

SART System™

Sartorius Aseptic Rapid Transfer System for Fluid Transfer



Description

The SART System™ is a connecting system that allows safe liquid transfer across a wall between two areas with different containment classifications. It consists of an external port and an internal port to be used in combination with a disposable connection device—the Gammasart ATD™.

Applying Rapid Transfer Port (RTP) technology at the port – as well as the disposable connector, the SART System™ utilizes a well-established aseptic connection technology. The principle of the connection is based on the alpha | beta concept using four V-shaped profiles matching exactly at the tip.

Key Features and Benefits

100% air leak test of the connector device and the port	Assurance of integrity
Extensive qualification of the systems	Proven robustness
Gammasart ATD™ allows up to 3 connections disconnections of traditional vessels or fully disposable systems.	Flexible operations
Integrated securities	Error-proof and safe aseptic transfer system

Applications

The SART System™ has been developed to facilitate liquid aseptic transfer from:

- a clean room into a protected environment (Isolator, Restricted Access Barrier System | RABS).
- a lower classified clean room or even corridor into a class A-ISO 5 clean room



Highest assurance of integrity

The SART System™ and the Gammasart ATD™ connector have been extensively qualified for use in critical applications within biopharmaceutical manufacturing processes. The Gammasart ATD™ connectors are 100 percent air leak-tested as a condition of lot release. In addition, the internal port of the SART System™ is tested at 350 mBars. A validation guide is available.

Proven robustness

The SART System™ is based on a well-established alpha | beta concept to allow for reliable and secured fluid transfer while maintaining the integrity of the critical area—isolators, RABS, and cleanrooms. The SART System™ has been extensively qualified to ensure the robustness of the connection and grade-A environment continuity for aseptic processing. The Gammasart ATD™ connector design has been reviewed to improve the tightness of the connection, providing increased robustness of the assembly.

Flexible, single-use devices

The disposable Gammasart ATD™ connector may be opened and re-used up to three times. Multiple transfers of different volumes can be processed without the need for re-sterilization. The disposable Gammasart ATD™ connector may be sterilized by autoclave or gamma irradiation, enabling use with either traditional vessels or fully disposable fluid-handling technologies.

Error-proof and safe aseptic transfer system

The port is equipped with mechanical interlock systems that prevent an accidental opening without the Gammasart ATD™ connector in place and accidental release of the outer part of the connector. In the event of a stoppage on a filling line, the Gammasart ATD™ connector allows the bulk to be safely removed from the line, stored, and reconnected later.

Technical Specifications

External Port

Port material	Stainless Steel 316L
Seal material	Silicone
Installation requirements	Wall thickness: 3 – 6 mm 0.08 – 0.23 in. Not suitable for installation on glass wall.
Dimensions	Diameter : 159 mm Thickness: 24 mm



SART System™ external port

Internal Port

Material	Stainless Steel 316L, PEEK
Dimensions	Length with straight arm: 200 mm Radius when arm claimed: 110 mm



SART System™ internal port

Gammasart ATD™ Connector

Connector body and cover material	PBT Celanex Grade 2404MT
Overmolded seal material	Thermoplastic Elastomer (TPE) THERMOLAST® M TM6ADT_S102

Sterilization

Supplied non-sterile or gamma-irradiated at 25 – 45kGy. Irradiated devices shall not be re-irradiated.

Gamma irradiation	≤ 50kGy
Autoclaving	Can afford sterilization by autoclave. Max recommended treatment: 30 min at 121°C (250°F)

Dimensions

External side	3/8" hose barb connection
Inner side	3/8" tube connection



Gammasart ATD™ connector

Ordering Information

Part Number	Description	Qty/Box
AN-CON-385025	SART System™, Sartorius Aseptic Rapid Transfer System, including external port, external port cover, internal port, internal stopper	1
AN-CON-122025	Gammasart ATD™, Non-sterile disposable connector for aseptic liquid transfer, including connector body and connector cover	50
AN-CON-121025	Gammasart ATD™, Gamma-irradiated disposable connector for aseptic liquid transfer, including connector body and connector cover	10

Library

Further documentation is available on request:

- SART System™ Validation Guide
- SART System™ User Manual

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Straße 11
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 For further contacts, visit
www.sartorius.com