



HiScreen™ Capto™ Core 400

HiScreen™ Capto™ Core 700

HiTrap™ Capto™ Core 400

HiTrap™ Capto™ Core 700

## Prepacked columns

### Instructions for Use

The Capto™ Core 400 and Capto Core 700 chromatography resins are designed for intermediate purification and polishing of viruses and other large biomolecules in flow-through mode. The products are based on the core bead concept. Each bead has a ligand-activated core and an inactive shell. The shell prevents larger molecules from entering the core, but smaller proteins and impurities can enter the core and bind to the hydrophobic and positively charged octylamine ligands.

The prepacked columns are used for optimization of methods and parameters, such as sample load and binding conditions, as well as for small-scale purifications. The columns provide fast and reproducible separations in convenient formats. The columns are optimally used with liquid chromatography systems such as ÄKTA™.

# Table of Contents

1	Product description .....	3
2	Optimization .....	7
3	Operation .....	10
4	Cleaning-In-Place (CIP) .....	15
5	Scale-up .....	16
6	Adjusting pressure limits .....	17
7	Storage .....	20
8	Troubleshooting .....	21
9	Ordering information .....	22
10	Related literature on the web .....	23

## Important

Read these instructions carefully before using the products.

## Safety

For use and handling of the products in a safe way, refer to the Safety Data Sheets.

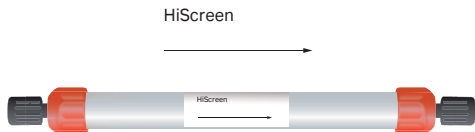
## Intended use

The products are intended for research use only, and shall not be used in any clinical or *in vitro* procedures for diagnostic purposes.

# 1 Product description

## HiScreen column characteristics

HiScreen™ columns are made of biocompatible polypropylene that does not interact with biomolecules. The arrow on the column label shows the recommended flow direction.



**Fig 1.** HiScreen column.

**Note:** Do not open or refill HiScreen columns.

**Note:** Make sure that the column is sealed properly to prevent leakage when stored.

If scale-up requires a larger bed height, two columns can be connected in series using a union to give a 20 cm bed height (see [Chapter 5 Scale-up, on page 16](#)).

**Table 1.** Characteristics of HiScreen column

Column volume (CV)	4.7 mL
Column dimensions	0.77 × 10 cm
Column hardware pressure limit	0.8 MPa (8 bar, 116 psi)

**Note:** The pressure over the packed bed varies depending on parameters such as resin characteristics, sample/liquid viscosity, and column tubing used.

## HiTrap column characteristics

HiTrap™ columns are made of biocompatible polypropylene that does not interact with biomolecules.

The columns are delivered with a stopper at the inlet and a snap-off end at the outlet. The first table below lists the characteristics of HiTrap columns.



**Fig 2.** HiTrap, 1 mL column.

**Note:** *HiTrap columns cannot be opened or refilled.*

**Note:** *Make sure that the connector is tight to prevent leakage.*

**Table 2.** Characteristics of HiTrap columns.

Column volume (CV)	1 mL
Column dimensions	0.7 × 2.5 cm
Column hardware pressure limit	0.5 MPa (5 bar, 72 psi)

**Note:** *The pressure over the packed bed varies depending on parameters such as resin characteristics, sample/liquid viscosity, and column tubing used.*

## Supplied Connector kit with HiTrap column

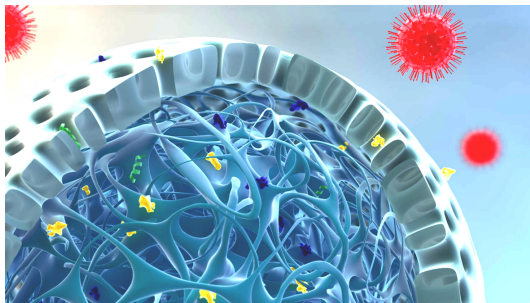
Connectors supplied	Usage	No. supplied
Union 1/16" male/luer female	For connection of syringe to HiTrap column	1
Stop plug female, 1/16"	For sealing bottom of HiTrap column	2, 5, or 7

## Resin properties

Capto Core 400 and Capto Core 700 are composed of a ligand-activated core and an inactive shell. The inactive shell excludes large molecules (average cut-offs:  $M_r \sim 400\,000$  for Capto Core 400, and  $M_r \sim 700\,000$  for Capto Core 700) from entering the core through the pores of the shell, see image below. These larger molecules are collected in the column flow-through, while smaller impurities bind to the internalized ligands.

