cm] (calculated*)	
	t = 2d	t = 7d	t = 14d	t = 14d (blank)	t = 14d	t = 14d (blank)	t = 14d
TuFlux® SIL (1/2" × 3/4")	5.4	19.0	38.0		3.5		190.1
Pharma 50 (1/2" × 3/4")	6.5	21.0	40.9	7.4	3.8	4.6	70.6

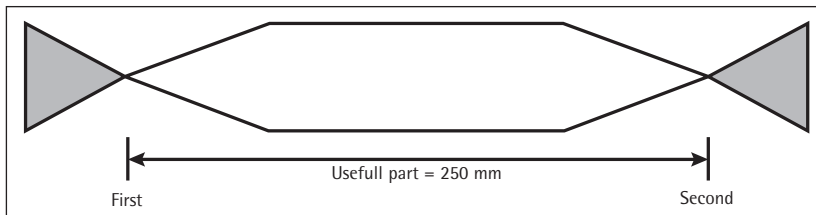
Note a : means that samples have been diluted 3 × prior to pH and conductivity measurements. Therefore, pH and conductivity have been extrapolated -"calculated"- to their initial values taking into account the dilution.

Barrier Properties to Water Vapor

The mass of each assembly is controlled at start and after 2, 7 and 14 days the samples being stored at 60 °C (140 °F). The percentage of mass loss is calculated from the weight at start of the test compared to the weight after storage time.

pH and Conductivity

pH and conductivity are measured at t = 0 and after 14 days of storage at 60 °C (140 °F).



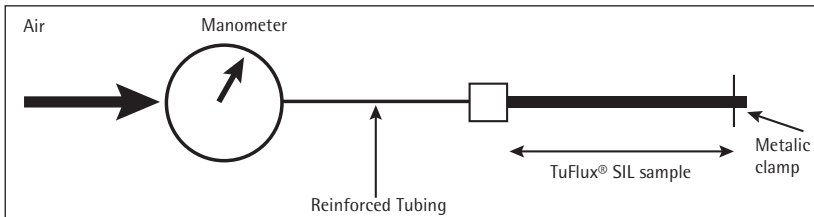
6. Pressure Resistance

Purpose

The goal of the burst pressure test is to assess the pressure resistance of the tubing depending on the tubing dimensions (inside and outside diameter).

Test Method

The test method is described in the scheme below. Three measurements of the pressure were taken for each tubing reference on Gamma sterilized samples at 50 kGy at burst.



Pressure Resistance Results:

pressure at which tubings inflate up to 10% of their initial dimensions

Dimension ["] (ID/OD)	Dimension [mm]	TuFlux® SIL Air Pressure Limit Resistance [bar]	Pharma 50 Air Pressure Limit Resistance [bar]
1/8" × 1/4"	3.2 × 6.4	4.5	3.8
1/4" × 3/8"	6.4 × 9.6	3.1	2.4
1/4" × 7/16"	6.4 × 11.1	3.5	3.3
3/8" × 5/8"	9.6 × 15.9	4.0	3.6
1/2" × 3/4"	12.7 × 19.1	4.6	2.4
3/4" × 1"	19.1 × 25.4	3.1	1.8
3/4" × 1 1/8"	19.1 × 28.6	4.7	2.4

7. Pumping Life Time

Purpose

The goal of the pumping life time test is to assess the mechanical resistance of the tubing under pumping conditions.

Test Method

The tubing is placed in a Watson Marlow serie 720 peristaltic pump and speed is set up to maximum (> 300 rpm). The tubing was pumping water at ambient temperature between 2 tanks mimicking recirculation conditions.

The test was stopped and time measured at tubing break resulting in leak.

Tubing	Pumping Life Time [h]
TuFlux® SIL	> 70
Pharma 50	> 55

8. Flow Rates

Purpose

The objective of this test was to assess the maximum flow rate of some TuFlux® SIL dimensions.

