



Lynx®S2S Connector

User Guide

The M logo and Lynx are registered trademarks of Merck KGaA, Darmstadt, Germany. All trademarks of third parties are the property of their respective owners. © 2007, 2015 EMD Millipore Corporation. All rights reserved. 00100285PU, Rev. B, 02/2015

Technical Assistance

For more information, contact the office nearest you or visit the Technical Service page at www.emdmillipore.com/techservice. Worldwide contact information is available at www.emdmillipore.com/offices.

Standard Warranty

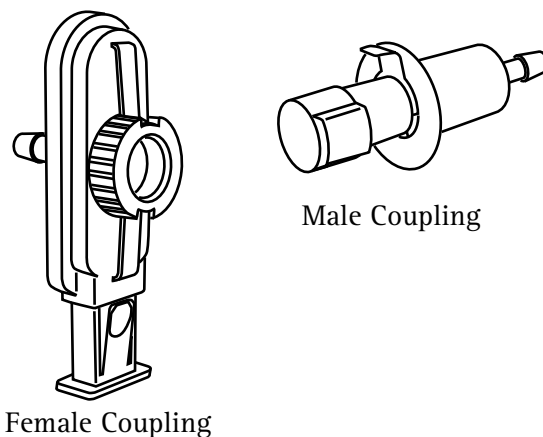
The applicable warranty for the products listed in this publication may be found at www.millipore.com/terms (within the “Terms and Conditions of Sale” applicable to your purchase transaction).

Introduction

The Lynx®S2S (sterile to sterile) connector is a single-use, single actuation, gamma and autoclave compatible disposable connector.

The Lynx®S2S connector includes a female coupling and a male coupling. Once these two components are assembled together, the sterile fluid path is enabled.

Lynx®S2S Connector components



Sterilization Instructions

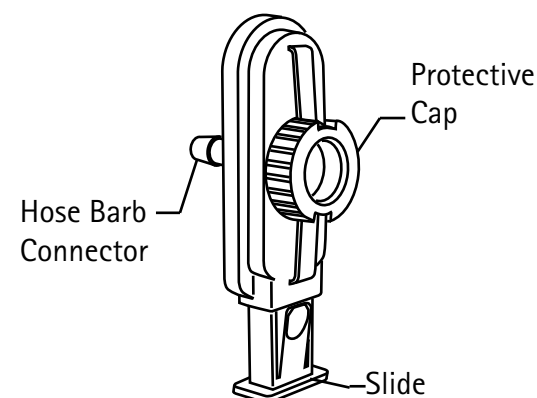
- The Lynx®S2S connector may be sterilized using gamma irradiation to a maximum dose of 45 kGy.
- The male and female connectors may be autoclaved at 121 to 130° C for up to 30 minutes.
- All system components installed on the Lynx®S2S connector and the packaging must be compatible with the chosen sterilization method.
- DO NOT STEAM-IN-PLACE.

Installation Instructions

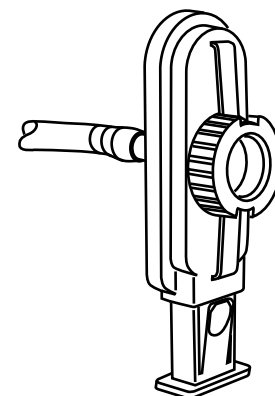
- Do not submerge components in liquids.
- Do not pressurize or attempt to flow liquid through couplings until actuated.

Installing the Female Coupling onto the Disposable Assembly

1. Remove coupling from outer packaging.
2. Remove from inner packaging.
DO NOT PUSH SLIDE INTO COUPLING BODY. DO NOT REMOVE CONNECTOR CAP.
3. Insert hose barb into tubing.



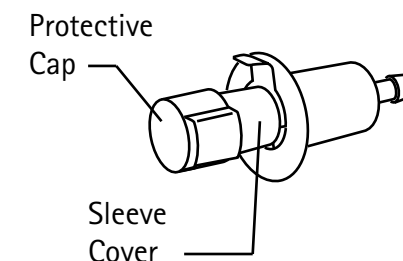
4. If required, install a user-supplied tubing fastener (clamp or tie wrap) over tubing to secure connection.



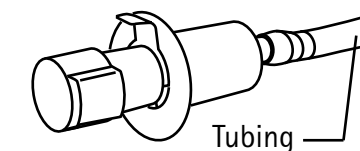
5. Verify that the protective cap is in place by twisting it clockwise
6. Package and sterilize the assembly by either gamma or autoclave.

Installing the Male Coupling onto the Disposable Assembly

1. Remove coupling from packaging.
DO NOT REMOVE SLEEVE COVER.
DO NOT REMOVE CONNECTOR CAP.



2. Insert hose barb into tubing.



3. If required, install a user-supplied tubing fastener (clamp or tie wrap) over tubing to secure connection.
4. Verify that the protective cap is in place by twisting it clockwise
5. Package and sterilize the assembly by either gamma or autoclave.

