

Clipster® Aseptic Disconnecter



Product Information

The Clipster® Aseptic Disconnecter is a single-use device developed by Sartorius Stedim Biotech that completes our range of products by performing aseptic disconnections of tubing. The Clipster® Aseptic Disconnecter may be sold as a standalone product or preassembled on our Fluid Management bag assemblies.

The Clipster® Aseptic Disconnecter is safe, quick and easy to use. The disconnection is performed with a hand-held tool which ensures easy execution in various space requirements.

Application

The Clipster® Aseptic Disconnecter is used after a fluid transfer to disconnect single-use transfer lines and bag assemblies used in biopharmaceutical applications. The Clipster® Aseptic Disconnecter allows aseptic disconnection in non-classified and classified environments while maintaining product sterility. It can be applied to multiple types and sizes of tubing.

Introduction

Features	Benefits
Mechanical disconnection	Could be performed on platinum-cured silicone and TPE tubings
Error-proof design	Prevents mistakes
Four-step operation	Easy, quick, robust and repeatable
Intensively qualified	Safe and robust
Available as standalone product or preassembled	Flexible
Hand-held tool	Easy to use
Three Clipster® sizes	Compatible with five tubing dimensions

Specifications

Reference	Designation	Material	Tube Size ID-OD*	Tube Material Silicone (Pt) Tuflux® SIL	Tube Material Tuflux® TPE	Tube Material Silicone (Pt) Pharma-50 Si	Tube Material TPE C-Flex®** 374
642XS-----D	Clipster® Aseptic	Polycarbonate	1/8" x 1/4"	■	■	■	■
642-XS-200-N	Disconnecter size XS		1/4" x 3/8"	■	■	■	■
642-S-----D	Clipster® Aseptic	Polycarbonate	1/4" x 7/16"	■	■	■	■
642-S-200-N	Disconnecter size S						
642-L-----D	Clipster® Aseptic	Polycarbonate	3/8" x 5/8"	■	■	■	■
642-L-100-N	Disconnecter size L		1/2" x 3/4"	■	■	■	■

* Note. For operating procedure and conditions, please refer to the instruction for use.

** C-Flex® is a registered trademark of Saint-Gobain Performance Plastics Corporation.

Sterilization Methods

Gamma Irradiation and X-Ray	≤ 50 kGy
Autoclaving	2 × 30 min at 125 °C

Validation

The Clipster® Aseptic Disconnecter has been qualified by applying the most stringent and innovative test regimes including Freeze and Thaw resistance. Biological and physical tests were combined to provide users of the Clipster® Aseptic Disconnecter with data representing a variety of process conditions. Methodologies and equipment are detailed in the Validation Guide.

Quality Assurance

Sartorius Stedim Biotech Quality Systems for single-use products follow applicable ISO and FDA regulations. Design, manufacturing and sterilization processes are conducted under conditions that mirror biopharmaceutical operations and meet cGMP requirements.

Security of Supply

Sartorius Stedim Biotech has established multiple manufacturing sites with consistent industrial processes. The expertise of designing single-use solutions based on collaborative supplier management and customer demand planning ensures a state-of-the-art and robust supply chain that can cope with strong market growth.

Ordering Information

Reference	Designation	Unit Box
642-XS-200-N	Clipster® Large Packaging XS	200
642-S-200-N	Clipster® Large Packaging S	200
642-L-100-N	Clipster® Large Packaging L	100
642-S--T	Hand-held tool for Clipster® XS/S	1
642-L--T	Hand-held tool for Clipster® L	1
642-S--B	Blades for Clipster® XS/S tool	20
642-L--B	Blades for Clipster® L tool	20
642-S-SP	Spare parts box for hand held tool for Clipster® XS S	1 kit
642-L-SP	Spare parts box for hand held tool for Clipster® L	1 kit
642XS-----D	Clipster® Aseptic Disconnecter size XS	10
642S-----D	Clipster® Aseptic Disconnecter size S	10
642L-----D	Clipster® Aseptic Disconnecter size L	10


Note. Make sure you have selected the right Clipster® Aseptic Disconnecter and appropriate Clipster® tool and blades according to your tubing material and dimensions.

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
[sartorius.com](https://www.sartorius.com)

Specifications subject to change without notice.

© 2024 Sartorius Stedim FMT S.A.S., Avenue de Jouques - Zone Industrielle Les Paluds - C.S. 91051, 13781 Aubagne Cedex, France

Status: 05 | 2024