



Fill-It

Consumables, Accessories
and Cryovial Compatibility Guide

Simplifying Progress

SARTORIUS

Consumables and Accessories for the Fill-It System

This brochure lists the consumables currently available for the Fill-It system. To order consumables please contact your local office, details on the back cover of this brochure, or email Royston-Orders@sartorius.com

Fill-It Tube Sets

Available in three versions:

- 4-way tube sets compatible with conventional cryovials held in Sartorius supplied 24 vial racks
- V8-way tube sets compatible with cryovials supplied in pre-assembled 48 vial racks
- 8-way tube sets compatible with cryovials supplied in pre-assembled 96 vial racks

Tube set specifications

Construction enables aseptic dispensing within GMP environments:	Supplied sterile and manufactured from animal-derived component free materials meeting USP class VI and FDA 21 CFR 177.2600 classifications
Certification	Supplied with a certificate of conformity
Packaging	Double bagged packaging provides reassurance on product integrity and includes a tear strip for aseptic unpacking
Dispensing nozzle construction	Manufactured from stainless steel to prevent corrosion and angled to reduce likelihood of splashing
Dispensing nozzle bore size (diameter)	0.68 mm
Dispensing nozzle protection	Fitted with a protective cover during transit to avoid damage to tube set and external packaging
Dispensed cell concentrations (typical)	1 to 100 million cells mL
Dead volume (typical after back-flushing)	<0.5 mL
Operating temperature range	+ 4°C to + 30°C
Weight (typical)	69 g
Shelf life	3 years from manufacture

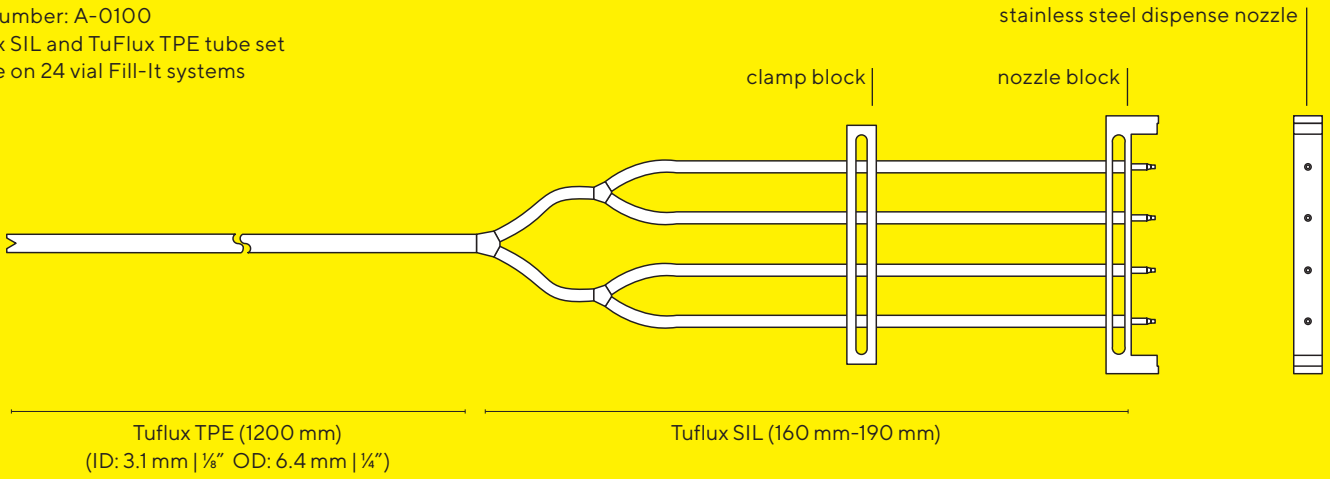


For more information, please visit

www.sartorius-stedim-tap.com/tap/cell_culture/Fill-It.htm

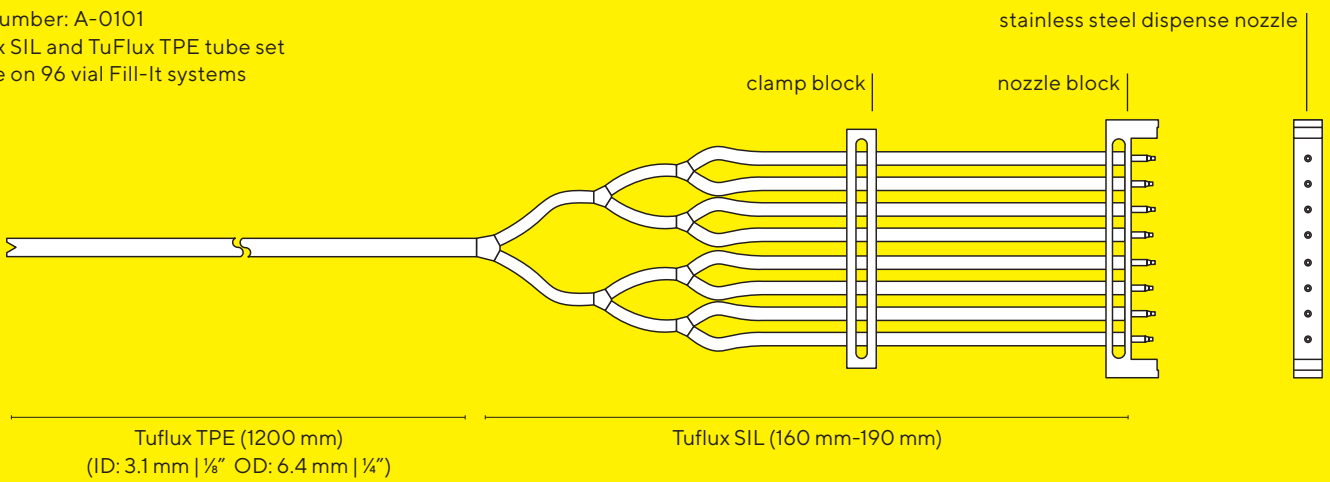
4-way tube set

Part number: A-0100
TuFlux SIL and TuFlux TPE tube set
for use on 24 vial Fill-It systems



