

Instructions for Use | Gebrauchsanleitung | Manuel d'utilisation |
Istruzioni per l'uso | Instrucciones de uso | Gebruiksaanwijzing

Clipster®

Aseptic Disconnecter | Sterile Trennvorrichtung | Déconnecteur aseptique |
Disconnettore aseptico | Desconectador aséptico | Steriele scheidingsinrichting



85032-538-41



SARTORIUS

English	page	3
Deutsch	Seite	17
Français	page	31
Italiano	pagina	45
Español	página	59
Nederlands	Pagina	73

Contents

1	Instruction for Use	4
1.1	Proper Use	4
1.2	Patents	4
2	Introduction	5
3	Operations.....	7
3.1	Clipster® Aseptic Disconnecter Components	7
3.2	Tool Description.....	8
3.3	Warnings.....	9
3.4	Operation Procedure	9
4	Maintenance.....	14
4.1	Procedure to Change the Blade	14
4.2	Cleaning of the Hand Tool	16

1 Instruction for Use

Please read this manual carefully. It includes important information for the safe and successful use of the Clipster® Aseptic Disconnecter.

1.1 Proper Use

Clipster® Aseptic Disconnecter was exclusively developed, constructed and built for the aseptic disconnection of tubing in biopharmaceutical applications.

 The system is exclusively designed for the purpose defined above. Any other application above or beyond this without the written permission of the manufacturer is not considered proper use.

 Sartorius Stedim Biotech cannot assume liability if Clipster® Aseptic Disconnecter is subjected to improper use or improper operation. Sartorius Stedim Biotech is not liable for any damages resulting there from. The operator bears the sole risk.

Proper use also includes compliance with operating, inspection and maintenance conditions prescribed by the manufacturer. Non-compliance with the warnings in the operating instructions constitutes improper use of the device.

1.2 Patents

Clipster® is a trademark of Sartorius Stedim Biotech. The manufacturer reserves the right to make any changes and improvements in the product and process described in this manual at any time without notice. All registered trademarks and patents are the property of their respective owners.

2 Introduction

The Clipster® Aseptic Disconnecter is a single-use device developed by Sartorius Stedim Biotech that completes our range of products by allowing tubing aseptic disconnection. The Clipster® Aseptic Disconnecter may be sold as a stand alone product or preassembled on our Fluid Management bag assemblies. The Clipster® Aseptic Disconnecter is safe, quick and easy to use. The aseptic disconnection is performed with a hand-held tool which ensures easy execution in various space requirements.

The Clipster® Aseptic Disconnecter is used after 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 by maintaining product sterility. It can be applied on multiple types and sizes of tubing.

Specifications

Reference	Designation	Material
642XS-----D	Clipster® Aseptic Disconnecter size XS	Polycarbonate
642-S-----D	Clipster® Aseptic Disconnecter size S	Polycarbonate
642-L-----D	Clipster® Aseptic Disconnecter size L	Polycarbonate

Reference	Tube Size ID – OD*	Tube Material	Tube Material
		Silicone (Pt) Tuflux® SIL	Silicone (Pt) Pharma-50 Si (Pt)
642XS-----D	1/8" × 1/4"	■	■
642XS-----D	1/4" × 3/8"	■	■
642-S-----D	1/4" × 7/16"		■
642-L-----D	3/8" × 5/8"	■	■
642-L-----D	1/2" × 3/4"		■

Reference	Tube Size ID – OD*	Tube Material
		TPE C-Flex®** 374
642XS-----D	1/8" × 1/4"	■
642XS-----D	1/4" × 3/8"	■
642-S-----D	1/4" × 7/16"	■
642-L-----D	3/8" × 5/8"	■
642-L-----D	1/2" × 3/4"	■

Reference	Designation
642-S--T	Hand-held tool for Clipster® XS S
642-S--B	Blades for Clipster® XS S tool
642-L--T	Hand-held tool for Clipster® L
642-L--B	Blades for Clipster® L tool

* Note: 300 mm is the minimum required length of tubing to use the Clipster® Aseptic Disconnecter

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

3 Operations

3.1 Clipster® Aseptic Disconnecter Components

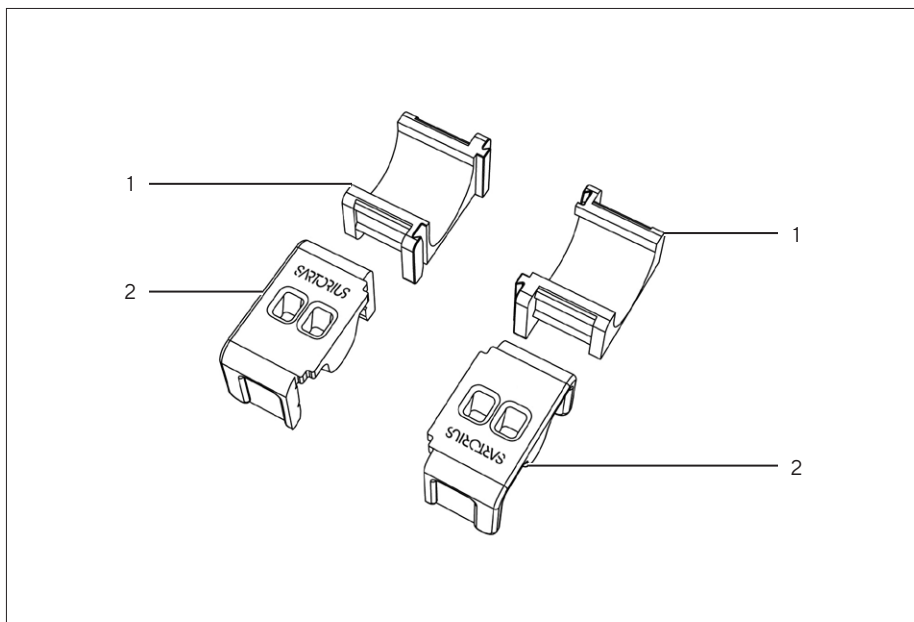


Fig.1

Pos.	Description
1	Clipster® Aseptic Disconnecter bottom component
2	Clipster® Aseptic Disconnecter top component

3.2 Tool Description

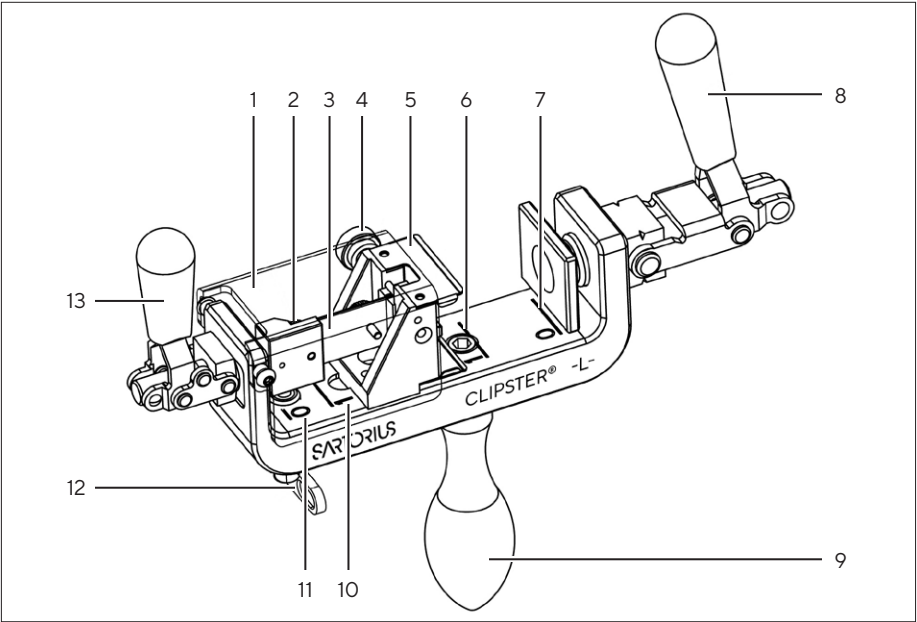


Fig.2

Pos.	Description	Pos.	Description
1	Protective cover	8	Clamping handle
2	Screw Head n° 1	9	Tool handle
3	Blade	10	Position 1 of the cutting handle
4	Screw Head n° 2	11	Position 0 of the cutting handle
5	Clipster® seat	12	Blade Safety Switch
6	Position 1 of the clamping handle	13	Cutting handle
7	Position 0 of the clamping handle		

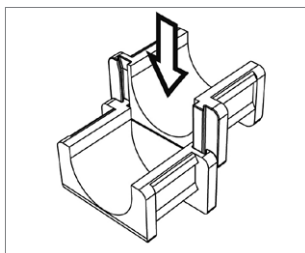
3.3 Warnings

⚠ Operators must be trained before using the Clipster® Aseptic Disconnecter and the handheld tool.

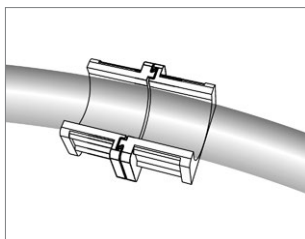
⚠ The Protective cover protects operators from potential injury caused by the blade. Sartorius Stedim Biotech is not liable for any injury if this protective cover is removed during use.

3.4 Operation Procedure

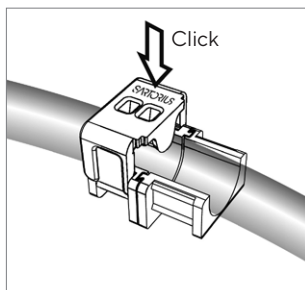
⚠ To guarantee the tubing aseptic disconnection, please follow carefully the operation procedure.



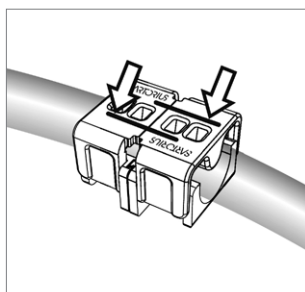
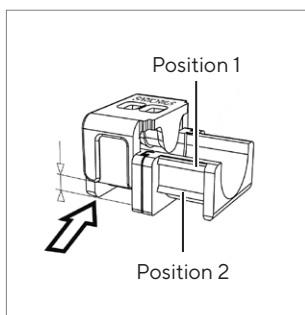
1. Pre-assemble the two bottom components by sliding together along the notched faces.



2. Place the tubing inside the two bottom components.

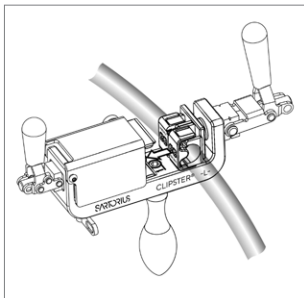


3. Snap one top component in position 1 of the bottom component.

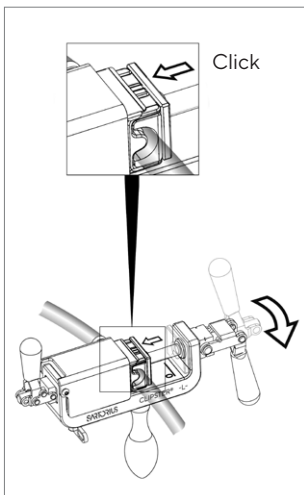
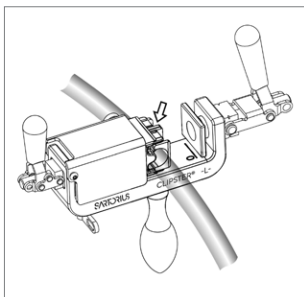


4. Snap the second top component in position 1 of the remaining bottom component and visually check that the Sartorius logo is aligned correctly.

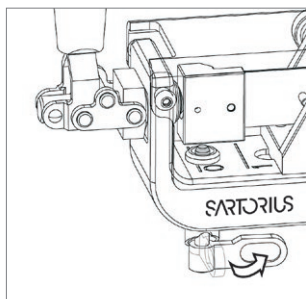
 Before using the hand-held tool, please check that the clamping and cutting handles are in the position 0 and that the security blade is engaged.



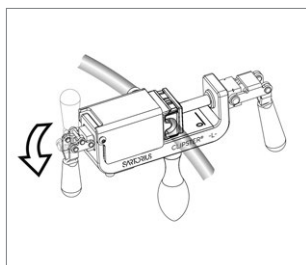
5. Place the assembled Clipster® Aseptic Disconnecter into the Clipster® seat by inserting the top components first (side with the Sartorius logo).



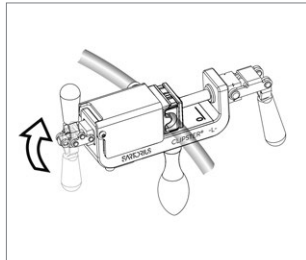
6. Turn the clamping handle to position 1 to clamp the Clipster® Aseptic Disconnecter.



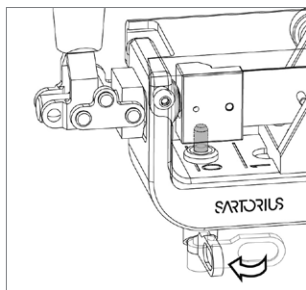
7. Turn the safety switch to release the blade.



8. Keep the clamping handle in position 1 and turn the cutting handle until the position 1 to cut the tubing.

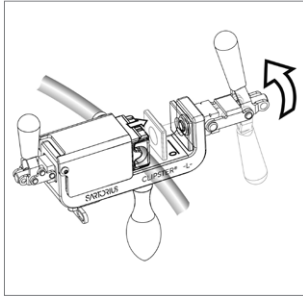


9. Return the cutting handle to the position 0.

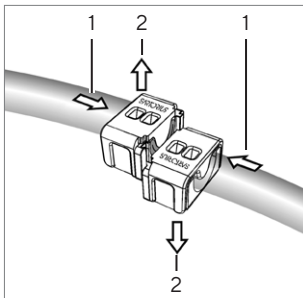


10. Return the blade safety switch to initial position.

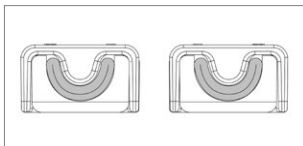
 It is important to return the blade safety switch to initial position for protecting operators from any injury and for using properly the hand-held tool.



11. Return the clamping handle to the position 0 and remove the Clipster® Aseptic Disconnecter from the hand-held tool.



12. Position the logo front of you and separate the two ends by pushing inward and sliding vertically.



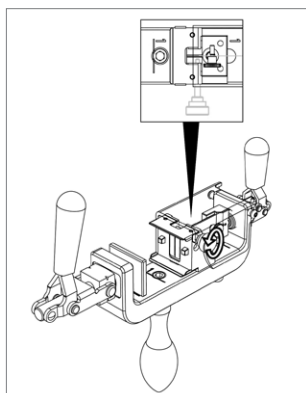
13. View of the two ends after disconnection.