The core of each bead is functionalized with ligands that are both hydrophobic and positively charged, resulting in a highly efficient multimodal binding of various impurities that are small enough to enter the core. The octylamine ligand ensures strong binding with most impurities over a wide range of pH and salt concentrations.

A schematic representation of the principle for Capto Core 400 and Capto Core 700 is given in the image below.



The products are based on a highly cross-linked agarose base matrix, which gives good flow properties.

The design of the bead, with an inactive shell that prevents binding of the large target molecules, allows for high resolution at high flow rates and short residence times. Further resin characteristics are found in the table below.

**Table 3.** Characteristics of Capto Core 400 and Capto Core 700

	Capto Core 400	Capto Core 700
<b>Matrix</b>	Highly cross-linked agarose	
<b>Median particle size, <math>d_{50v}</math><sup>1</sup></b>	90 $\mu\text{m}$	85 $\mu\text{m}$
<b>Ligand</b>	Octylamine	
<b>Dynamic binding capacity, <math>Q_{B10}</math><sup>2</sup></b>	~22 mg ovalbumin/mL resin	~13 mg ovalbumin/mL resin
<b>Exclusion limit, <math>M_r</math></b>	~ 400 000	~ 700 000
<b>Recommended operating flow velocity, maximum<sup>3</sup></b>	700 cm/h	500 cm/h
<b>pH stability</b>		
<b>Operational<sup>4</sup></b>	3 to 13	
<b>CIP<sup>5</sup></b>	3 to 14	
<b>Working temperature</b>	4°C to 30°C	
<b>Chemical stability</b>	Stable in commonly used aqueous buffers, 1 M NaOH <sup>6</sup> , 6 M guanidine hydrochloride, 30% isopropanol, 70% ethanol	
<b>Avoid</b>	Oxidizing agents, anionic detergents	

<sup>1</sup> Median particle size of the cumulative volume distribution.

<sup>2</sup> Dynamic binding capacity at 10% breakthrough with a residence time of 3 minutes (1.6 mL/min = 200 cm/h) in HiScreen columns. Buffer: 20 mM Tris-HCl, 0.15 M NaCl, pH 7.5.

<sup>3</sup> At 20 cm bed height, at room temperature using buffers and samples with the same viscosity as water. < 2 bar (0.2 MPa). Recommended operating flow velocity, maximum must be optimized. Start with a lower flow rate to prevent exceeding pressure limits.

<sup>4</sup> pH range where resin can be operated without significant change in function.

<sup>5</sup> pH range where resin can be subjected to cleaning- or sanitization-in-place without significant change in function.

<sup>6</sup> 1 M NaOH should only be used for cleaning purposes.

## 2 Optimization

### General

The aim of designing and optimizing a separation process is to identify conditions that promote binding of the highest amount of impurities, in the shortest possible time, with highest possible product recovery and purity.

Already from the start of the process development it is necessary to consider process cost, cleaning of the resin, and environmental constraints.

The HiScreen column format is suitable for method optimization when developing a new purification process. The small column volume (CV) of 4.7 mL, and the 10 cm bed height allow for scalable experiments at relevant process flow rates. Two columns can be connected in series with a union to give a 20 cm bed height (see [Chapter 5 Scale-up, on page 16](#)).

The HiTrap column format is suitable for initial screening of binding and elution conditions.

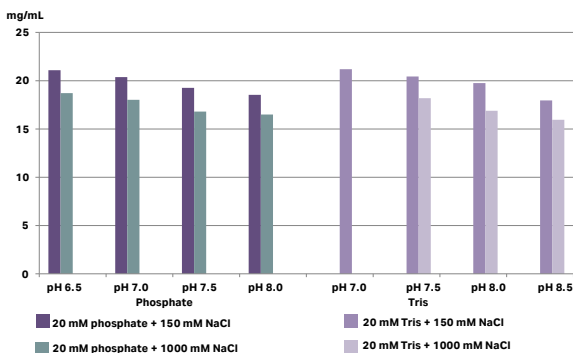
### Buffer and pH

The octylamine ligand is multimodal (both hydrophobic and positively charged), giving a broad window of operation. The function of the ligand has been verified in buffers containing up to 1 M NaCl, see [Fig. 3, on page 8](#).

Capto Core 400 and Capto Core 700 are compatible with most commonly used buffers for size exclusion chromatography and ion exchange chromatography, such as phosphate and Tris. The use of citrate buffer might result in decreased binding capacity.

Most host cell proteins are negatively charged above pH 7. Since the ligands are positively charged, it is advised to use a pH of 7 to 9 to provide good binding of the host cell proteins to the ligands.