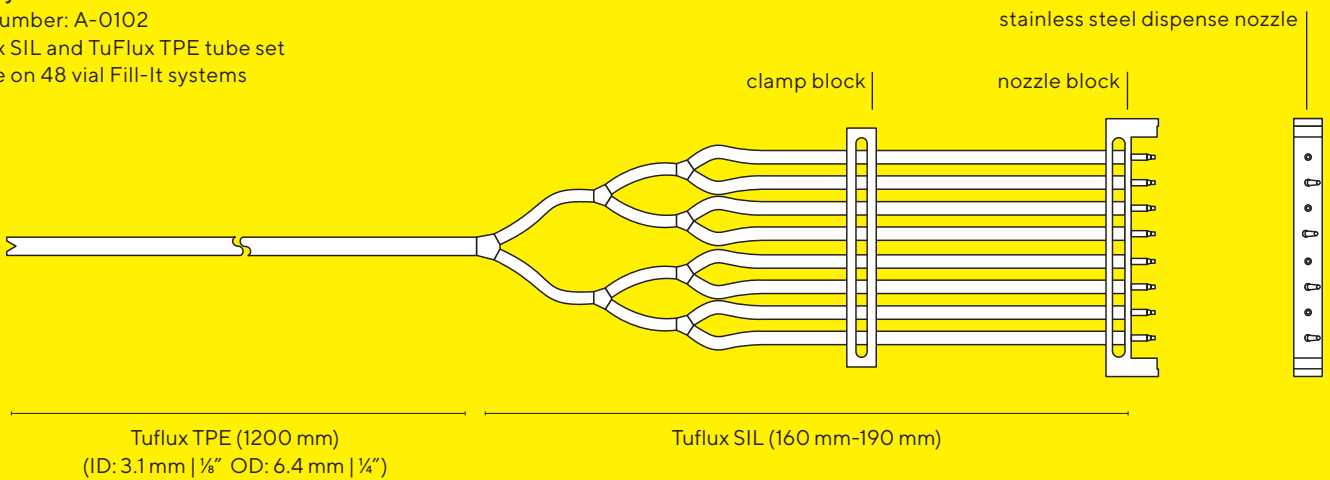
8-way tube set

Part number: A-0101
TuFlux SIL and TuFlux TPE tube set
for use on 96 vial Fill-It systems



V8-way tube set

Part number: A-0102
TuFlux SIL and TuFlux TPE tube set
for use on 48 vial Fill-It systems



Note: TuFlux Tube Sets – due for launch January 2021

Dispensing Pots



Mini-Stacker
Trolley



GMP Priming vessel



Fill-It Priming Tubs

24-way priming tub, stainless steel

Part no. 063-8D14

48-way and 96-way filling system priming tub, stainless steel

Part no. 063-8D15

Fill-It Dispensing Pots

Large dispensing pot, 1.5 L

Part no. 063-8D12

Glass Dispensing pot suitable for GMP laboratories

Small dispensing pot, 650 mL

Glass Dispensing pot suitable for GMP laboratories

Part no. 063-8D11

Handling Devices

Mini-stacker trolley

For safe handling and transport of Fill-It

Part no. 063-8D13



Fill-It Cryovial Racks

Specially designed 24 vial racks are available to hold loose cryovial from specific manufacturers.

O63-8D20 24 vial rack for Nunc and Corning cryovials 1-2 mL

O63-8D16 24 vial rack for Nunc and Corning cryovials 4-5 mL

O63-8D17 24 vial rack for Nalgene cryovials 1 mL

O63-8D18 24 vial rack for Nalgene cryovials 1.5 mL

Fill-It Compatible Cryovial Types

The standard fill-it system is built configured to fit up to 3 sizes of one of the following cryovial types.

Nunc: 96 vial rack system	Manufacturer catalogue number
Internal thread	
0.5 mL	374082, 374083, 374086, 374087
1.0 mL	374085, 374088, 374089

Nunc: 48 vial rack system	
External thread	
1.8 mL (2D coded)	374500, 374500TS, 374500-BR latch rack
1.8 mL (uncoded)	374501 latch rack
Internal thread	
2.0 mL (2D coded)	374221 latch rack
2.0 mL (uncoded)	374223 latch rack
5.0 mL (2D coded)	374220LOCK black lock-in rack
5.0 mL (uncoded)	374222LOCK black lock-in rack

Nunc: Loose cryovials for 24 vial systems	
External thread	
1.8 mL	347627, 375418
4.5 mL	347643, 337516
Internal thread	
1.8 mL	377267
4.5 mL	379146

Matrix: 96 vial rack system	Manufacturer catalogue number
Internal thread	
0.5 mL	3744 (Gre, Yel, Blu, Red, Whi, Purp)
1.0 mL	3741 (Gre, Yel, Blu, Red, Whi, Purp)

Corning: Loose cryovials for 24 vial systems	
External thread	
2.0 mL	430659
4.0 mL	430662
5.0 mL	430663
Internal thread	
1.2 mL	430487
2.0 mL	430488

Nalgene: Loose cryovials for 24 vial systems	
External thread	
1.0 mL	5000-1012
1.5 mL	5000-1020



SARTORIUS


FILL-IT

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