4 Maintenance

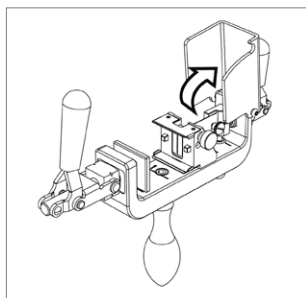
⚠ Operators must be trained before using the Clipster® Aseptic Disconnect and the hand-held tool.

For technical assistance contact your Sartorius Stedim Biotech Service Representative.

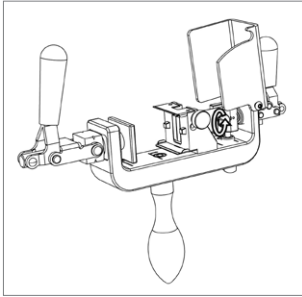
4.1 Procedure to Change the Blade



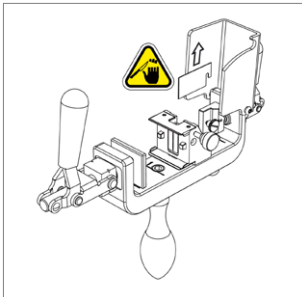
1. Partially unscrew Screw Head n°2 (see figure 2) on the protective cover until the blade is accessible.



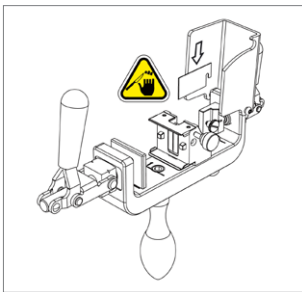
2. Open the protective cover.



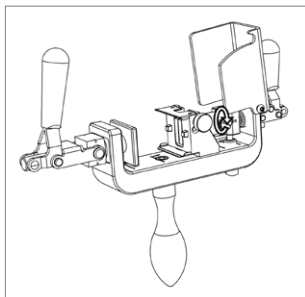
3. Unscrew Screw Head n°1 (see figure 2) on the blade until the blade is released.



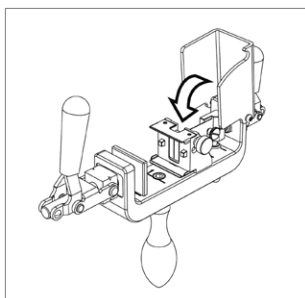
4. Carefully remove the blade.



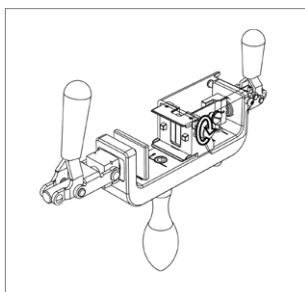
5. Insert a new blade.



6. Tighten Screw Head n°1 to secure the blade.



7. Close the protective cover.



8. Tighten Screw Head n°2 to the protective cover.

4.2 Cleaning of the Hand Tool

The Clipster® Aseptic Disconnector can be cleaned with most disinfection solutions such as isopropyl alcohol.

Inhalt

1	Gebrauchsanweisung	18
1.1	Sachgemäße Verwendung	18
1.2	Patente	18
2	Einleitung	19
3	Gebrauch	21
3.1	Bestandteile der sterilen Clipster®-Trennvorrichtung	21
3.2	Beschreibung der Vorrichtung	22
3.3	Warnhinweise	23
3.4	Arbeitsschritte	23
4	Wartung	28
4.1	Verfahren zum Austausch der Klinge	28
4.2	Reinigen des Handgerätes	30

1 Gebrauchsanweisung

Lesen Sie diese Anweisung sorgfältig durch. Sie enthält wichtige Informationen zur sicheren und erfolgreichen Verwendung der sterilen Clipster®-Trennvorrichtung.

1.1 Sachgemäße Verwendung

Die sterile Clipster®-Trennvorrichtung wurde speziell für das sterile Trennen von Schläuchen in biopharmazeutischen Anwendungen entwickelt, konstruiert und gebaut.

 Die Vorrichtung ist ausschließlich für den oben genannten Zweck vorgesehen. Jede andere, darüber hinausgehende Anwendung wird ohne die schriftliche Genehmigung des Herstellers als nicht sachgemäße Verwendung betrachtet.

 Sartorius Stedim Biotech kann keine Haftung übernehmen, wenn die sterile Clipster®-Trennvorrichtung unsachgemäß eingesetzt oder verwendet wurde. Sartorius Stedim Biotech übernimmt keine Haftung für daraus entstehende Schäden. Der Benutzer trägt das alleinige Risiko.

Die sachgemäße Verwendung umfasst außerdem die Einhaltung der Arbeits-, Inspektions- und Wartungsvorschriften des Herstellers. Die Nichteinhaltung der Warnhinweise in der Betriebsanleitung stellt eine unsachgemäße Verwendung des Gerätes dar.

1.2 Patente

Clipster® ist ein Warenzeichen von Sartorius Stedim Biotech. Der Hersteller behält sich das Recht vor, jederzeit und ohne vorherige Mitteilung Änderungen und Verbesserungen am Produkt und am in dieser Anweisung beschriebenen Verfahren vorzunehmen. Alle eingetragenen Warenzeichen und Patente sind Eigentum der jeweiligen Besitzer.

2 Einleitung

Die sterile Clipster®-Trennvorrichtung ist eine Einweg-Vorrichtung von Sartorius Stedim Biotech, die unsere Produktpalette durch das sterile Trennen von Schläuchen ergänzt. Die sterile Clipster®-Trennvorrichtung kann als Einzelprodukt oder zusammen mit unserer Fluidmanagement-Beuteleinheit verkauft werden. Die sterile Clipster®-Trennvorrichtung ist sicher, schnell und einfach in der Handhabung. Die sterile Trennung erfolgt mit einem Handgerät, das eine einfache Ausführung unter verschiedenen Platzbedingungen gewährleistet.

Die sterile Clipster®-Trennvorrichtung wird nach der Übertragung des Fluids verwendet, um Einweg-Transferleitungen und -Beuteleinheiten in biopharmazeutischen Anwendungen zu trennen. Die sterile Clipster®-Trennvorrichtung ermöglicht eine sterile Trennung in nicht klassifizierten und klassifizierten Umgebungen und gewährleistet Produktsterilität. Sie kann für verschiedene Arten und Größen von Schläuchen verwendet werden.

Technische Daten

Referenz	Bezeichnung	Material
642XS-----D	Clipster® Sterile Trennvorrichtung Größe XS	Polykarbonat
642-S-----D	Clipster® Sterile Trennvorrichtung Größe S	Polykarbonat
642-L-----D	Clipster® Sterile Trennvorrichtung Größe L	Polykarbonat

Referenz	Schlauchgröße ID – OD*	Schlauchmaterial	Schlauchmaterial
		Silikon (Pt) Tuflux® SIL	Silikon (Pt) Pharma-50 Si (Pt)
642XS-----D	1/8" × 1/4"	■	■
642XS-----D	1/4" × 3/8"	■	■
642-S-----D	1/4" × 7/16"		■
642-L-----D	3/8" × 5/8"	■	■
642-L-----D	1/2" × 3/4"		■

Referenz	Schlauchgröße ID – OD*	Schlauchmaterial
		TPE C-Flex®** 374
642XS-----D	1/8" × 1/4"	■
642XS-----D	1/4" × 3/8"	■
642-S-----D	1/4" × 7/16"	■
642-L-----D	3/8" × 5/8"	■
642-L-----D	1/2" × 3/4"	■

Referenz	Bezeichnung
642-S--T	Handwerkzeuggerät für Clipster® XS S
642-S--B	Klingen für Clipster® XS S Werkzeug
642-L--T	Handwerkzeuggerät für Clipster® L
642-L--B	Klingen für Clipster® L Werkzeug

* Hinweis: Für die sterile Clipster®-Trennvorrichtung sind Schläuche mit einer Mindestlänge von 300 mm notwendig.

** C-Flex® ist eine eingetragene Marke der Saint-Gobain Performance Plastics Corporation.

3 Gebrauch

3.1 Bestandteile der sterilen Clipster®-Trennvorrichtung

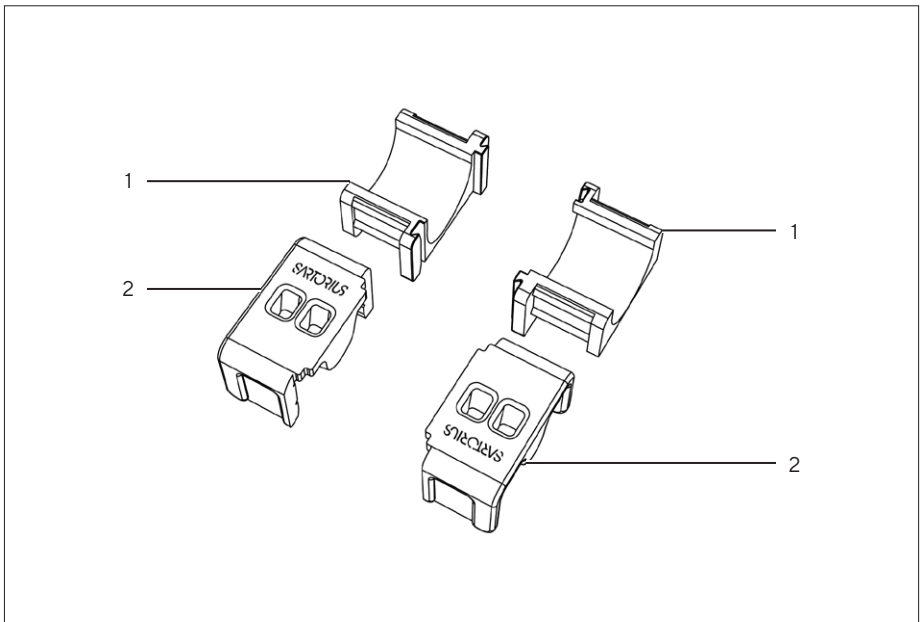


Abb.1

Pos.	Beschreibung
1	Unterteil der sterilen Clipster®-Trennvorrichtung
2	Oberteil der sterilen Clipster®-Trennvorrichtung

3.2 Beschreibung der Vorrichtung

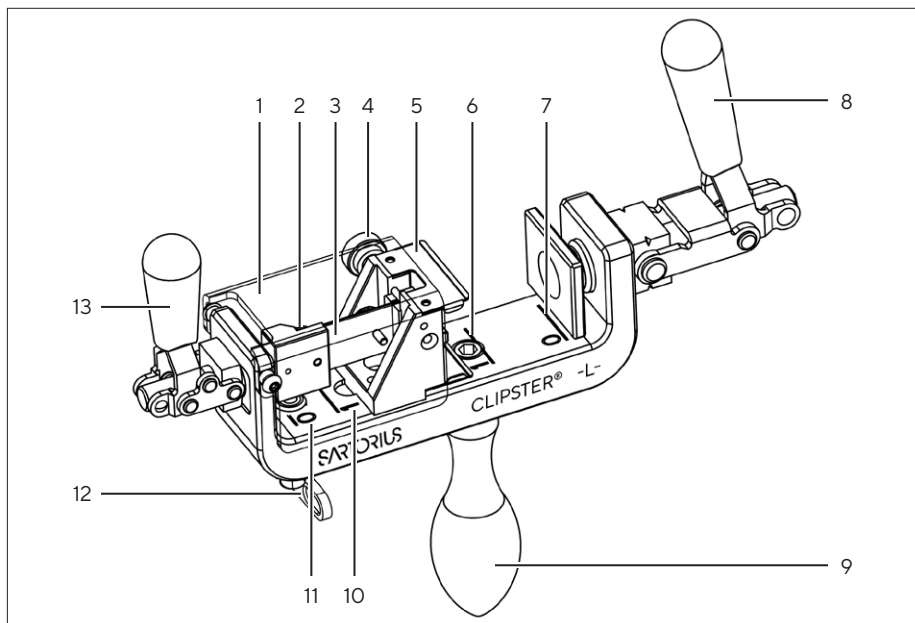


Abb.2

Pos.	Beschreibung
1	Schutzabdeckung
2	Schraubenkopf Nr. 1
3	Klinge
4	Schraubenkopf Nr. 2
5	Clipster®-Sitz
6	Position 1 des Klemmgriffes
7	Position 0 des Klemmgriffes

Pos.	Beschreibung
8	Klemmgriff
9	Griff der Vorrichtung
10	Position 1 des Schneidegriffes
11	Position 0 des Schneidegriffes
12	Klingensicherheitsschalter
13	Schneidegriff

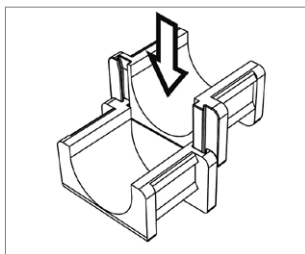
3.3 Warnhinweise

⚠ Die Benutzer müssen vor der Verwendung der sterilen Clipster®-Trennvorrichtung und des Handgerätes geschult werden.

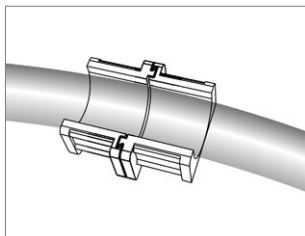
⚠ Die Schutzabdeckung schützt den Benutzer vor möglichen Verletzungen durch die Klinge. Sartorius Stedim Biotech haftet nicht für Verletzungen, wenn diese Schutzabdeckung während der Verwendung abgenommen wurde.

3.4 Arbeitsschritte

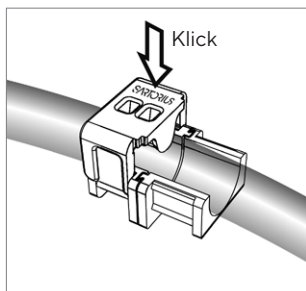
⚠ Eine sterile Trennung von Schläuchen ist nur möglich, wenn folgende Schritte sorgfältig eingehalten werden.



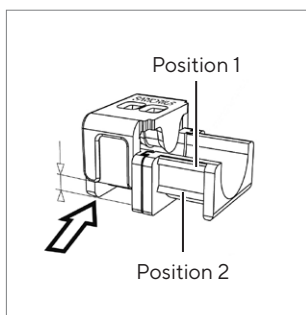
1. Die beiden Unterteile an den eingekerbten Flächen zusammenschieben.



