

StreamLink® CC 15

Automated, High-Throughput,
and Efficient mAb Sample
Preparation

Product Information

Introducing StreamLink® CC 15: The next generation in cell line and process development. Simplify mAb sample preparation without compromising quality with an automated high-throughput system that offers fast clarification and purification of 5 – 15 mL samples ready for critical quality attribute (CQA) analysis.

Process 24 – 48 samples directly from Ambr® 15 | Ambr® 250 bioreactors to cut processing time from days to hours and reduce manual input from 6 hours to just 10 minutes.



Features and Benefits

- **High throughput:** Clarification and | or purification of 24 × 10 ± 5 mL culture samples in 2 hours.
- **All-in-one system:** Achieve a streamlined, controlled, and reproducible clarification and purification process in one step.
- **Automation:** Automated walk-away process with integrated error recoveries to save time and reduced full time equivalent (FTE).

Introduction

Relevant Applications

- Cell line development
(harvest | clarification and purification)
- Process development (sample prep for CQAs)
- Process characterization (sample prep for CQAs)
- Molecule identification | transient-expressing cell lines
(material generation for molecule testing)

Relevant Process Steps

- mAb sample preparation in cell line development
- Clarification | harvest
- Chromatography | capture and polish

Technical Specifications

StreamLink® CC 15

Dimension W×D×H [mm]	530×1,220×660
Weight [kg]	50
Nominal input volume [mL]	5 - 15
Maximum input volume [mL]	30
Minimum input volume [mL]	3
Maximum sample throughput	12 samples per hour
Purification consumable	Sartobind® Rapid A Nano
Clarification consumable	Sartoclear® Disc
Maximum input samples	48
Maximum system pressure [bar]	4
Maximum pumped flow rate [mL/min]	75
Number of labware holder positions	6
Maximum clarification filter holder capacity (per filter station)	30
Liquid handler deck bottles	5
Liquid handler in-line supply bottles	3
Filter station supply bottles (excl. unused air)	6



Ordering Information

Item	Description	Order Number
StreamLink® CC 15 System	<p>One complete StreamLink® CC 15 system with all included components packaged in one outer box on a pallet</p> <p>Inside the system box will be the following:</p> <ul style="list-style-type: none"> ▪ 2× Filter station (each FS will have a unique serial number) ▪ Liquid handler (each LH will have a unique serial number) ▪ Controller ▪ Accessories: <ul style="list-style-type: none"> ▪ 061-2B20 Feed Waste bottle set ▪ 061-2B21 Feed Waste bottle lid set ▪ 061-2B22 Feed connection tubing set ▪ 061-2B23 Liquid handler bottle set ▪ 061-2B24 Liquid handler bottle lids ▪ 061-2B26 Syringe pump pH probe assembly <p>User Manual in six languages (English, German, Spanish, French, Chinese and Italian)</p>	061-8B01-P

Peripherals and Accessories

Product Name	Order Number
Ambr® 15 vessel holder including lid	061-8B50
Microplate holder	061-8B51
50 mL centrifuge tube holder including lid	061-8B52
15 mL centrifuge tube holder including lid	061-8B53
Lid for centrifuge tube holder	061-8B61
Lid for Ambr® 15 vessel holder	061-8B60
Feed Waste bottle set	061-2B20
Feed Waste bottle lid set	061-2B21
Feed Connection tubing set	061-2B22
Liquid handler bottle set	061-2B23
Liquid handler bottle lids	061-2B24
Custom bottle and tube set	061-2B25
Syringe pump pH probe assembly	061-2B26
Levelling Platform	061-8B30
Power Cable set - UK plug type	061-8B10
Power Cable set - EU plug type	061-8B11
Power Cable set - US plug type	061-8B12

Consumables	Description	Order Number
Sartoclear® Disc (Box of 24) for clarification	Sartoclear® Disc is a single-use filter, with 20 cm ² filtration area. Each StreamLink® CC 15 system can filter up to 24 samples in 2 hours, which requires 24 Sartoclear® Discs. Sartoclear® Discs are provided in boxes of 24 pieces.	29DF020-DIA----2
Sartobind® Protein A Nano (Box of 2) for purification	Sartobind® Rapid A Nano is a disposable membrane adsorber, with 1.2 mL membrane volume. Each StreamLink® CC 15 system can filter up to 24 samples in 2 - 4 hours, which requires 2 Sartobind® Rapid A devices. Sartobind® Rapid A Nano are provided in boxes of 2 pieces	96R-PA19D-T11--V
Sartobind® Protein A Nano (Box of 4) for purification	Sartobind® Rapid A Nano are provided in boxes of 2 pieces or 4 pieces.	96R-PA19D-T11--A

United Kingdom

Sartorius Stedim TAP
Grantham CI
York Way
Royston SG8 5WY

 **For more information, visit**
www.sartorius.com