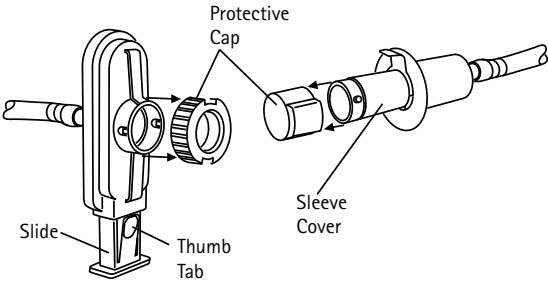
Actuating the Lynx®S2S Connector

The Lynx®S2S Connector is a single use unit. Once the female and male couplings are assembled they cannot be disassembled. No special clean room equipment is required to ensure a sterile connection.

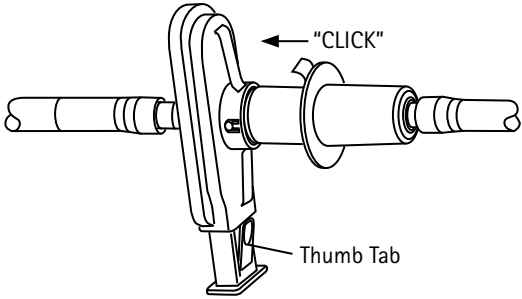
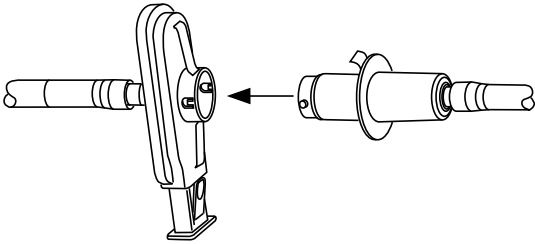
NOTE

Do not initiate fluid flow to the couplings until male and female couplings are actuated. If necessary, use a tubing clamp to keep fluid and pressure away from the couplings until actuation steps are complete.

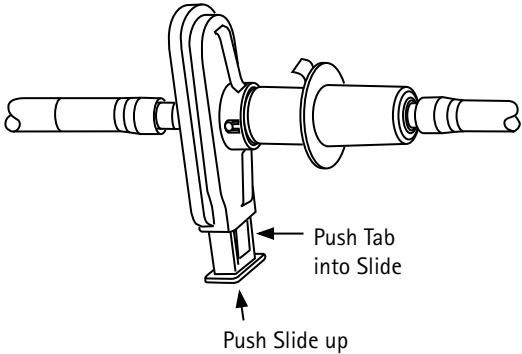
1. Remove protective caps from the female and male couplings by turning the caps counter clockwise. (The protective sleeve cover remains on the male coupling.)



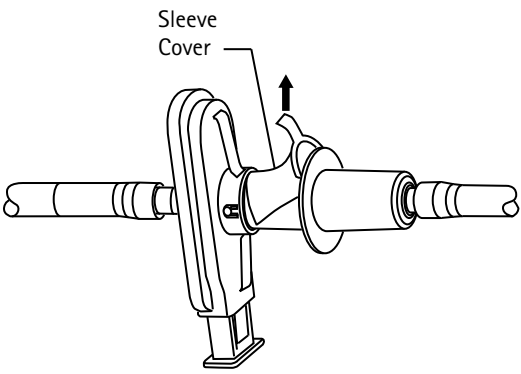
2. Insert the male coupling into the female coupling aligning the pins on the male coupling with the holes in the female coupling. A “click” will be heard when the components are properly connected.



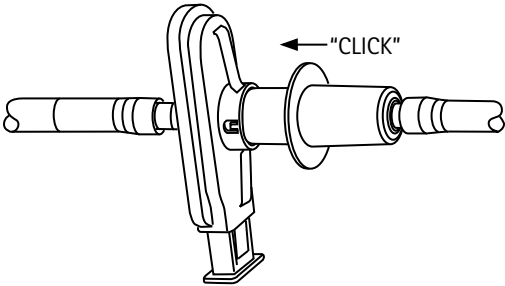
3. Press the thumb tab on the slide, then push the slide into the body of the female coupling until it stops and a “click” is heard.



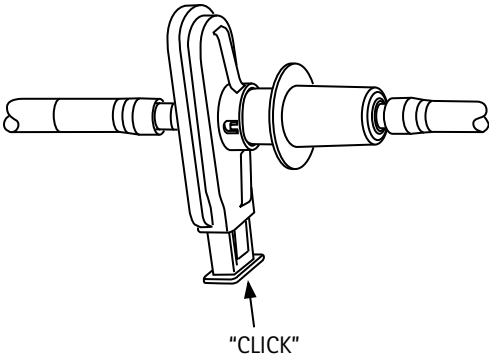
4. Pull the tab on the sleeve cover to remove the cover.



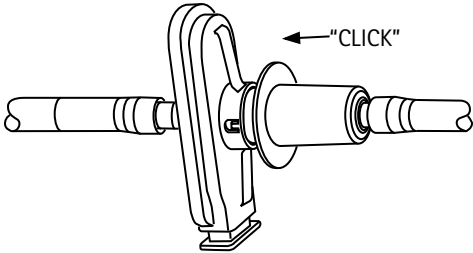
5. Push male coupling into body of female coupling until it stops and a “click” is heard.



6. Push slide completely into body of female coupling until it stops and a “click” is heard.



7. Push male coupling into female coupling until a “click” is heard. If a “click” is not heard the connection is not actuated and leakage or breach of sterility may occur. If a “click” is not heard, push male coupling into female coupling again until a “click” is heard. If unit is still not actuated, do not use.



8. Sterile fluid path is now ready for fluid flow.