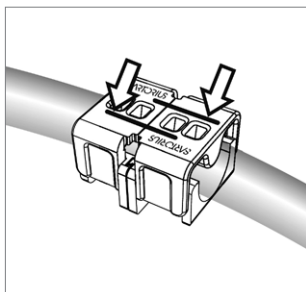
2. Den Schlauch in die beiden Unterteile legen.



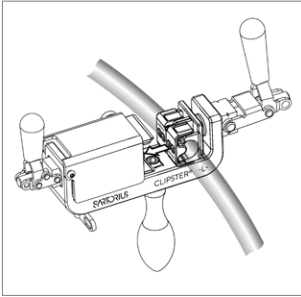
3. Ein Oberteil in Position 1 des Unterteils einrasten.



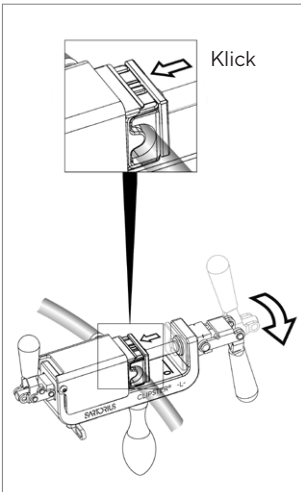
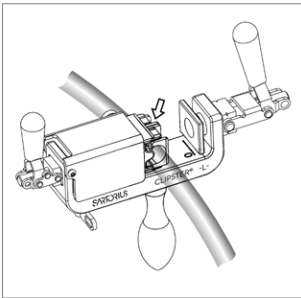
4. Das zweite Oberteil in Position 1 des anderen Unterteils einrasten und visuell überprüfen, ob das Sartorius Logo richtig ausgerichtet ist.



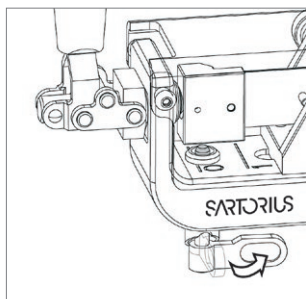
⚠ Vor der Verwendung des Handgeräts müssen sich die Klemm- und Schneidegriffe in Position 0 befinden und die Sicherheitsklinge muss eingearastet sein.



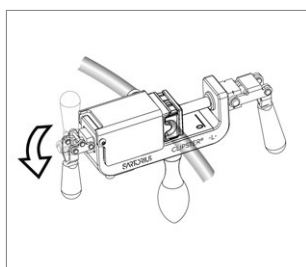
5. Die montierte sterile Clipster®- Trenvorrichtung mit den Oberteilen (Seite mit Sartorius Logo) zuerst in den Clipster®- Sitz einsetzen.



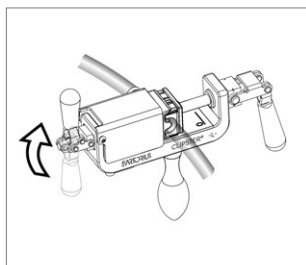
6. Den Klemmgriff in Position 1 drehen, um die sterile Clipster®-Trenvorrichtung festzuklemmen.



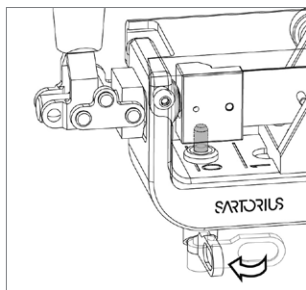
7. Den Sicherheitsschalter drehen, um die Klinge freizugeben.



8. Den Klemmgriff in Position 1 halten und den Schneidegriff bis zur Position 1 drehen, um den Schlauch durchzutrennen.

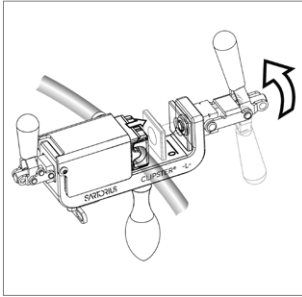


9. Den Schneidegriff wieder in Position 0 zurückbringen.

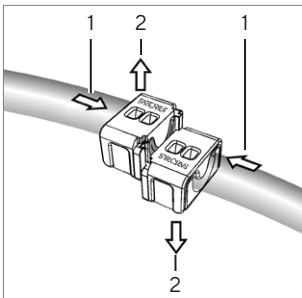


10. Den Klingensicherheitsschalter in die Ausgangsposition stellen.

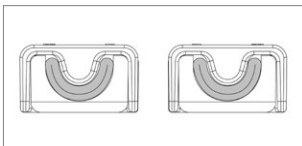
⚠ Es ist wichtig, den Klingensicherheitsschalter in die Ausgangsposition zurückzubringen, um die Benutzer vor Verletzungen zu schützen und das Handgerät sachgemäß zu verwenden.



11. Den Klemmgriff in die Position 0 zurückstellen und die sterile Clipster®-Trennvorrichtung aus dem Handgerät herausnehmen.



12. Das Logo in Ihre Richtung drehen und die beiden Enden auseinanderziehen, indem Sie nach innen drücken und nach oben schieben.



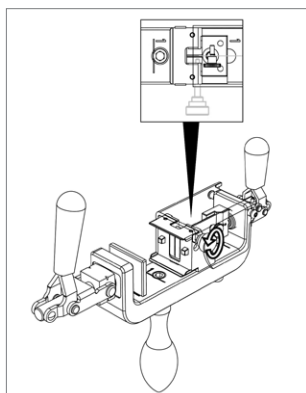
13. Die beiden Enden nach dem Trennen überprüfen.

4 Wartung

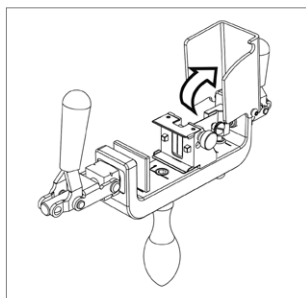
⚠ Die Benutzer müssen vor der Verwendung der sterilen Clipster®-Trennvorrichtung und des Handgerätes geschult werden.

Technische Unterstützung erhalten Sie von Ihrem Ansprechpartner bei Sartorius Stedim Biotech.

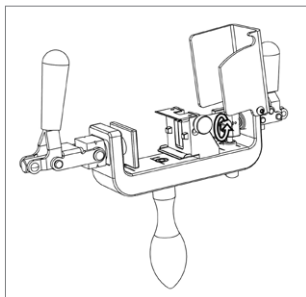
4.1 Verfahren zum Austausch der Klinge



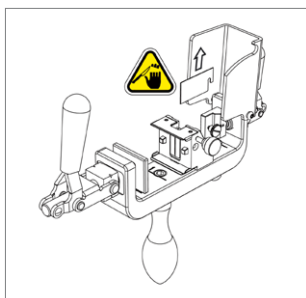
1. Den Schraubenkopf Nr. 2 (siehe Abbildung 2) auf der Schutzabdeckung etwas lösen, bis die Klinge freigegeben wird.



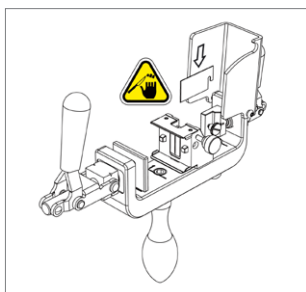
2. Die Schutzabdeckung öffnen.



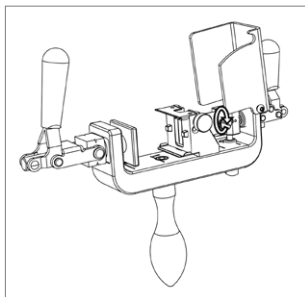
3. Den Schraubenkopf Nr. 1 (siehe Abbildung 2) auf der Klinge etwas lösen, bis die Klinge freigegeben wird.



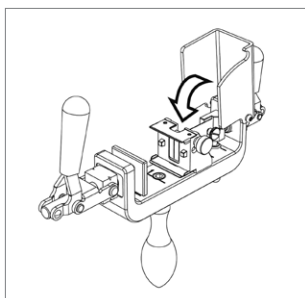
4. Die Klinge vorsichtig herausnehmen.



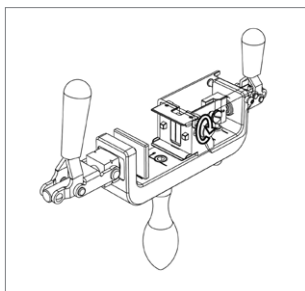
5. Eine neue Klinge einsetzen.



6. Den Schraubenkopf Nr. 1 festziehen, um die Klinge zu sichern.



7. Die Schutzabdeckung schließen.



8. Den Schraubenkopf Nr. 2 auf der Schutzabdeckung festziehen.

4.2 Reinigen des Handgerätes

Die sterile Clipster®-Trennvorrichtung kann mit den meisten Desinfektionslösungen wie Isopropylalkohol gereinigt werden.

Table des matières

1	Conseils d'utilisation	32
1.1	Utilisation conforme	32
1.2	Brevets	32
2	Introduction	33
3	Fonctionnement	35
3.1	Composants du déconnecteur aseptique Clipster®	35
3.2	Description de l'outil	36
3.3	Avertissements	37
3.4	Procédure de fonctionnement	37
4	Maintenance	42
4.1	Remplacement de la lame	42
4.2	Nettoyage de l'outil à main	44

1 Conseils d'utilisation

Veuillez lire attentivement ce manuel. Il contient des informations importantes qui permettent une utilisation sûre et correcte du déconnecteur aseptique Clipster®.

1.1 Utilisation conforme

Le déconnecteur aseptique Clipster® a été développé, conçu et construit exclusivement pour la déconnexion aseptique de tuyaux dans des applications biopharmaceutiques.

 Le système est destiné exclusivement à l'usage défini ci-dessus. Toute utilisation différente de celle définie ci-dessus et n'ayant pas reçu l'autorisation écrite du fabricant est considérée comme non conforme.

 Sartorius Stedim Biotech ne peut pas être considéré comme responsable si le déconnecteur aseptique Clipster® est soumis à une utilisation ou à fonctionnement impropres. Sartorius Stedim Biotech n'est pas tenu responsable des dommages pouvant en résulter. Seul l'utilisateur en assume les risques.

Pour utiliser l'appareil de manière conforme, il faut également respecter les conditions de fonctionnement, d'inspection et de maintenance prévues par le fabricant. La non-observation des avertissements contenus dans le mode d'emploi est considérée comme une utilisation impropre du dispositif.

1.2 Brevets

Clipster® est une marque déposée de Sartorius Stedim Biotech. Le fabricant se réserve le droit de modifier et d'améliorer à tout moment et sans préavis le produit et le processus décrits dans ce manuel. Toutes les marques et brevets déposés sont la propriété de leurs propriétaires respectifs.

2 Introduction

Le déconnecteur aseptique Clipster® est un dispositif à usage unique développé par Sartorius Stedim Biotech qui complète notre gamme de produits en permettant la déconnexion aseptique de tuyaux. Le déconnecteur aseptique Clipster® peut être vendu séparément ou préassemblé sur nos systèmes de poches à usage unique. Le déconnecteur aseptique Clipster® est sûr, rapide et facile d'utilisation. La déconnexion aseptique s'effectue grâce à un outil qui permet une utilisation sûre et facile quelque soit l'encombrement spacial.

Le déconnecteur aseptique Clipster® est utilisé après un transfert de fluides pour déconnecter des lignes de transfert à usage unique ou des systèmes de poches utilisés dans des applications biopharmaceutiques. Le déconnecteur aseptique Clipster® permet la déconnexion aseptique dans des environnements non classifiés et classifiés en maintenant la stérilité du produit. Il peut être utilisé avec des tubes de différentes dimensions.

Caractéristiques techniques

Référence	Désignation	Matière
642XS-----D	Déconnecteur aseptique Clipster® dimensions XS	Polycarbonate
642-S-----D	Déconnecteur aseptique Clipster® dimensions S	Polycarbonate
642-L-----D	Déconnecteur aseptique Clipster® dimensions L	Polycarbonate

Référence	Dimensions du tube ID – OD*	Matière du tube Silicone (Pt) TuFlux® SIL	Matière du tube Silicone (Pt) Pharma-50 Si (Pt)
642XS-----D	1/8" × 1/4"	■	■
642XS-----D	1/4" × 3/8"	■	■
642-S-----D	1/4" × 7/16"		■
642-L-----D	3/8" × 5/8"	■	■
642-L-----D	1/2" × 3/4"		■

Référence	Dimensions du tube ID – OD*	Matière du tube TPE C-Flex®** 374
642XS-----D	1/8" × 1/4"	■
642XS-----D	1/4" × 3/8"	■
642-S-----D	1/4" × 7/16"	■
642-L-----D	3/8" × 5/8"	■
642-L-----D	1/2" × 3/4"	■

Référence	Désignation
642-S--T	Outil à main pour Clipster® XS S
642-S--B	Lames pour Clipster® XS S outil
642-L--T	Outil à main pour Clipster® L
642-L--B	Lames pour Clipster® L outil

* Remarque : pour utiliser le déconnecteur aseptique Clipster®, les tubes doivent avoir une longueur minimum de 300 mm.