DNA is negatively charged over a wider pH range and the efficiency of DNA removal is thus less dependent on pH. To remove DNA with Capto Core 400 and Capto Core 700, it is strongly recommended to treat the sample with a DNase/RNase mixture first, see next section.



**Fig 3.** Static binding capacity (SBC) of Capto Core 700 for ovalbumin in sodium phosphate and Tris buffers with different NaCl concentrations and pH values.

## Reduction of DNA and RNA

In some cases, high levels of DNA and RNA can affect the performance of the Capto Core resins. It is therefore recommended to reduce DNA and RNA levels before the purification step, for example, by using an anion exchange step, or by treatment with a DNase/RNase mixture.



The DNase/RNase mixture degrades all nucleic acids down to oligonucleotides of approximately 3 to 5 base pairs in length. These small fragments and the DNase/RNase mixture itself can enter the core of the resins where they are bound to the octylamine ligands.

## Sample load

Capto Core 400 and Capto Core 700 work as scavengers, meaning that the target passes in the flow-through fraction while the impurities bind to the interior of the beads.

To determine the maximum sample volume that can be loaded onto the column while maintaining the desired purity of target, it is recommended to collect fractions during sample application and analyze the fractions for recovery and purity of target. Typically, 5 to 20 column volumes can be loaded but this is highly dependent on sample composition.

Make sure to start with a high load during method optimization and determine the correct load by analysis of the fractions. After washing the column with binding buffer, the bound impurities will be removed during the cleaning-in-place (CIP) of the resin, see [Chapter 4 Cleaning-In-Place \(CIP\)](#), on page 15.

## 3 Operation

### Prepare the sample

Step	Action
------	--------

- |   |   |
|---|---|
| 1 | Adjust the sample to the composition of the binding buffer, using one of the following methods: <ul style="list-style-type: none"><li>• Dilute the sample with binding buffer.</li><li>• Exchange the buffer using a prepacked column for desalting, see the table below.</li></ul> |
| 2 | Filter the sample through a 0.45 µm filter or centrifuge immediately before loading it to the column. This prevents clogging and increases the longevity of the column when loading large sample volumes.   |

**Note:**

*Water and chemicals used for buffer preparation must be of high purity. Filter buffers through a 0.22 µm or a 0.45 µm filter before use.*

## Prepacked columns for desalting and buffer exchange

**Table 4.** Prepacked columns for desalting and buffer exchange

Column	Loading volume	Elution volume
HiPrep™ 26/10 Desalting <sup>1</sup>	2.5 to 15 mL	7.5 to 20 mL
HiTrap Desalting <sup>2</sup>	0.25 to 1.5 mL	1.0 to 2.0 mL
PD-10 Desalting <sup>3</sup>	1.0 to 2.5 mL <sup>4</sup> 1.75 to 2.5 mL <sup>5</sup>	3.5 mL Up to 2.5 mL
PD MiniTrap™ G-25	0.1 to 2.5 mL <sup>4</sup> 0.2 to 0.5 mL <sup>5</sup>	1.0 mL Up to 0.5 mL
PD MidiTrap™ G-25	0.5 to 1 mL <sup>4</sup> 0.75 to 1 mL <sup>5</sup>	1.5 mL Up to 1 mL

<sup>1</sup> Prepacked with Sephadex™ G-25 Fine and requires a pump or a chromatography system to run.

<sup>2</sup> Prepacked with Sephadex G-25 Superfine and requires a syringe or pump to run.

<sup>3</sup> Prepacked with Sephadex G-25 and can be run by the gravity flow or centrifugation.

<sup>4</sup> Volumes with gravity elution.

<sup>5</sup> Volumes with centrifugation.

## Recommended flow rates

To allow proteins and fragmented DNA to bind properly, the flow rate during sample application must not be too high. A good starting point is 1.2 mL/min for HiScreen and 1.0 mL/min for HiTrap (150 cm/h). During column equilibration and wash steps, higher flow rates can be used, see tables below.

## Flow rates for Capto Core 400

**Table 5.** Recommended flow rates for different operations

Type of operation	Flow rate (mL/min)		Flow velocity (cm/h)	
	HiScreen	HiTrap	HiScreen	HiTrap
Equilibration <sup>1</sup>	up to 5.4	up to 4.0	up to 700	up to 600
Wash <sup>1</sup>	up to 5.4	up to 4.0	up to 700	up to 600
Load of sample	0.8 to 2.3	0.6 to 1.9	100 to 300	100 to 300
Cleaning-in-place <sup>2</sup>	0.16 to 0.31	0.13 to 0.26	20 to 40	20 to 40

<sup>1</sup> Stated flow rates are for buffers with the same viscosity as water at 20°C. For solutions with higher viscosities, for example, 20% ethanol, use lower flow rates.