Test Method

The time to transfer 100 L of water at room temperature with non sterile tubing using a peristaltic pump set up at the maximum speed (310 rpm) was measured in duplicates.

Tubing Size (ID × OD)	Tubing Material	
	TuFlux® SIL	Pharma 50
1/8" × 1/4" 3.2 × 6.4 mm (WT 1.6 mm)	> 0.3 L/min	-
3/8" × 5/8" 9.5 × 15.9 mm (WT 3.2 mm)	> 6.0 L/min	> 4.9 L/min
1/2" × 3/4" 12.7 × 19.05 mm (WT 3.2 mm)	> 8.9 L/min	> 8.0 L/min

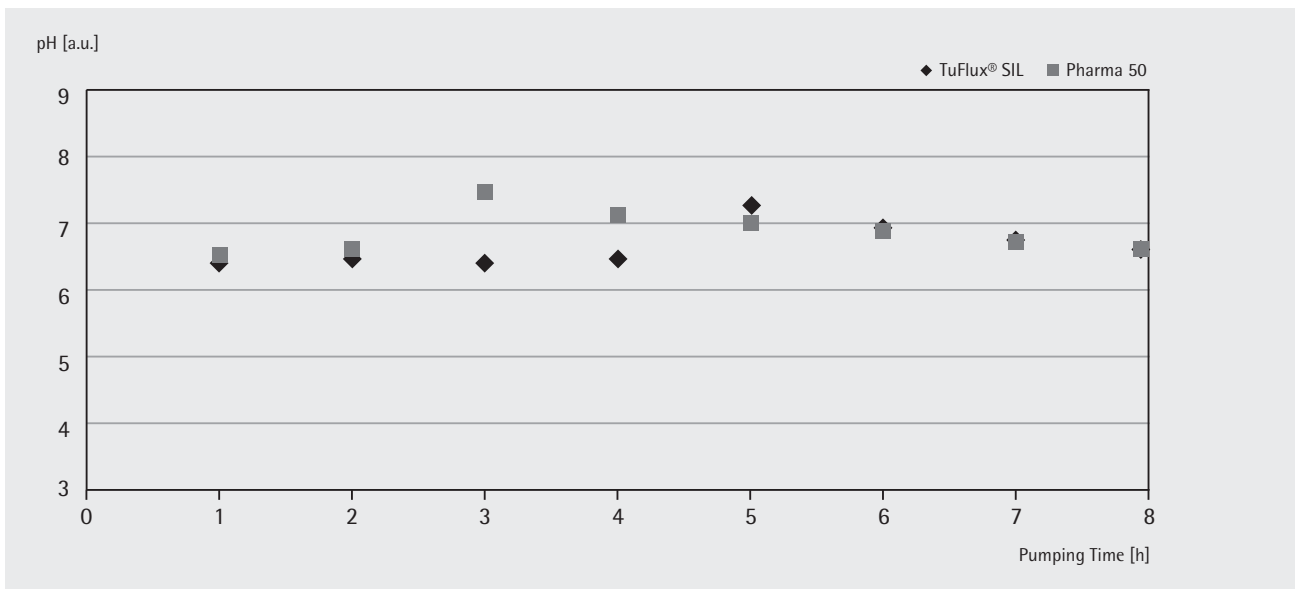
9. Tubing Benchmarking after Recirculation

The goal of this test is to compare tubes irradiated at 50 kGy about the influence of the pumping time on pH, conductivity, particulates, UV|Vis and TOC.