** C-Flex® est une marque déposée de Saint-Gobain Performance Plastics Corporation.

3 Fonctionnement

3.1 Composants du déconnecteur aseptique Clipster®

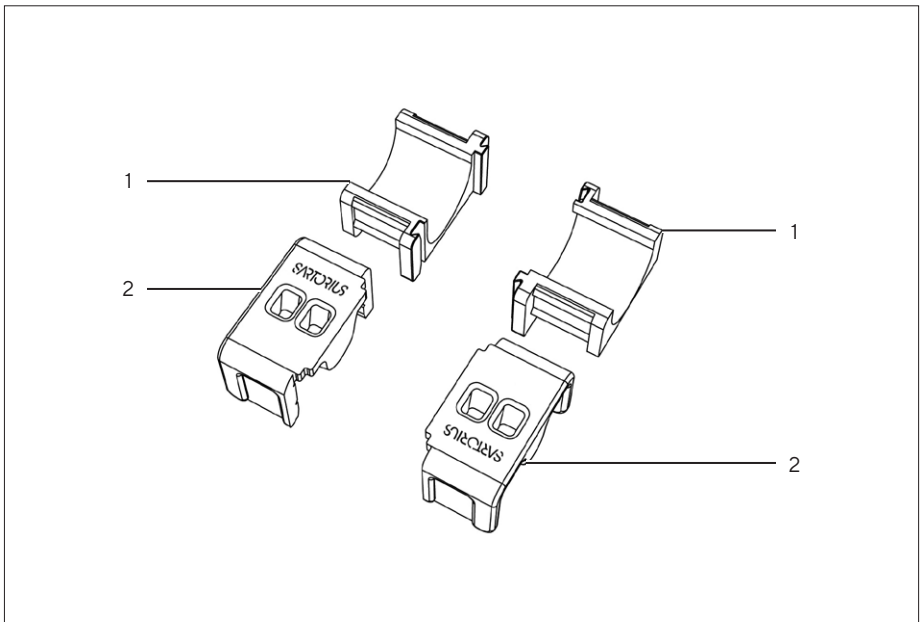


Fig.1

Pos.	Description
1	Composant inférieur du déconnecteur aseptique Clipster®
2	Composant supérieur du déconnecteur aseptique Clipster®

3.2 Description de l'outil

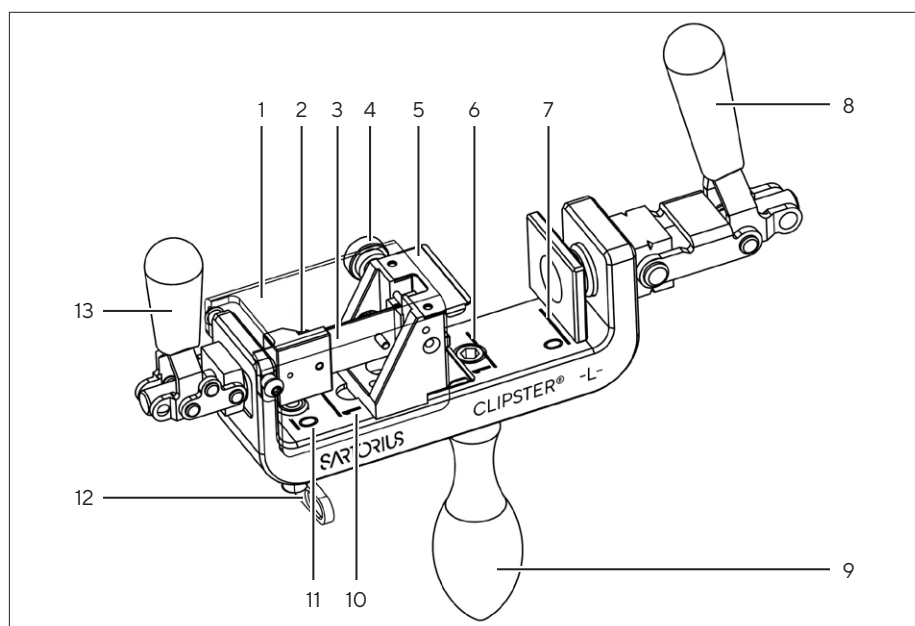


Fig.2

Pos.	Description
1	Cache de protection
2	Tête de vis n° 1
3	Lame
4	Tête de vis n° 2
5	Emplacement du Clipster®
6	Position 1 de la poignée de serrage
7	Position 0 de la poignée de serrage

Pos.	Description
8	Poignée de serrage
9	Poignée de l'outil
10	Position 1 de la poignée de coupe
11	Position 0 de la poignée de coupe
12	Dispositif de sécurité de la lame
13	Poignée de coupe

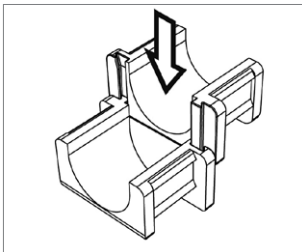
3.3 Avertissements

⚠ Les opérateurs doivent être formés avant d'utiliser le déconnecteur aseptique Clipster® et son outil associé.

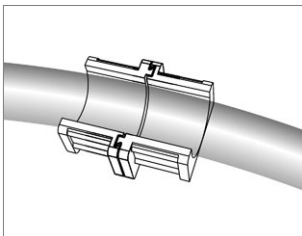
⚠ Le cache de protection protège les opérateurs de blessures potentielles causées par la lame. Sartorius Stedim Biotech décline toute responsabilité en cas de blessures si le cache de protection est enlevé pendant l'utilisation du dispositif.

3.4 Procédure de fonctionnement

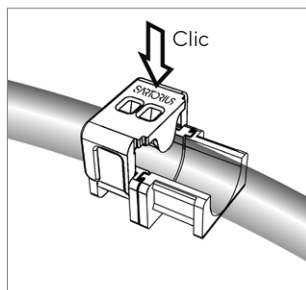
⚠ Pour garantir la déconnexion aseptique des tubes, veuillez suivre scrupuleusement la procédure de fonctionnement.



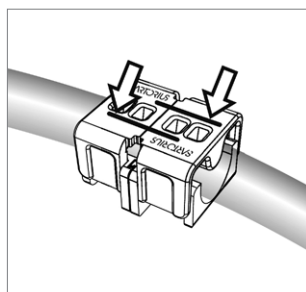
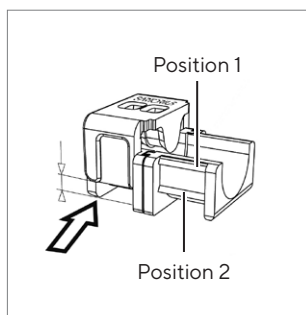
1. Préassemblez les deux composants inférieurs en les faisant coulisser l'un dans l'autre le long des rainures.



2. Mettez le tube à l'intérieur des deux composants inférieurs.

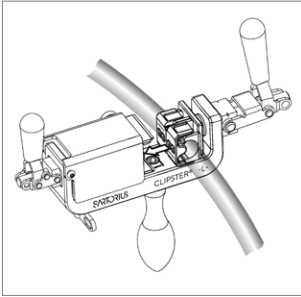


3. Assemblez l'un des composants supérieurs dans la position 1 du composant inférieur.

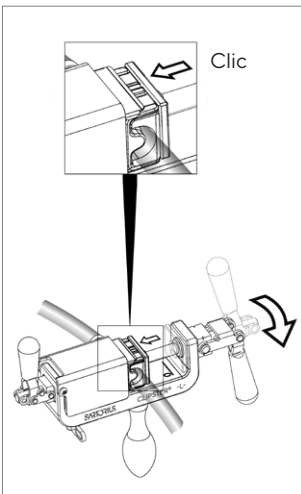
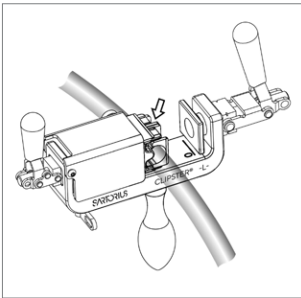


4. Assemblez le deuxième composant supérieur dans la position 1 du composant inférieur restant et vérifiez que le logo Sartorius est correctement écrit.

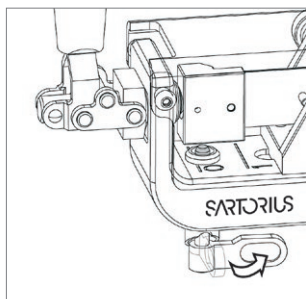
⚠ Avant d'utiliser l'outil à main, vérifiez que les poignées de serrage et de coupe sont dans la position 0 et que la lame de sécurité est engagée.



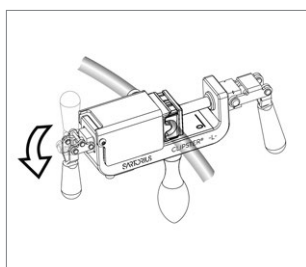
5. Mettez le déconnecteur aseptique Clipster® assemblé dans l'emplacement prévu en insérant d'abord les composants supérieurs (côté avec le logo Sartorius).



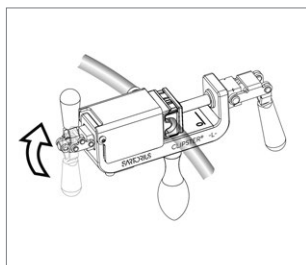
6. Tournez la poignée de serrage dans la position 1 pour clamper | fermer le déconnecteur aseptique Clipster®.



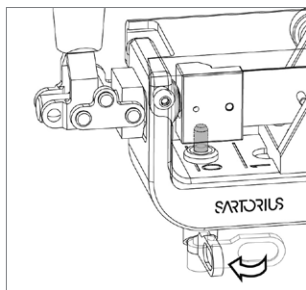
7. Tournez le dispositif de sécurité pour libérer la lame.



8. Maintenez la poignée de serrage dans la position 1 et tournez la poignée de coupe jusqu'à la position 1 pour couper le tube.

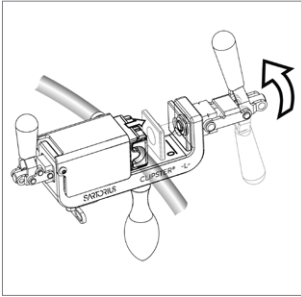


9. Remettez la poignée de coupe dans la position 0.

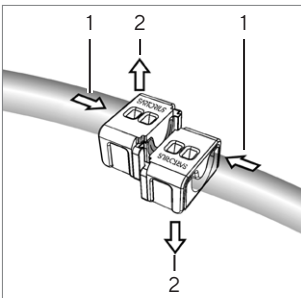


10. Remettez le dispositif de sécurité de la lame dans sa position initiale.

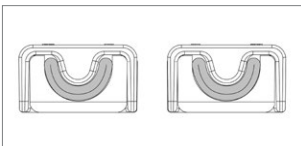
⚠ Il est important de remettre le dispositif de sécurité de la lame dans sa position initiale pour protéger les opérateurs de tout risque de blessure et pour utiliser correctement l'outil.



11. Remettez la poignée de serrage dans la position 0 et enlevez le déconnecteur aseptique Clipster® de l'outil à main.



12. Placez le logo en face de vous et séparez les deux extrémités en les poussant vers l'intérieur tout en les couissant verticalement.



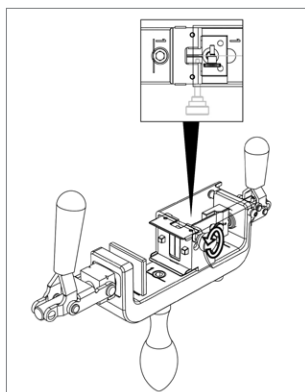
13. Vue des deux extrémités après la déconnexion.

4 Maintenance

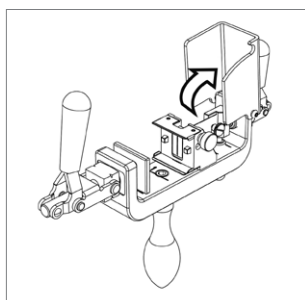
⚠ Les opérateurs doivent avoir été formés avant d'utiliser le déconnecteur aseptique Clipster® et l'outil à main.

Pour obtenir de l'aide, contactez le service après-vente de Sartorius Stedim Biotech.

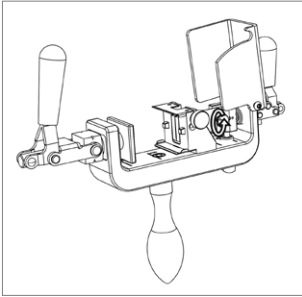
4.1 Remplacement de la lame



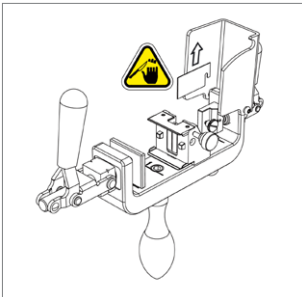
1. Dévissez en partie la tête de vis n°2 (voir figure 2) sur le cache de protection jusqu'à ce que la lame soit libérée.



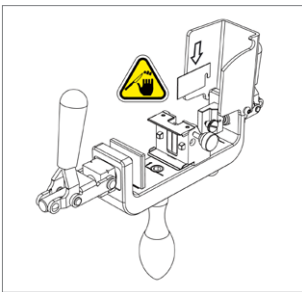
2. Ouvrez le cache de protection.



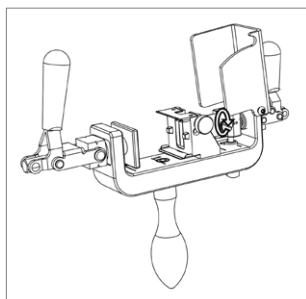
3. Dévissez en partie la tête de vis n°1 (voir figure 2) sur la lame jusqu'à ce que la lame soit libérée.



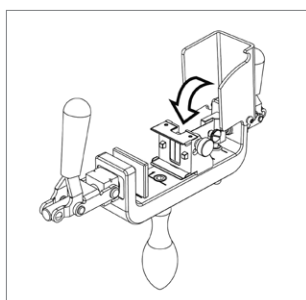
4. Enlevez la lame avec précaution.



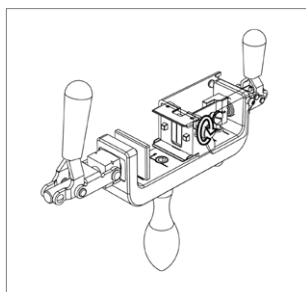
5. Insérez une nouvelle lame.



6. Vissez la tête de vis n°1 pour fixer la lame.



7. Fermez le cache de protection.



8. Vissez la tête de vis n°2 sur le cache de protection.

4.2 Nettoyage de l'outil à main

Le déconnecteur aseptique Clipster® peut être nettoyé avec la plupart des solutions désinfectantes telles que l'alcool isopropylique.

Indice

1 Istruzioni per l'uso.....	46
1.1 Uso previsto	46
1.2 Brevetti	46
2 Introduzione	47
3 Funzionamento	49
3.1 Componenti del disconnettore asettico Clipster®	49
3.2 Descrizione dell'utensile	50
3.3 Avvertenze	51
3.4 Procedura di funzionamento.....	51
4 Manutenzione	56
4.1 Procedura per la sostituzione della lama	56
4.2 Pulizia dell'utensile a mano	58

1 Istruzioni per l'uso

Si prega di leggere attentamente questo manuale di istruzioni. Contiene informazioni importanti per un utilizzo sicuro e corretto del disconnettore aseptico Clipster®.

1.1 Uso previsto

Il disconnettore aseptico Clipster® è stato sviluppato, progettato e costruito esclusivamente per lo scollegamento in aseptico dei tubi nelle applicazioni biofarmaceutiche.

 Il sistema è concepito esclusivamente per lo scopo definito sopra. Qualsiasi altra applicazione diversa da quella sopra menzionata, senza autorizzazione scritta da parte del costruttore, è da considerarsi non conforme all'uso previsto.