<sup>2</sup> See also [Chapter 4 Cleaning-In-Place \(CIP\)](#), on page 15.

## Flow rates for Capto Core 700

**Table 6.** Recommended flow rates for different operations

Type of operation	Flow rate (mL/min)		Flow velocity (cm/h)	
	HiScreen	HiTrap	HiScreen	HiTrap
Equilibration <sup>1</sup>	up to 3.9	up to 3.2	up to 500	up to 500
Wash <sup>1</sup>	up to 3.9	up to 3.2	up to 500	up to 500
Load of sample	0.8 to 2.3	0.6 to 1.9	100 to 300	100 to 300
Cleaning-in-place <sup>2</sup>	0.16 to 0.31	0.13 to 0.26	20 to 40	20 to 40

<sup>1</sup> Stated flow rates are for buffers with the same viscosity as water at 20°C. For solutions with higher viscosities, for example, 20% ethanol, use lower flow rates.

<sup>2</sup> See also [Chapter 4 Cleaning-In-Place \(CIP\)](#), on page 15.

## Purification

**Flow rate:** See the preceding two sections for the recommended flow rates for different operations. Collect fractions throughout the separation.

**Column tubing:** Choose the optimal tubing kit (i.d. 0.25, 0.50, or 0.75 mm) for the column and for the application that you intend to use. Tubing with a wider inner diameter gives broader peaks, whereas tubing with a smaller inner diameter gives a higher back pressure.

Step	Action
1	Before connecting the column to a system: <ul style="list-style-type: none"><li>a. Remove the two stoppers from the HiScreen column. Or:</li><li>b. Remove the stopper from the inlet and the snap-off end at the HiTrap column outlet.</li></ul>
2	Connect the column to the system with fingertight 1/16" connectors (28401081).  <b>Note:</b> <i>Make a drop-to-drop connection to prevent air from entering the column.</i>  <b>Note:</b> <i>Make sure that the connectors are tight to prevent leakage.</i>

Step	Action
------	--------

- |   |  |
|---|--|
| 3 | Wash with 1 CV of distilled water to remove the ethanol. This prevents precipitation of buffer salts at exposure to ethanol. |
|---|--|

**Note:**

*The viscosity for 20% ethanol is higher than that for water. For this step, do not use a higher flow rate than mentioned below:*

- 1.9 mL/min for HiScreen (250 cm/h)
- 1.6 mL/min for HiTrap (250 cm/h)

- |   |  |
|---|--|
| 4 | Equilibrate the column with at least 5 CV binding buffer or until the UV baseline, eluent pH, and conductivity are stable. |
|---|--|

- |   |  |
|---|--|
| 5 | Adjust the sample to the chosen starting pH and conductivity, see <a href="#">Prepare the sample, on page 10</a> . |
|---|--|

- |   |   |
|---|---|
| 6 | Load approximately 5 to 20 CV of the sample onto the column. Collect fractions for analysis of purity and recovery of target, see <a href="#">Chapter 2 Optimization, on page 7</a> . |
|---|---|

- |   |  |
|---|--|
| 7 | Wash with 5 to 10 CV binding buffer or until the UV trace of the flowthrough returns to near baseline. |
|---|--|

**Note:**

*The target is in the flowthrough.*

- |   |  |
|---|--|
| 8 | Perform a CIP to elute and clean the column, see <a href="#">Chapter 4 Cleaning-In-Place (CIP), on page 15</a> . |
|---|--|

Step	Action
------	--------

- |   |  |
|---|--|
| 9 | Equilibrate with 5 to 10 CV binding buffer or until the UV baseline, pH, and conductivity reach the required values. |
|---|--|

**Note:**

*Equilibration can be shortened by washing with a high concentration buffer first to roughly obtain the desired pH value. Then wash with binding buffer until the conductivity and pH values are stable.*

**Note:**

*Do not exceed the maximum flow rate (see [Table 5, on page 12](#) and [Table 6, on page 12](#)) or the column hardware pressure limit (see [Table 1, on page 3](#) and [Table 2, on page 4](#)).*

## 4 Cleaning-In-Place (CIP)

### General description

Regular CIP prevents the buildup of contaminants and helps maintaining the capacity, flow properties, and performance of the chromatography resin and prepacked columns.

It is recommended to perform a CIP:

- after each run with live feed
- when an increase in back pressure is seen
- if reduced column performance is observed
- before first-time use or after long term storage

## CIP protocol

For increased contact time and due to the viscosity of the CIP solutions, it is recommended to use a lower flow rate than during purification, see [Table 5, on page 12](#) and [Table 6, on page 12](#).