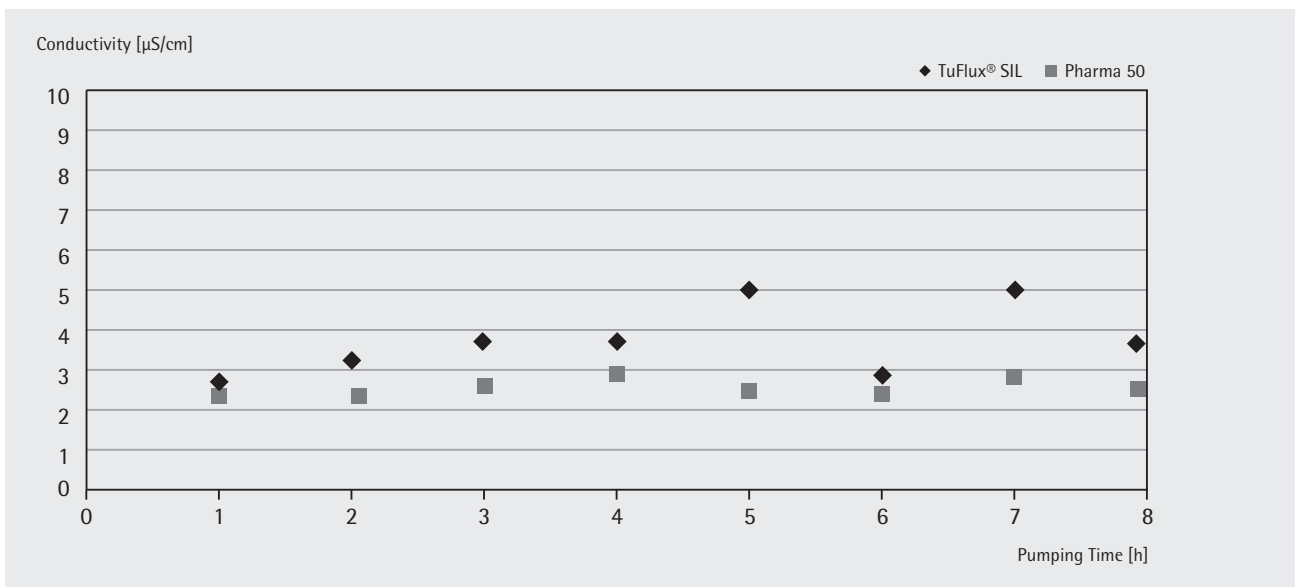
Tests Conditions:

- Tube length: 2 m of $\frac{1}{2} \times \frac{3}{4}$ " Tubing
- Volume WFI: 3 L
- Glass flask + plastic tube
- Rotation speed of the pump: 310 rpm
- Samples are taken at $t = 0$ and after x hours of pumping (x = 1, 2, 3, 4, 6, 8, 10).