 Sartorius Stedim Biotech non si assume alcuna responsabilità relativa a uso o funzionamento improprio del disconnettore aseptico Clipster®. Sartorius Stedim Biotech declina ogni responsabilità per eventuali danni da ciò derivanti. L'utilizzo avviene a rischio esclusivo dell'operatore.

Per uso previsto s'intende anche il rispetto delle condizioni di funzionamento, ispezione e manutenzione previste dal costruttore. La non osservanza delle avvertenze contenute nelle istruzioni per l'uso costituisce un uso improprio del dispositivo.

1.2 Brevetti

Clipster® è un marchio registrato di Sartorius Stedim Biotech. Il costruttore si riserva il diritto di apportare modifiche o miglioramenti al prodotto e al processo descritto nel presente manuale in ogni momento e senza preavviso alcuno. Tutti i marchi registrati e i brevetti sono di proprietà dei legittimi proprietari.

2 Introduzione

Il disconnettore asettico Clipster® è un dispositivo monouso, sviluppato da Sartorius Stedim Biotech a completamento della gamma di prodotti, che permette lo scollegamento in asettico dei tubi. Il disconnettore asettico Clipster® può essere venduto come prodotto stand-alone oppure preassemblato sui sistemi di sacche per il trattamento dei fluidi. Il disconnettore asettico Clipster® è sicuro, rapido e facile da usare. Lo scollegamento in asettico viene eseguito con un utensile a mano che permette una facile esecuzione a seconda delle diverse esigenze di spazio.

Il disconnettore asettico Clipster® viene usato dopo il trasferimento del fluido per scollegare le linee di trasferimento monouso e i sistemi di sacche usati nelle applicazioni biofarmaceutiche. Il disconnettore asettico Clipster® consente lo scollegamento in asettico in ambienti classificati e non classificati mantenendo la sterilità del prodotto. Questo dispositivo può essere usato per tubi di svariati modelli e dimensioni.

Specifiche tecniche

Codice	Denominazione	Materiale
642XS-----D	Disconnettore asettico Clipster® dimensione XS	Policarbonato
642-S-----D	Disconnettore asettico Clipster® dimensione S	Policarbonato
642-L-----D	Disconnettore asettico Clipster® dimensione L	Policarbonato

Codice	Dimensioni del tubo ID – OD*	Materiale del tubo	Materiale del tubo
		Silicone (Pt) Tuflux® SIL	Silicone (Pt) Pharma-50 Si (Pt)
642XS-----D	$\frac{1}{8}" \times \frac{1}{4}"$	■	■
642XS-----D	$\frac{1}{4}" \times \frac{3}{8}"$	■	■
642-S-----D	$\frac{1}{4}" \times \frac{7}{16}"$		■
642-L-----D	$\frac{3}{8}" \times \frac{5}{8}"$	■	■
642-L-----D	$\frac{1}{2}" \times \frac{3}{4}"$		■

Codice	Dimensions du tube ID – OD*	Materiale del tubo
		TPE C-Flex®** 374
642XS-----D	$\frac{1}{8}" \times \frac{1}{4}"$	■
642XS-----D	$\frac{1}{4}" \times \frac{3}{8}"$	■
642-S-----D	$\frac{1}{4}" \times \frac{7}{16}"$	■
642-L-----D	$\frac{3}{8}" \times \frac{5}{8}"$	■
642-L-----D	$\frac{1}{2}" \times \frac{3}{4}"$	■

Codice	Denominazione
642-S--T	Strumento tenuto in mano per Clipster® XS S
642-S--B	Lame per Clipster® XS S strumento
642-L--T	Strumento tenuto in mano per Clipster® L
642-L--B	Lame per Clipster® L strumento

* Nota: 300 mm è la lunghezza minima del tubo richiesta per l'uso del disconnettore aseptico Clipster®

** C-Flex® è un marchio registrato di Saint-Gobain Performance Plastics Corporation

3 Funzionamento

3.1 Componenti del disconnettore asettico Clipster®

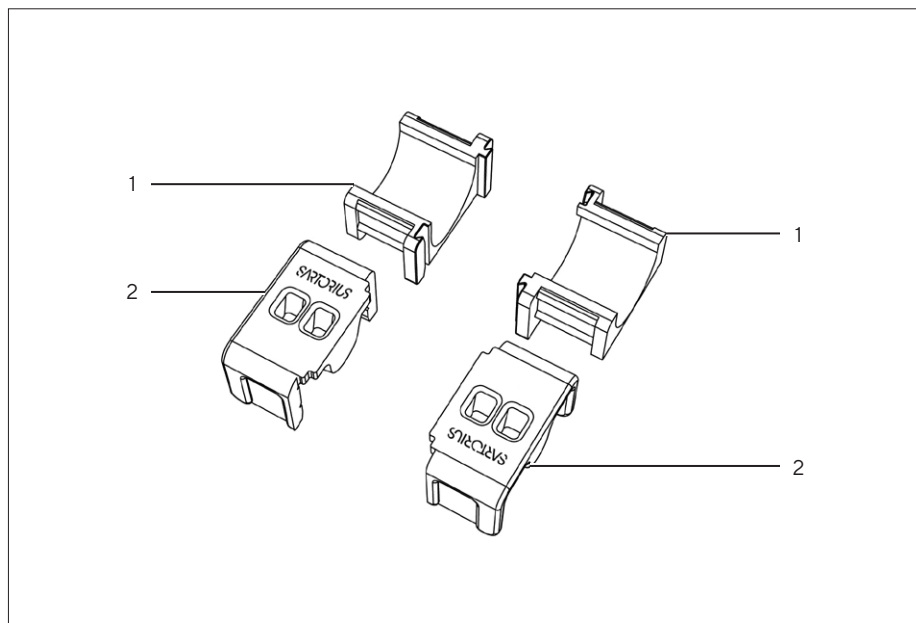


Fig.1

Pos.	Descrizione
1	Componente inferiore del disconnettore asettico Clipster®
2	Componente superiore del disconnettore asettico Clipster®

3.2 Descrizione dell'utensile

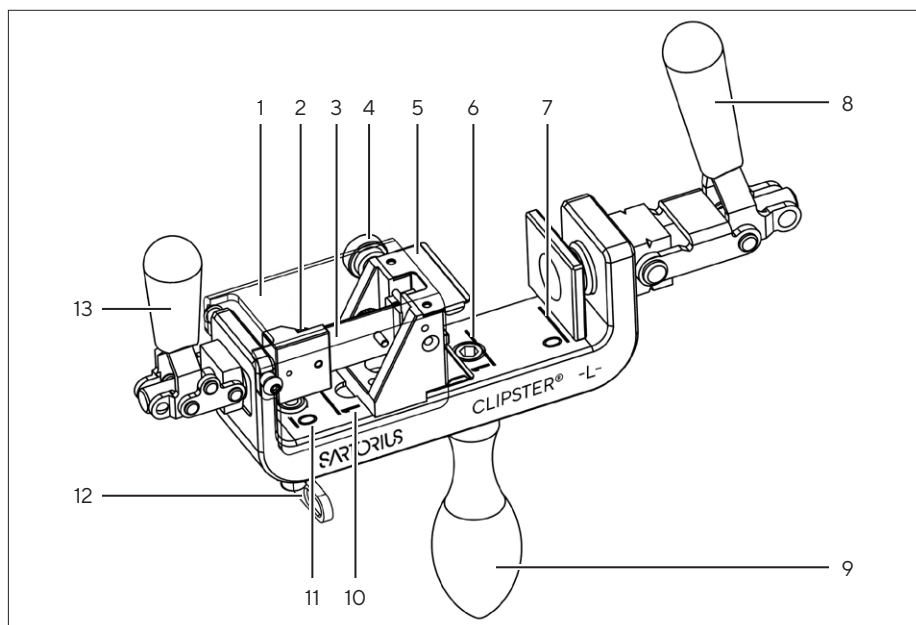


Fig.2

Pos.	Descrizione
1	Copertura di protezione
2	Testa della vite n° 1
3	Lama
4	Testa della vite n° 2
5	Alloggiamento del Clipster®
6	Posizione 1 della maniglia di serraggio
7	Posizione 0 della maniglia di serraggio

Pos.	Descrizione
8	Maniglia di serraggio
9	Manico dell'utensile
10	Posizione 1 della maniglia di taglio
11	Posizione 0 della maniglia di taglio
12	Dispositivo di sicurezza della lama
13	Maniglia di taglio

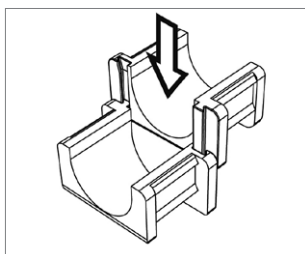
3.3 Avvertenze

⚠ Gli operatori devono essere addestrati prima di usare il disconnettore asettico Clipster® e l'utensile a mano.

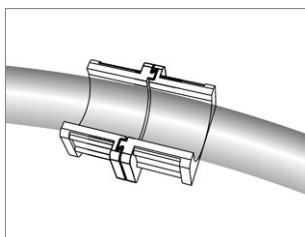
⚠ La copertura protegge gli operatori da potenziali lesioni derivanti dalla lama. Sartorius Stedim Biotech declina ogni responsabilità per eventuali lesioni se durante l'uso è stata tolta la copertura di protezione.

3.4 Procedura di funzionamento

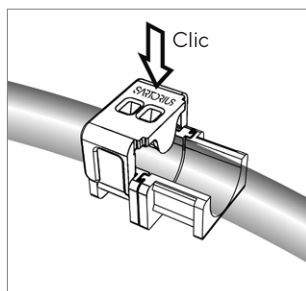
⚠ Per garantire lo scollegamento in asettico del tubo, seguire attentamente la procedura di funzionamento.



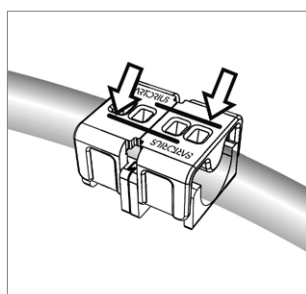
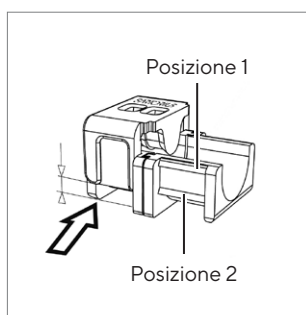
1. Preassemblare i due componenti inferiori facendoli scivolare insieme lungo i lati incavati.



2. Posizionare il tubo all'interno dei due componenti inferiori.

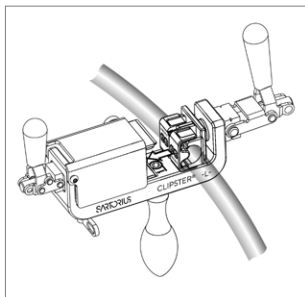


3. Far scattare uno dei componenti superiori nella posizione 1 del componente inferiore.

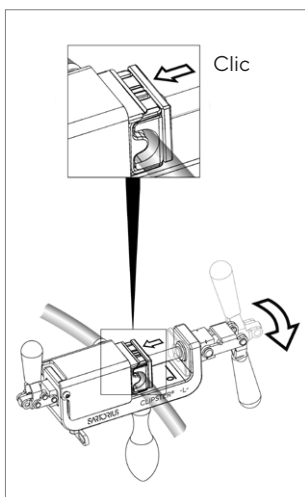
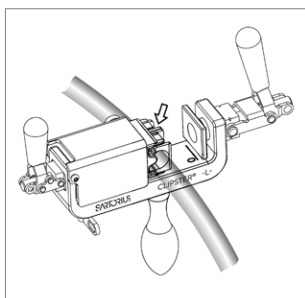


4. Far scattare il secondo componente superiore nella posizione 1 del componente inferiore restante e controllare che il logo Sartorius sia allineato correttamente.

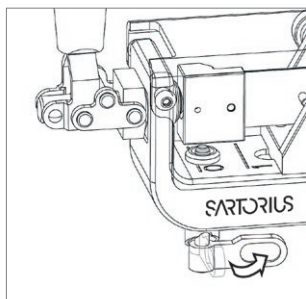
⚠ Prima di usare l'utensile a mano, controllare che la maniglia di serraggio e quella di taglio siano nella posizione 0 e che la lama di sicurezza sia innestata.



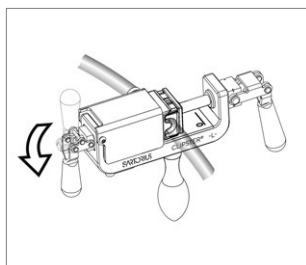
5. Posizionare il disconnettore asettico Clipster® assemblato nell'apposito alloggiamento inserendo per primi i componenti superiori (lato con il logo Sartorius).



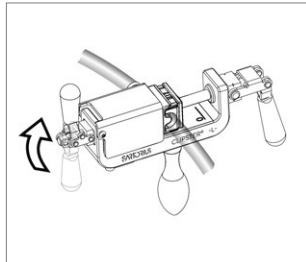
6. Girare la maniglia di serraggio nella posizione 1 per fissare il disconnettore asettico Clipster®.



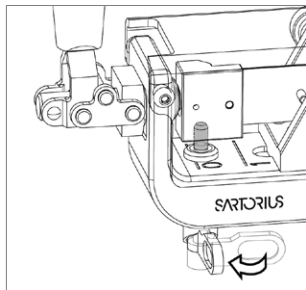
7. Girare il dispositivo di sicurezza per liberare la lama.



8. Tenere la maniglia di serraggio nella posizione 1 e girare la maniglia di taglio fino alla posizione 1 per tagliare il tubo.

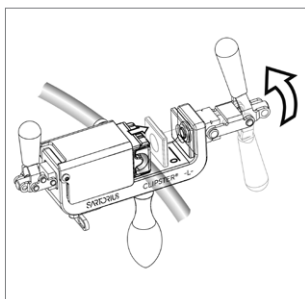


9. Riportare la maniglia di taglio nella posizione 0.

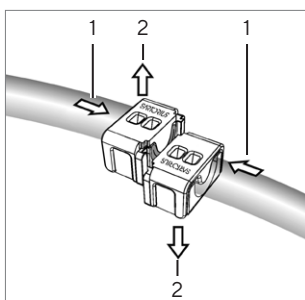


10. Riportare il dispositivo di sicurezza della lama nella posizione iniziale.

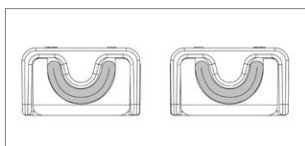
⚠ E' importante riportare il dispositivo di sicurezza della lama nella posizione iniziale sia per proteggere gli operatori da eventuali lesioni sia per usare in modo corretto l'utensile a mano.



11. Riportare la maniglia di serraggio nella posizione 0 e togliere il disconnettore asettico Clipster® dall'utensile a mano.



12. Con il logo rivolto verso l'alto, separare le due estremità premendo verso l'interno e facendole scivolare verticalmente.



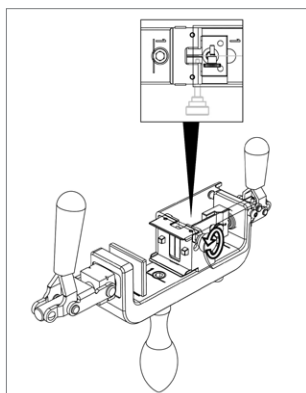
13. Vista delle due estremità dopo lo scollegamento.

4 Manutenzione

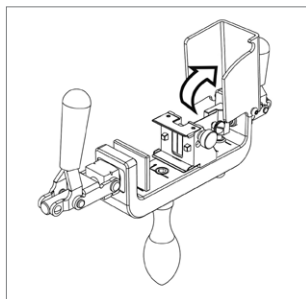
⚠ Gli operatori devono essere addestrati prima di usare il disconnettore aseptico Clipster® e l'utensile a mano.

Per richiedere assistenza tecnica, contattare il proprio rappresentante Sartorius Stedim Biotech.

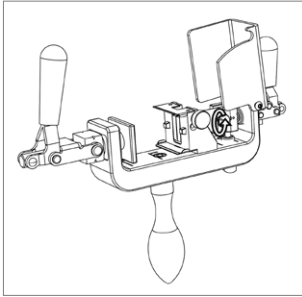
4.1 Procedura per la sostituzione della lama



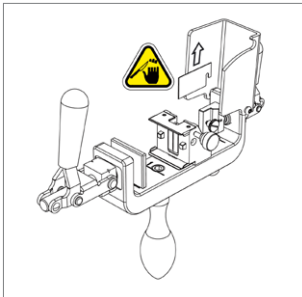
1. Svitare parzialmente la testa della vite n° 2 (vedi Figura 2) presente sulla copertura di protezione, fino allo sganciamento della lama.



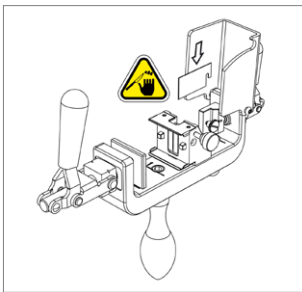
2. Aprire la copertura di protezione.



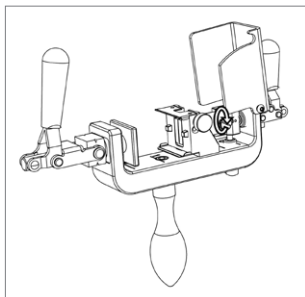
3. Svitare parzialmente la testa della vite n° 1 (vedi Figura 2) presente sulla lama, fino allo sganciamento della lama.



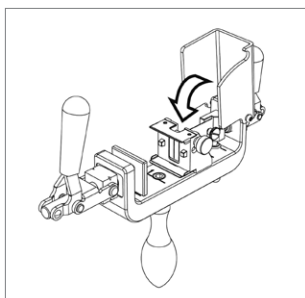
4. Togliere con attenzione la lama.



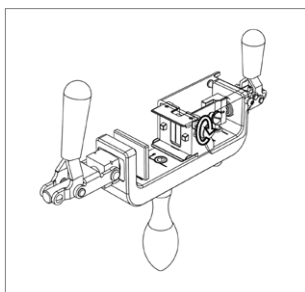
5. Inserire la nuova lama.



6. Stringere la testa della vite n° 1 per fissare la lama.



7. Chiudere la copertura di protezione.



8. Stringere la testa della vite n° 2 sulla copertura di protezione.

4.2 Pulizia dell'utensile a mano

Per la pulizia del disconnettore asettico Clipster® si può usare la maggior parte delle soluzioni disinfettanti come alcool isopropilico.

Índice

1	Instrucciones de uso	60
1.1	Uso previsto	60
1.2	Patentes	60
2	Introducción	61
3	Funcionamiento	63
3.1	Componentes del desconector aséptico Clipster®	63
3.2	Descripción de la herramienta	64
3.3	Advertencias	65
3.4	Procedimiento	65
4	Mantenimiento	70
4.1	Procedimiento de sustitución de la cuchilla	70
4.2	Limpieza de la herramienta de mano	72

1 Instrucciones de uso

Lea este manual atentamente. Contiene información importante para el uso seguro y correcto del desconector aséptico Clipster®.

1.1 Uso previsto

El desconector aséptico Clipster® ha sido desarrollado y fabricado exclusivamente para la desconexión aséptica de tubos en aplicaciones biofarmacéuticas.

 El sistema está diseñado exclusivamente para el uso previsto descrito anteriormente. Cualquier otro uso, fin o aplicación deberá contar previamente con el permiso por escrito del fabricante, o se considerará uso no previsto e inadecuado.

 Sartorius Stedim Biotech no asumirá responsabilidad alguna si el desconector aséptico Clipster® se utiliza de forma indebida o inadecuada, o si se destina a un fin distinto del previsto. Sartorius Stedim Biotech no será responsable de ningún daño que pueda derivarse de dicho uso no previsto. El usuario será enteramente responsable.

El uso previsto abarca también el cumplimiento de las condiciones de funcionamiento, inspección y mantenimiento prescritas por el fabricante. Cualquier incumplimiento de las advertencias contenidas en el manual de instrucciones constituye un uso inadecuado del dispositivo.

1.2 Patentes

Clipster® es una marca comercial de Sartorius Stedim Biotech. El fabricante se reserva el derecho a incorporar, en cualquier momento y sin previo aviso, cualquier cambio o mejora que considere oportunos, tanto en el producto como en los procesos descritos en este manual. Todas las marcas registradas y patentes son propiedad de sus respectivos propietarios.

2 Introducción

El desconectador aséptico Clipster® es un dispositivo de un solo uso desarrollado por Sartorius Stedim Biotech para complementar nuestra gama de productos con la posibilidad de la desconexión aséptica de tubos. El desconectador aséptico Clipster® se puede vender como producto independiente o preensamblado como parte de nuestros conjuntos de bolsas de gestión de fluidos. El desconectador aséptico Clipster® es seguro, rápido y fácil de usar. La desconexión aséptica se realiza con una herramienta de mano que garantiza una ejecución fácil en situaciones con distintas disponibilidades de espacio.

El desconectador aséptico Clipster® se utiliza después de la transferencia de fluidos para desconectar los conductos de transferencia y los conjuntos de bolsa desechables que suelen emplearse en aplicaciones biofarmacéuticas. El desconectador aséptico Clipster® permite la desconexión aséptica en entornos clasificados y no clasificados, manteniendo la esterilidad del producto. Puede emplearse para múltiples tipos y tamaños de tubos.

Especificaciones

Referencia	Denominación	Material
642XS-----D	Desconectador aséptico Clipster® tamaño XS	Polycarbonato
642-S-----D	Desconectador aséptico Clipster® tamaño S	Polycarbonato
642-L-----D	Desconectador aséptico Clipster® tamaño L	Polycarbonato

Referencia	Dimensiones del tubo diámetro int. ext.*	Material del tubo	Material del tubo
		Silicone (Pt) Tuflux® SIL	Silicone (Pt) Pharma-50 Si (Pt)
642XS-----D	1/8" x 1/4"	■	■
642XS-----D	1/4" x 3/8"	■	■
642-S-----D	1/4" x 7/16"		■
642-L-----D	3/8" x 5/8"	■	■
642-L-----D	1/2" x 3/4"		■

Referencia	Dimensiones del tubo diámetro int. ext.*	Material del tubo
		TPE C-Flex®** 374
642XS-----D	1/8" x 1/4"	■
642XS-----D	1/4" x 3/8"	■
642-S-----D	1/4" x 7/16"	■
642-L-----D	3/8" x 5/8"	■
642-L-----D	1/2" x 3/4"	■

Referencia	Denominación
642-S--T	Herramienta de mano para Clipster® XS S
642-S--B	Cuchillas para Clipster® XS S herramienta
642-L--T	Herramienta de mano para Clipster® L
642-L--B	Cuchillas para Clipster® L herramienta

* Nota: El desconector aséptico Clipster® requiere tubos con una longitud mínima de 300 mm.

** C-Flex® es un marca comercial registrada de Saint-Gobain Performance Plastics Corporation.

3 Funcionamiento

3.1 Componentes del desconector aséptico Clipster®

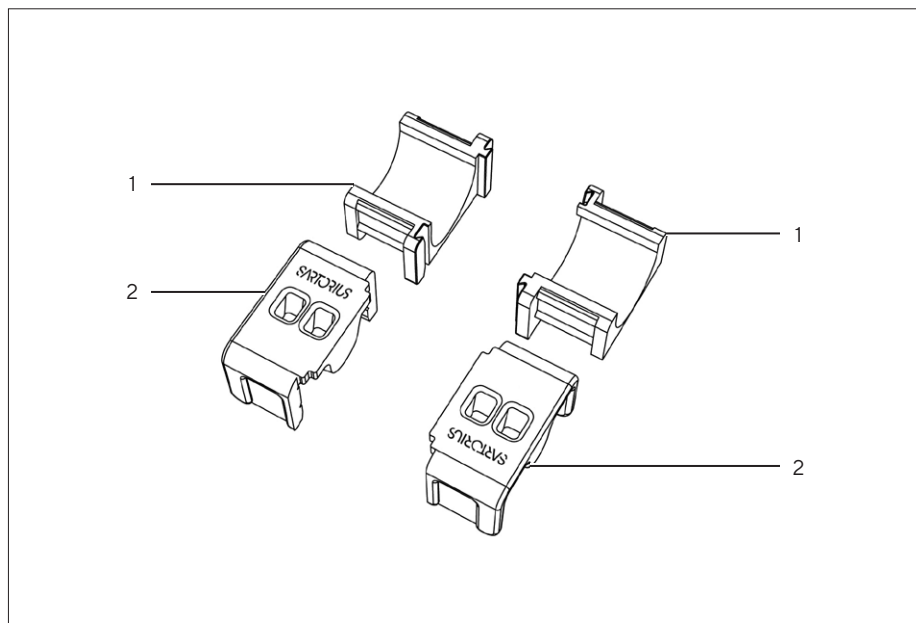


Fig.1

Pos.	Descripción
1	Componente inferior del desconector aséptico Clipster®
2	Componente superior del desconector aséptico Clipster®

3.2 Descripción de la herramienta

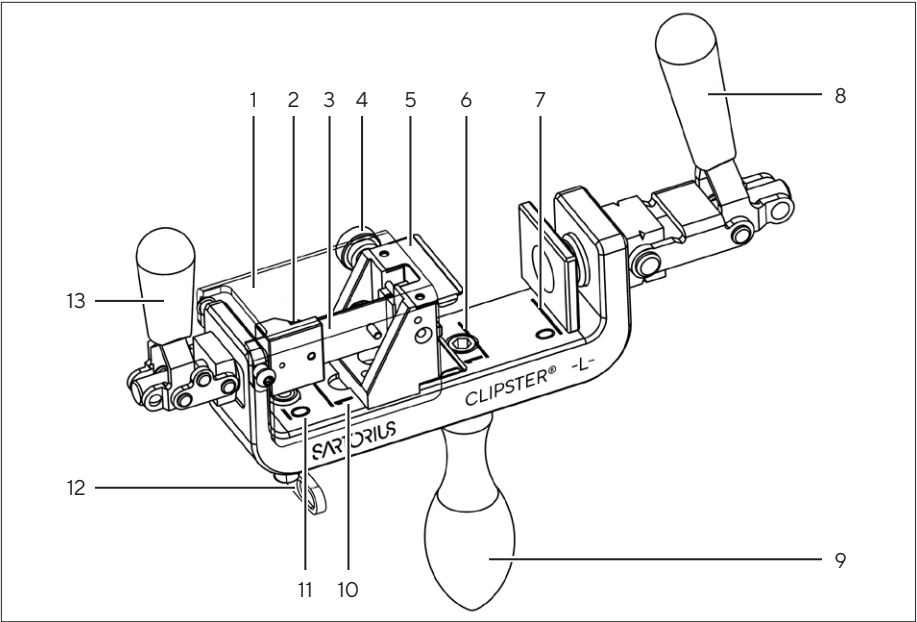


Fig.2

Pos.	Descripción
1	Cubierta protectora
2	Cabeza de tornillo nº 1
3	Cuchilla
4	Cabeza de tornillo nº 2
5	Asiento del Clipster®
6	Posición 1 del mango de abrazadera
7	Posición 0 del mango de abrazadera

Pos.	Descripción
8	Mango de abrazadera
9	Mango de la herramienta
10	Posición 1 del mango de corte
11	Posición 0 del mango de corte
12	Interruptor de seguridad de la cuchilla
13	Mango de corte

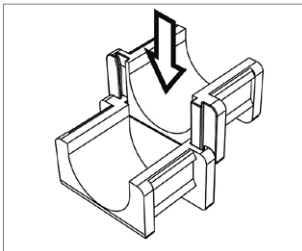
3.3 Advertencias

⚠ Los usuarios deben disponer de los conocimientos adecuados antes de usar el desconector aséptico Clipster® y la herramienta manual.

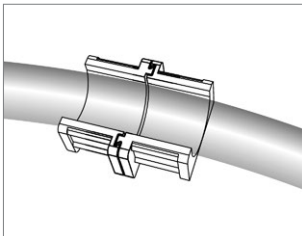
⚠ La cubierta protectora protege a los usuarios contra posibles lesiones causadas por la cuchilla. Sartorius Stedim Biotech no será responsable de ninguna lesión causada por haber retirado la cubierta protectora durante el uso.

3.4 Procedimiento

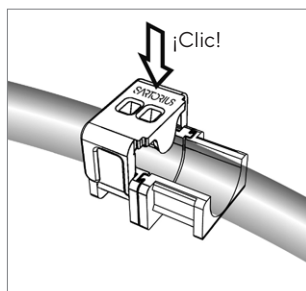
⚠ Para garantizar la desconexión aséptica del tubo, siga meticulosamente los pasos de este procedimiento.