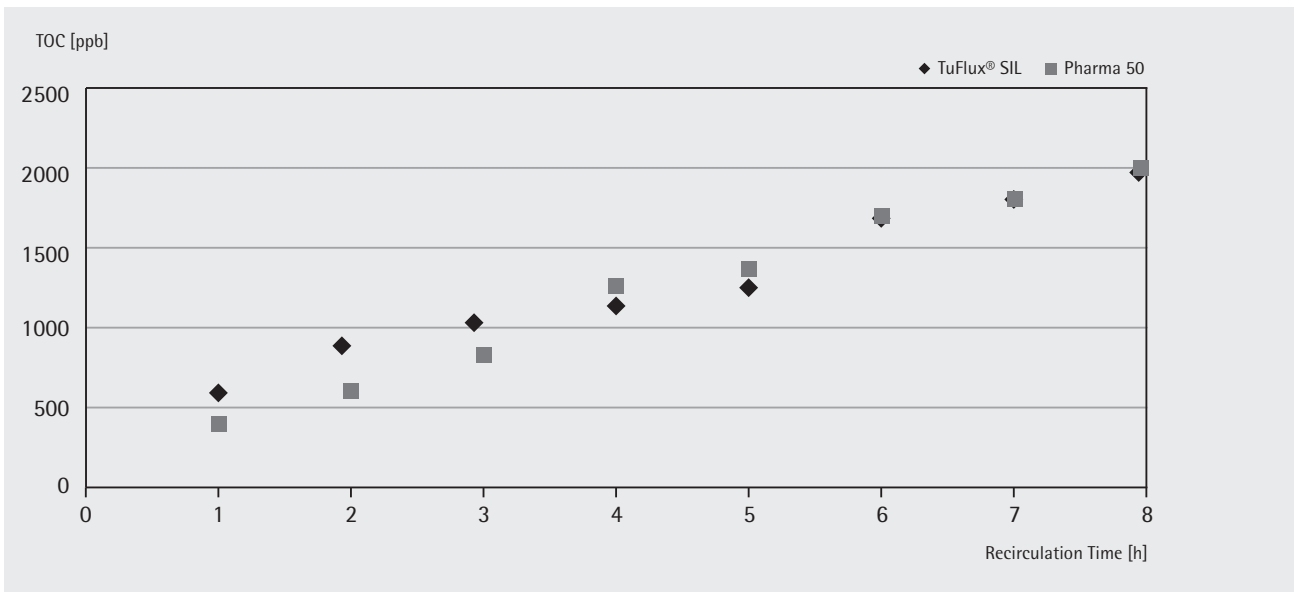
9.1 Pumping Time vs pH



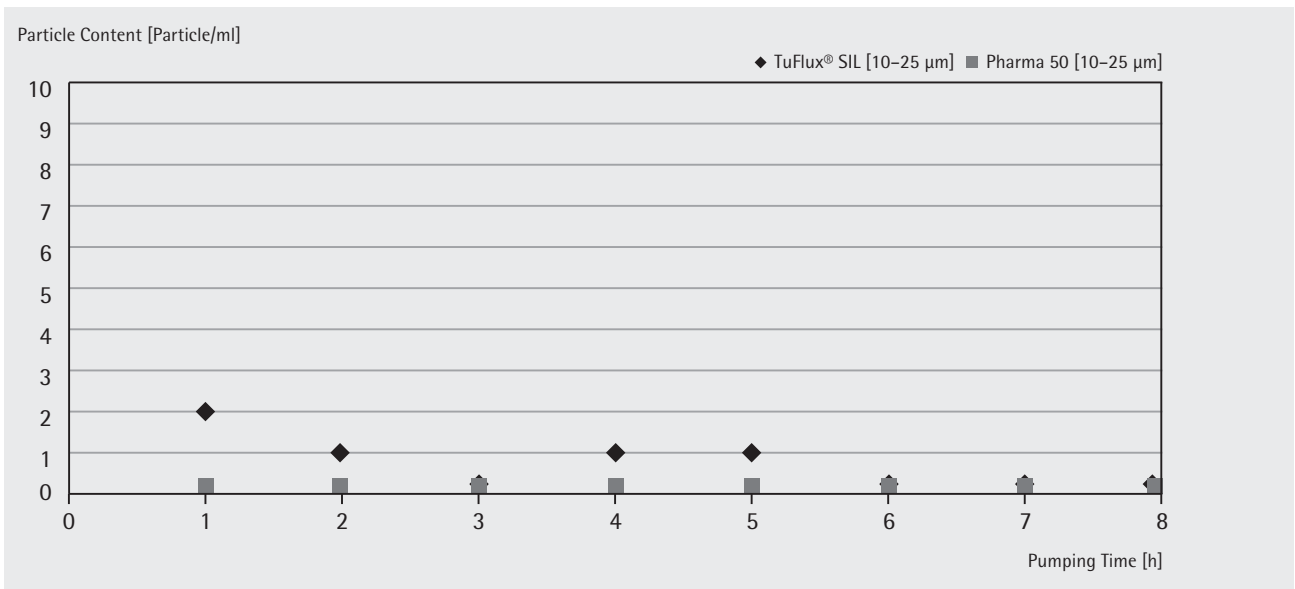
9.2 Pumping Time vs Conductivity



9.3 Pumping Time vs TOC Content



9.4 Pumping Time vs Particle Content

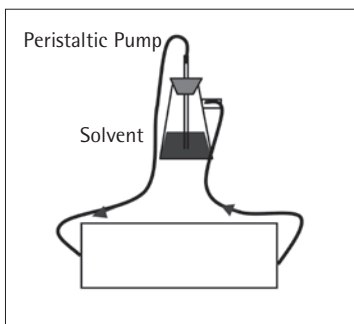


10. Extractable Studies

The comparison of tubing materials has been performed by solvent recirculation in tubing at room temperature:

- with EtOH for 2H-
- with 1M-HCl, 1MNaOH, WFI for 10H

→ Experiments as follow:



Purpose

Extractables are substances that can be extracted from a bioprocessing containment, such as a tubing, using extraction solvents and conditions that are more aggressive than the usual conditions of use. The goal of this extractable profile test is to supply worst case extractable data to support process developers and toxicologists in their validation studies.

Reporting Limits for the Different Analytical Methods:

- Volatile Compounds (HS-GC|MS):
Reporting limit: 0.01 µg/mL
- Non Volatile Compounds (LC|MS):
Reporting limit: 0.01 µg/mL
(for WFI only)
- Metal Analysis (ICP-MS|OES):
Reporting limit: 0.1 µg/mL
- Semi Volatile Compounds (GC|MS) :
Reporting limit: 0.05 µg/mL

The Data shown below represents cleaned Data.

Only values which are three fold above the blank are reported.

10.1 Results from Exposure of Silicone Tubing to WFI Solution

10.1.1 Volatiles by Headspace-GC|MS

Analytes	Quantification [$\mu\text{g/mL}$]	
	TuFlux® SIL	Pharma 50
Ethanol	0–0.05	0–0.05
Dichloromethane	0.05–0.1	0–0.05

Ethanol and dichloromethane are surely cross contaminant chemicals during sample preparation.

10.1.2 Semi Volatile Compounds (GC|MS)

Analytes	Quantification [$\mu\text{g/L}$]	
	TuFlux® SIL	Pharma 50
(None)		

10.1.3 Metal Analysis (ICP-MS|OES)

Solvent	Analytes	Quantification [$\mu\text{g/L}$]	
		TuFlux® SIL	Pharma 50
WFI	Phosphorus	0.1–0.5	0.1–0.5
	Silicone	0.1–0.5	0.1–0.5
	Sodium	1–5	1–5

10.2 Results from Exposure of Silicone Tubing to NaOH 1M Solution

10.2.1 Semi Volatile Compounds (GC|MS)

Analytes	Quantification [$\mu\text{g/L}$]	
	TuFlux® SIL	Pharma 50
Aromatic hydrocarbon	0	0

10.2.2 Metal Analysis (ICP-MS|OES)

Solvent	Analytes	Quantification [$\mu\text{g/L}$]	
		TuFlux® SIL	Pharma 50
NaOH 1M	Aluminium	1–5	n.d.
	Boron	5–10	n.d.
	Silicone	50–100	50–100

n.d. = not detected

10.3 Results from Exposure of Silicone Tubing to HCl 1M Solution

10.3.1 Semi Volatile Compounds (GC|MS)

Analytes	Quantification [$\mu\text{g/L}$]	
	TuFlux® SIL	Pharma 50
None	-	-

10.3.2 Metal Analysis (ICP-MS|OES)

Solvent	Analytes	Quantification [$\mu\text{g/L}$]	
		TuFlux® SIL	Pharma 50
HCl 1M	Calcium	0.1–0.5	0.1–0.5
	Magnesium	0.1–0.5	0.5–1
	Silicone	1–5	1–5
	Sodium	0.5–1	n.d.
	Zinc	1–5	5–10

n.d. = not detected

10.4 Results from Exposure of Silicone Tubing to EtOH Solution

10.4.1 Semi Volatile Compounds (GC|MS)

Analytes	CAS Number	Quantification [µg/L]	
		TuFlux® SIL	Pharma 50
Octamethyl Cyclotetrasiloxane NaOH 1M	[566-67-2]	0.5-1	n.d.
Decamethyl Cyclopentasiloxane	[541-02-6]	1-5	5-10
Dodecamethyl Cyclohexasiloxane	[540-97-6]	1-5	10-50
Tetradecamethyl Cycloheptasiloxane	[107-50-6]	1-5	10-50
Hexadecamethyl Cyclooctasiloxane	[556-68-3]	10-50	10-50
Octadecamethyl Cyclononasiloxane	[556-71-8]	1-5	10-50
Other siloxanes	-	1-50	100-250