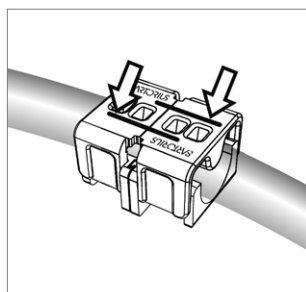
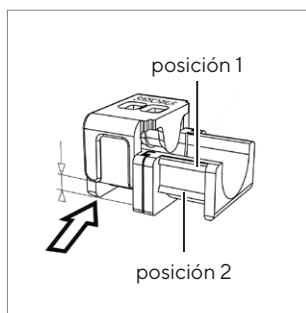
1. Preensamble los dos componentes inferiores deslizándolos uno contra el otro a lo largo de las caras con muescas.



2. Coloque el tubo en ambos componentes inferiores.

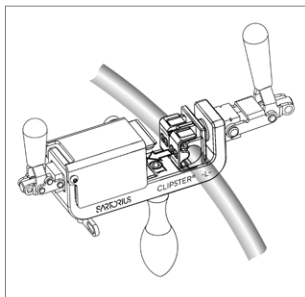


3. Encaje un componente superior en la posición 1 del componente inferior.

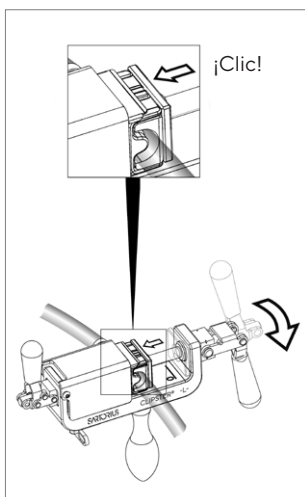
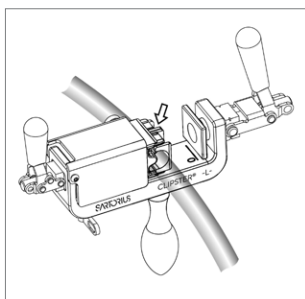


4. Encaje el segundo componente superior en la posición 1 del otro componente inferior y compruebe si el logotipo de Sartorius está alineado correctamente.

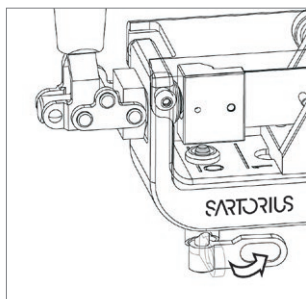
⚠ Antes de usar la herramienta de mano, asegúrese de que los mangos de abrazadera y de corte se encuentren en la posición 0 y de que el interruptor de seguridad de la cuchilla esté accionado.



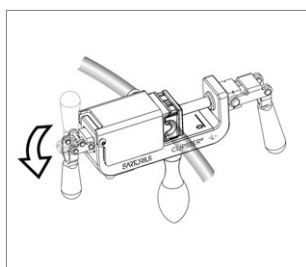
5. Coloque el desconectador aséptico Clipster® preensamblado en el asiento del Clipster® insertando primero los componentes superiores (lado a lado con el logotipo de Sartorius).



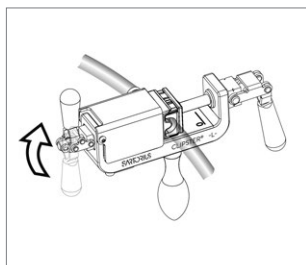
6. Gire el mango de abrazadera a la posición 1 para inmovilizar el desconectador aséptico Clipster®.



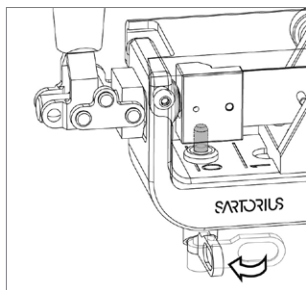
7. Gire el interruptor de seguridad para liberar la cuchilla.



8. Mantenga el mango de abrazadera en la posición 1 y gire el mango de corte a la posición 1 para cortar el tubo.

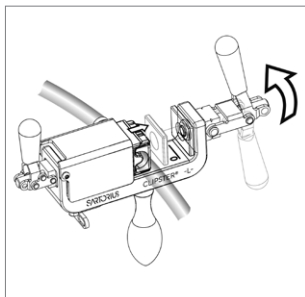


9. Devuelva el mango de corte a la posición 0.

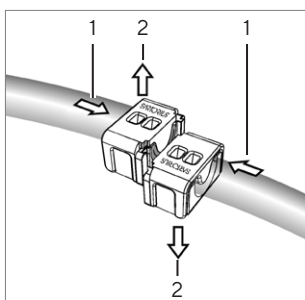


10. Devuelva el interruptor de seguridad de la cuchilla a la posición inicial.

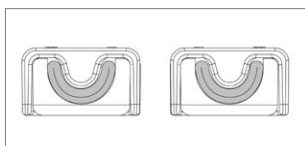
⚠ Es importante devolver el interruptor de seguridad de la cuchilla a la posición inicial para proteger al usuario de posibles lesiones y para emplear correctamente la herramienta de mano.



11. Devuelva el mango de abrazadera a la posición 0 y retire el desconectador aséptico Clipster® de la herramienta de mano.



12. Coloque el logotipo frente a usted y separe los dos extremos empujando por dentro y deslizando verticalmente.



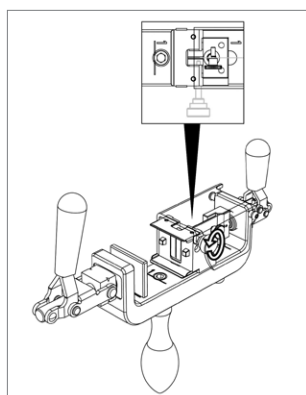
13. Vista de ambos extremos después de la desconexión.

4 Mantenimiento

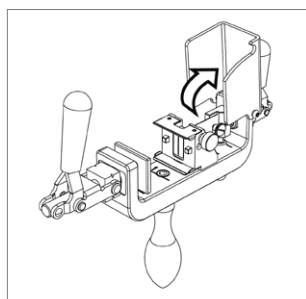
⚠ Los usuarios deben recibir la formación adecuada antes de usar el desconector aséptico Clipster® y la herramienta de mano.

Para obtener asistencia técnica, póngase en contacto con su representante de mantenimiento de Sartorius Stedim Biotech.

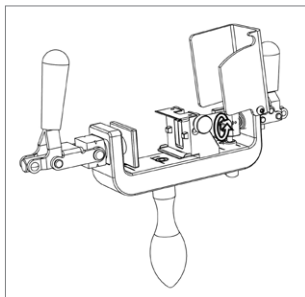
4.1 Procedimiento de sustitución de la cuchilla



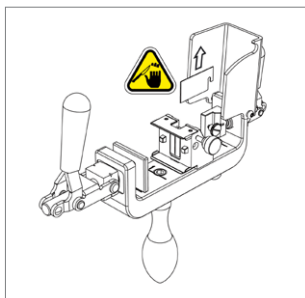
1. Desenrosque parcialmente la cabeza del tornillo n° 2 (ver la figura 2) de la cubierta protectora hasta liberar la cuchilla.



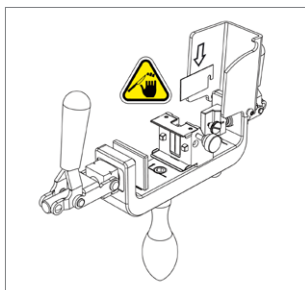
2. Abra la cubierta protectora.



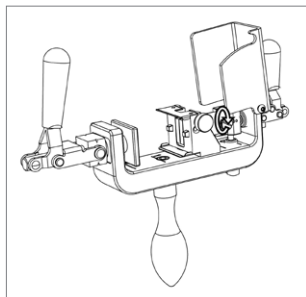
3. Desenrosque parcialmente la cabeza de tornillo n° 1 (ver la figura 2) de la cuchilla hasta liberar ésta.



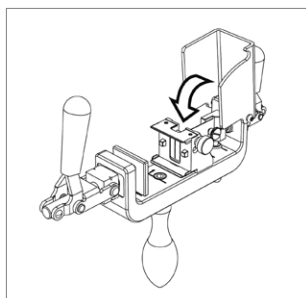
4. Extraiga cuidadosamente la cuchilla.



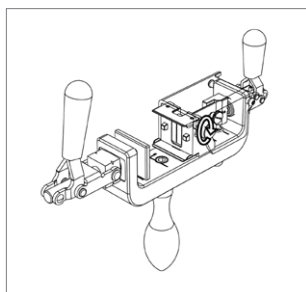
5. Inserte una nueva cuchilla.



6. Apriete la cabeza del tornillo n°1 para asegurar la cuchilla.



7. Cierre la cubierta protectora.



8. Apriete la cabeza del tornillo n°2 en la cubierta protectora.

4.2 Limpieza de la herramienta de mano

El desconectador aséptico Clipster® se puede limpiar con la mayoría de soluciones desinfectantes, como alcohol isopropílico.

Inhoud

1	Gebruiksaanwijzing	74
1.1	Deskundig gebruik	74
1.2	Patenten	74
2	Inleiding	75
3	Gebruik	77
3.1	Onderdelen van de steriele Clipster®-scheidingsinrichting	77
3.2	Beschrijving van de inrichting	78
3.3	Waarschuwingen	79
3.4	Werkstappen	79
4	Onderhoud	84
4.1	Procedure voor vervanging van het mesblad	84
4.2	Handapparaat reinigen	86

1 Gebruiksaanwijzing

Lees deze gebruiksaanwijzing goed door. Deze bevat belangrijke informatie over veilig en succesvol gebruik van de steriele Clipster®-scheidingsinrichting.

1.1 Deskundig gebruik

De steriele Clipster®-scheidingsinrichting werd speciaal ontwikkeld, geconstrueerd en vervaardigd voor het steriele scheiden van slangen in biofarmaceutische toepassingen.

 De inrichting is uitsluitend bestemd voor het bovengenoemde doel. Elke andere toepassing buiten deze wordt zonder schriftelijke toestemming van de fabrikant beschouwd als oneigenlijk gebruik.

 Sartorius Stedim Biotech aanvaardt geen aansprakelijkheid wanneer de steriele Clipster®-scheidingsinrichting ondeskundig is ingezet of gebruikt. Sartorius Stedim Biotech aanvaardt geen aansprakelijkheid voor hieruit voortvloeiende schade. Het risico ligt uitsluitend bij de gebruiker.

Deskundig gebruik houdt ook in dat de werk-, inspectie- en onderhoudsvoorschriften van de fabrikant worden opgevolgd. Het niet in acht nemen van de waarschuwingen in de handleiding geldt als oneigenlijk gebruik van het apparaat.

1.2 Patenten

Clipster® is een handelsmerk van Sartorius Stedim Biotech. De fabrikant behoudt zich het recht voor om te allen tijde en zonder voorafgaande kennisgeving wijzigingen en verbeteringen aan het product en de in deze gebruiksaanwijzing beschreven procedures aan te brengen. Alle geregistreerde handelsmerken en patenten zijn eigendom van hun respectievelijke eigenaars.

2 Inleiding

De steriele Clipster®-scheidingsinrichting is een eenweg-inrichting van Sartorius Stedim Biotech, die een aanvulling vormt op ons productassortiment door het steriele scheiden van slangen. De steriele Clipster®-scheidingsinrichting kan als afzonderlijk product of samen met onze Fluidmanagement-zakeenheid worden verkocht. De steriele Clipster®-scheidingsinrichting is veilig, snel en eenvoudig in gebruik. De steriele scheiding vindt plaats met een handapparaat, dat zorgt voor een eenvoudige uitvoering bij uiteenlopende ruimte-omstandigheden.

De steriele Clipster®-scheidingsinrichting wordt na de overdracht van de fluid gebruikt om eenweg-transferleidingen en -zakeenheden in biofarmaceutische toepassingen te scheiden. De steriele Clipster®-scheidingsinrichting maakt een steriele scheiding in niet-geclassificeerde en geclassificeerde omgevingen mogelijk en waarborgt productsteriliteit. Hij kan worden gebruikt voor verschillende soorten en maten slangen.

Technische gegevens

Referentie	Aanduiding	Materiaal
642XS-----D	Clipster® steriele scheidingsinrichting maat XS	polycarbonaat
642-S-----D	Clipster® steriele scheidingsinrichting maat S	polycarbonaat
642-L-----D	Clipster® steriele scheidingsinrichting maat L	polycarbonaat

Referentie	Slangmaat ID – OD*	Slangmateriaal	Slangmateriaal
		Siliconen (Pt) TuFlux® SIL	Siliconen (Pt) Pharma-50 Si (Pt)
642XS-----D	1/8" x 1/4"	■	■
642XS-----D	1/4" x 3/8"	■	■
642-S-----D	1/4" x 7/16"		■
642-L-----D	3/8" x 5/8"	■	■
642-L-----D	1/2" x 3/4"		■

Referentie	Slangmaat ID – OD*	Slangmateriaal
		TPE C-Flex®** 374
642XS-----D	1/8" x 1/4"	■
642XS-----D	1/4" x 3/8"	■
642-S-----D	1/4" x 7/16"	■
642-L-----D	3/8" x 5/8"	■
642-L-----D	1/2" x 3/4"	■

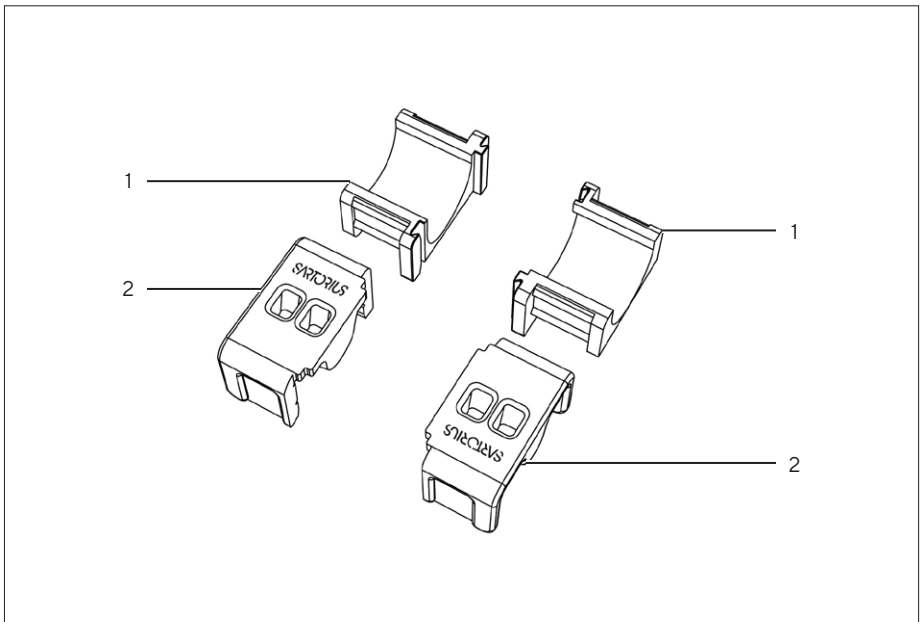
Referentie	Aanduiding
642-S--T	Handgereedschapapparaat voor Clipster® XS S
642-S--B	Mesbladen voor Clipster® XS S gereedschap
642-L--T	Handgereedschapapparaat voor Clipster® L
642-L--B	Mesbladen voor Clipster® L gereedschap

* Aanwijzing: Voor de steriele Clipster®-scheidingsinrichting zijn slangen met een minimumlengte van 300 mm nodig.

** C-Flex® is een gedeponeerd handelsmerk van Saint-Gobain Performance Plastics Corporation.

3 Gebruik

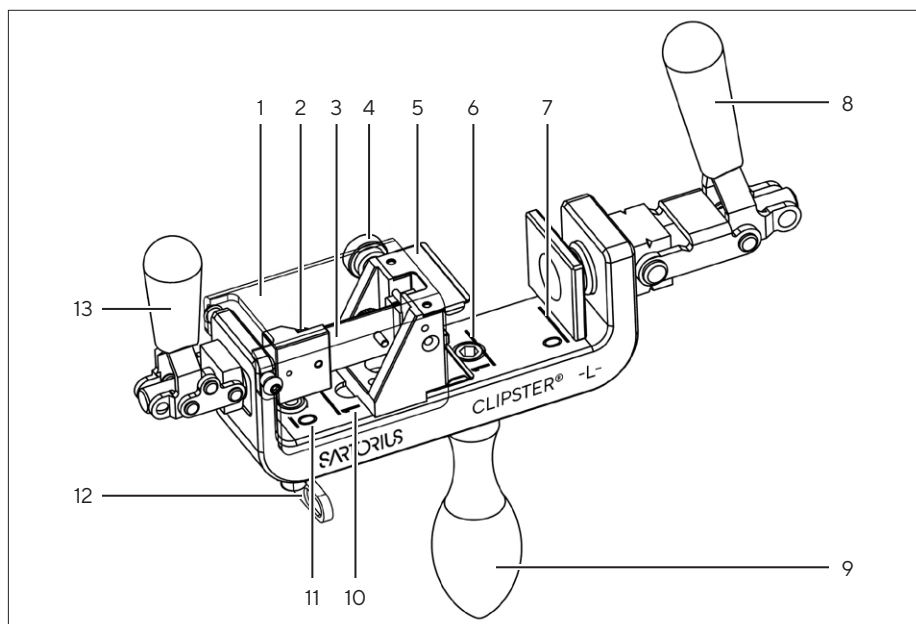
3.1 Onderdelen van de steriele Clipster®-scheidingsinrichting



Afb.1

Pos.	Beschrijving
1	Onderste deel van de steriele Clipster®-scheidingsinrichting
2	Bovendeel van de steriele Clipster®-scheidingsinrichting

3.2 Beschrijving van de inrichting



Afb.2

Pos.	Beschrijving
1	Beschermingsfadedkking
2	Schroefkop nr. 1
3	Mesblad
4	Schroefkop nr. 2
5	Clipster®-zitting
6	Positie 1 van de klemgreep
7	Positie 0 van de klemgreep

Pos.	Beschrijving
8	Klemgreep
9	Greep van de inrichting
10	Positie 1 van de snijgreep
11	Positie 0 van de snijgreep
12	Mesblad-veiligheidsschakelaar
13	Snijgreep

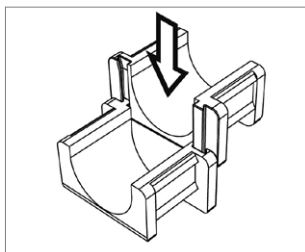
3.3 Waarschuwingen

⚠ De gebruikers moeten worden getraind voorafgaand aan het gebruik van de steriele Clipster®-scheidingsinrichting en het handapparaat.

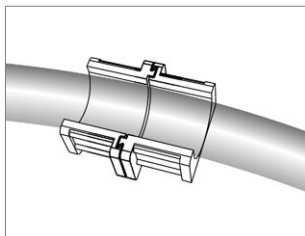
⚠ De beschermingsafdekking beschermt de gebruiker tegen mogelijk letsel door het mesblad. Sartorius Stedim Biotech aanvaardt geen aansprakelijkheid wanneer deze beschermingsafdekking tijdens het gebruik is verwijderd.

3.4 Werkstappen

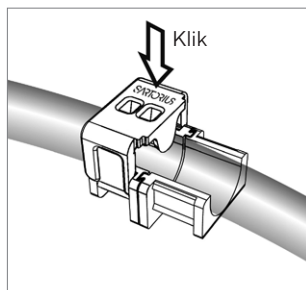
⚠ Steriele scheiding van slangen is alleen mogelijk als de volgende stappen zorgvuldig worden gevolgd.



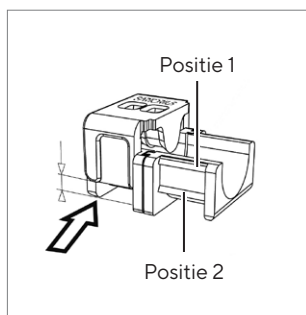
1. Schuif de beide onderste delen bij de inkepingen in elkaar.



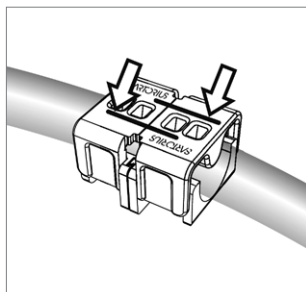
2. Leg de slang in de beide onderste delen.



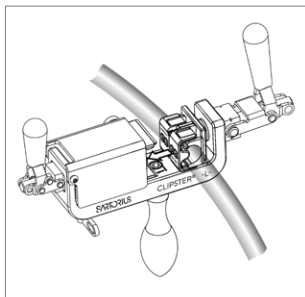
3. Klik een bovendeel vast in positie 1 van het onderste deel.



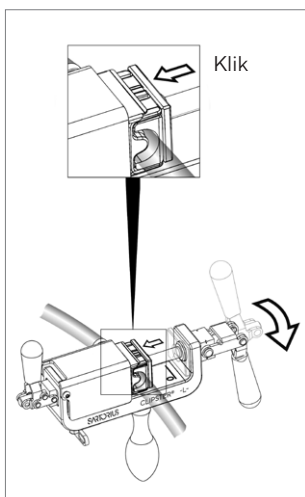
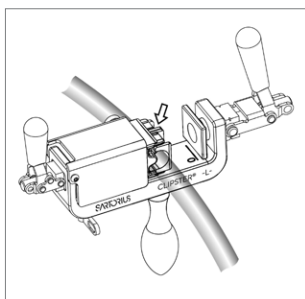
4. Klik het tweede bovendeel vast op positie 1 van het andere onderste deel en controleer visueel of het logo van Sartorius correct is uitgelijnd.



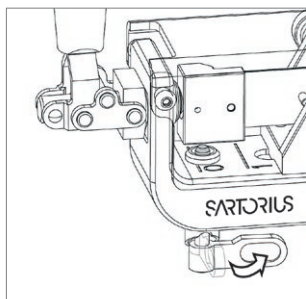
⚠ Vóór gebruik van het handapparaat moeten de klem- en de snijgreep in positie 0 staan en moet het veiligheidsmesblad zijn vastgeklit.



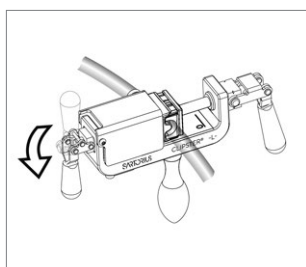
5. Plaats de gemonteerde steriele Clipster®-scheidingsinrichting met de bovendelen (kant met het logo van Sartorius) eerst in de Clipster®-zitting.



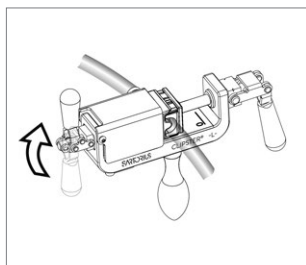
6. Draai de klemgreep naar positie 1 om de steriele Clipster®-scheidingsinrichting vast te klemmen.



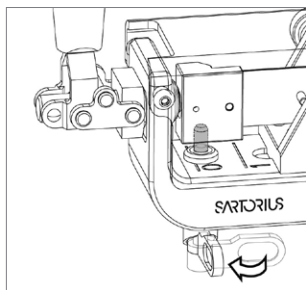
7. Draai de veiligheidsschakelaar om het mesblad los te maken.



8. Houd de klemgreep in positie 1 en draai de snijgreep naar stand 1 om de slang door te snijden.

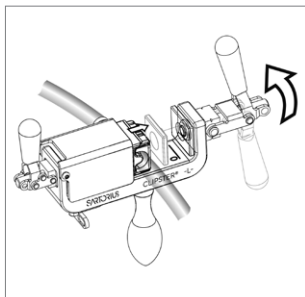


9. Zet de snijgreep terug in positie 0.

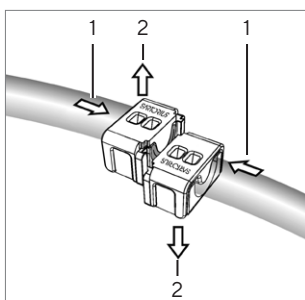


10. Zet de veiligheidsschakelaar van het mesblad in de uitgangspositie.

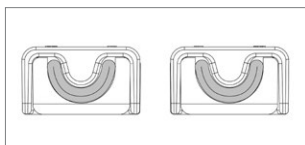
 Het is belangrijk om de veiligheidsschakelaar van het mesblad terug te zetten in zijn uitgangspositie, om gebruikers te beschermen tegen letsel en om het handapparaat op de juiste manier te gebruiken.



11. Zet de klemgreep terug in de positie 0 en haal de steriele Clipster®-scheidingsinrichting uit het handapparaat.



12. Draai het logo in uw richting en trek de beide uiteinden uit elkaar door ze naar binnen te drukken en omhoog te schuiven.



13. Controleer beide uiteinden na het scheiden.

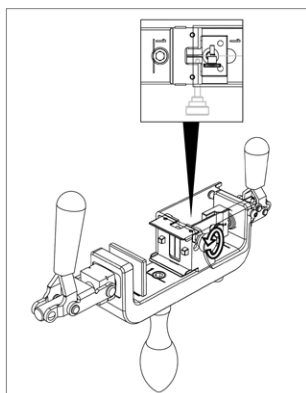
4 Onderhoud

 De gebruikers moeten worden getraind voorafgaand aan het gebruik van de steriele Clipster®-scheidingsinrichting en het handapparaat.

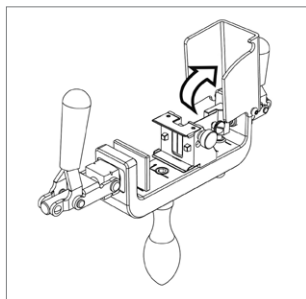
Technische ondersteuning krijgt u van uw contactpersoon bij Sartorius Stedim Biotech.

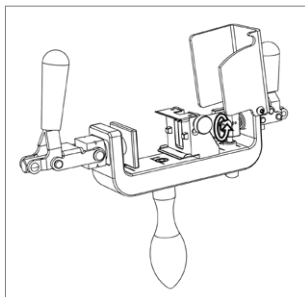
4.1 Procedure voor vervanging van het mesblad

1. Draai de schroefkop nr. 2 (zie afbeelding 2) op de beschermingsafdekking iets los, totdat het mesblad vrijkomt.

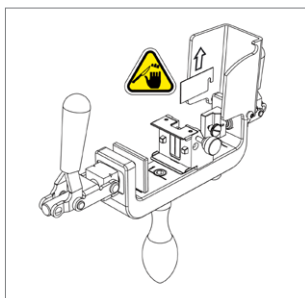


2. Open de beschermingsafdekking.

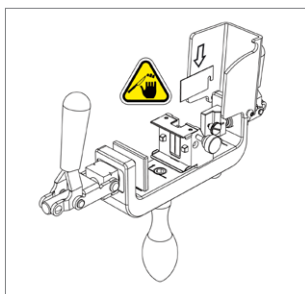




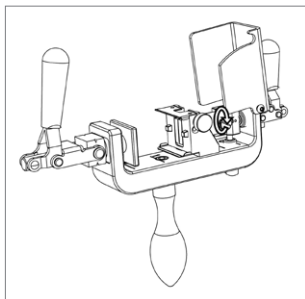
3. Draai de schroefkop nr. 1 (zie afbeelding 2) op het mesblad iets los, totdat het mesblad vrijkomt.



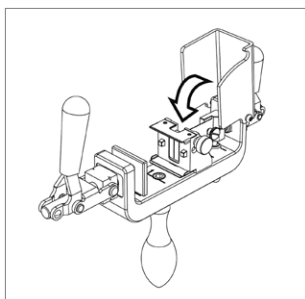
4. Verwijder het mesblad voorzichtig.



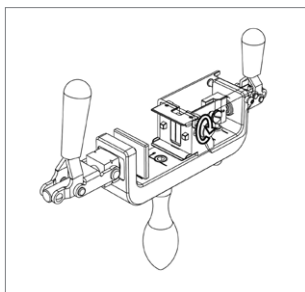
5. Plaats een nieuw mesblad.



6. Draai de schroefkop nr. 1 vast om het mesblad vast te zetten.



7. Sluit de beschermingsafdekking.



8. Draai de schroefkop nr. 2 vast op de beschermingsafdekking.

4.2 Handapparaat reinigen

De steriele Clipster®-scheidingsinrichting kan worden gereinigd met de meeste desinfecterende oplossingen, zoals isopropylalcohol.

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone: +49 551 308 0
www.sartorius.com

The information and figures contained in these instructions correspond to the version date specified below.

Sartorius reserves the right to make changes to the technology, features, specifications and design of the equipment without notice.

Masculine or feminine forms are used to facilitate legibility in these instructions and always simultaneously denote the other gender as well.

Copyright notice:

This instruction manual, including all of its components, is protected by copyright.

Any use beyond the limits of the copyright law is not permitted without our approval.

This applies in particular to reprinting, translation and editing irrespective of the type of media used.

© Sartorius Germany

Last updated:
12 | 2020