10.4.2 Metal Analysis (ICP-MS|OES)

Solvents	Analytes	Quantification [µg/L]	
		TuFlux® SIL	Pharma 50
Ethanol	Silicone	50-100	100-250
	Zinc	0.1-0.5	n.d.

Solvents = Ethanol
n.d. = not detected

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
Fax +49.551.308.3289

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

Phone +49.5665.407.0
Fax +49.5665.407.2200

France

Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques – CS 91051
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.845619

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques – CS 71058
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.846545

Austria

Sartorius Stedim Austria GmbH
Franzosengraben 12
1030 Vienna

Phone +43.1.7965763.18
Fax +43.1.796576344

Belgium

Sartorius Stedim Belgium N.V.
Leuvensesteenweg, 248/B
1800 Vilvoorde

Phone +32.2.756.06.80
Fax +32.2.756.06.81

Hungary

Sartorius Stedim Hungária Kft.
Kagyló u. 5
2092 Budakeszi

Phone +36.23.457.227
Fax +36.23.457.147

Italy

Sartorius Stedim Italy S.p.A.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)

Phone +39.055.63.40.41
Fax +39.055.63.40.526

Netherlands

Sartorius Stedim Netherlands B.V.

Phone +31.30.60.25.080
Fax +31.30.60.25.099

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
Fax +48.61.879.25.04

Russian Federation

LLC "Sartorius ICR"
Uralskaya str. 4, Lit. B
199155, Saint-Petersburg

Phone +7.812.327.5.327
Fax +7.812.327.5.323

Spain

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

Phone +34.902.110.935
Fax +34.91.358.96.23

Switzerland

Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen

Phone +41.52.354.36.36
Fax +41.52.354.36.46

U.K.

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

Phone +44.1372.737159
Fax +44.1372.726171

America

USA

Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716

Toll-Free +1.800.368.7178
Fax +1.631.254.4253

Argentina

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

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

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

Phone +55.11.4362.8900
Fax +55.11.4362.8901

Mexico

Sartorius de México S.A. de C.V.
Circuito Circunvalación Poniente
No. 149
Ciudad Satélite
53100, Estado de México
México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

Asia|Pacific

Australia

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

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius Stedim Biotech (Beijing) Co. Ltd.
No. 33 Yu'an Road
Airport Industrial Zone B
Shunyi District, Beijing 101300

Phone +86.10.80426516
Fax +86.10.80426580

Sartorius Stedim Biotech (Beijing) Co. Ltd.

Shanghai Branch Office
3rd Floor, North Wing, Tower 1
No. 4560 Jin ke Road
Pudong District, Shanghai, 201210

Phone +86.21.68782300
Fax +86.21.68782332|68782882

Sartorius Stedim Biotech (Beijing) Co. Ltd.

Guangzhou Representative Office
Unit K, Building 23
Huihua Commerce Et Trade Building
No. 80 Xianlie Middle Road
Room 704, Broadway Plaza,
Guangzhou 510070

Phone +86.20.37618687|37618651
Fax +86.20.37619051

India

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

Phone +91.80.4350.5250
Fax +91.80.4350.5253

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
Fax +81.3.4331.4301

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
Fax +60.3.8996.0755

Singapore

Sartorius Stedim Singapore Pte. Ltd.
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park II
Singapore 117528

Phone +65.6872.3966
Fax +65.6778.2494

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
Fax +82.31.622.5799



◀▶ www.sartorius-stedim.com