Wash with a solution of 1 M NaOH in 30% isopropanol or in 27% 1-propanol with reversed flow direction.

Use a total contact time of 30 to 60 minutes, depending on the sample. For more effective cleaning, a pause of 15 to 30 minutes can be included.

The characteristics of the sample determine the final CIP protocol, so the protocol might require optimization. The main parameters to vary during CIP optimization are:

- NaOH and solvent concentrations
- contact time and volume of the CIP solution

## 5 Scale-up

After method optimization at laboratory scale, the process is ready for scale-up. For quick and small scale-up, two HiTrap or HiScreen columns can be connected in series with a union (18112093).

**Note:** *The back pressure increases when columns are connected in series. This can be addressed by decreasing the flow rate.*

Factors such as clearance of critical impurities might change when column bed height is modified and must be validated using the final bed height.



Scale-up is typically performed by keeping bed height and flow velocity (cm/h) constant while increasing bed diameter and flow rate (mL/min or L/h).

For further scale-up, bulk resins are available, see [Chapter 9 Ordering information, on page 22](#).

## Procedure

Step	Action
1	Select bed volume according to required sample load. Keep sample concentration constant.
2	Select column diameter to obtain the desired bed height. The excellent rigidity of the high flow base matrix allows for flexibility in choice of bed heights.
3	Large equipment used during scale-up can cause deviations from the method optimized at small-scale. In such cases, check the buffer delivery and monitoring systems for time delays or volume changes.

## 6 Adjusting pressure limits

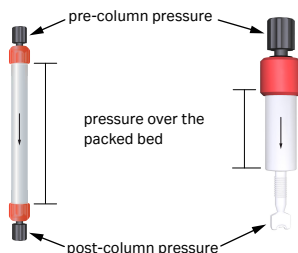
### Introduction

The pressure generated by the flow through a column affects the packed bed and the column hardware, see the following image. Increased pressure is generated when running the column using one or more of the following:

- high flow rates
- high-viscosity buffers or samples
- low temperatures

- a flow restrictor
- long and narrow tubing

**Note:** Exceeding the flow limit can damage the column. See column-specific limit in [Recommended flow rates, on page 11](#).



## ÄKTA avant and ÄKTA pure chromatography systems

The system monitors the pressures automatically. The following table describes which pressures are monitored by each system.

System	Pressures monitored
ÄKTA pure™ without column valve <b>V9-C</b>	<ul style="list-style-type: none"> <li>• system pressure</li> <li>• pre-column pressure</li> </ul>
ÄKTA™ avant and ÄKTA pure with column valve <b>V9-C</b>	<ul style="list-style-type: none"> <li>• system pressure</li> <li>• pre-column pressure</li> <li>• pressure over the packed bed, <math>\Delta p</math></li> </ul>

The pre-column pressure limit is the column hardware pressure limit. The limits are described in [Chapter 1 Product description, on page 3](#).

The maximum pressure for the packed bed depends on resin characteristics and sample or liquid viscosity. The measured value also depends on the tubing that is used to connect the column to the system.

## Systems without multiple pressure sensors

Systems without multiple pressure sensors only measure the system pressure. For optimal system functionality, adjust the pressure limit in the software as follows:

Step	Action
1	<p><b>a.</b> Replace the column with either a piece of tubing with a large inner diameter or a connector with zero dead-volume. Keep all tubing connected to the instrument, including the tubing running to and from the column.</p> <p><b>b.</b> Run the pump at the maximum intended flow rate.</p> <p><b>c.</b> Record the pressure as total system pressure <b>P1</b>.</p> <p><b>Note:</b> <i>The actual pressure over the packed bed (<math>\Delta p</math>) during a run is equal to the measured pressure minus the total system pressure, <b>P1</b>.</i></p>
2	<p><b>a.</b> Disconnect the tubing and run the pump at the maximum intended flow rate.</p> <p><b>Note:</b> <i>The column valve will drip.</i></p>

Step	Action
	<b>b.</b> Record the pressure as <b>P2</b> .
<b>3</b>	<b>a.</b> Calculate the new pressure limit as the sum of <b>P2</b> and the column hardware pressure limit. See <a href="#">Chapter 1 Product description, on page 3</a> for the hardware pressure limit for the given column.
	<b>b.</b> Replace the pressure limit in the software with the calculated value.

**Note:** Repeat the procedure each time parameters are changed.

## 7 Storage

Store HiScreen and HiTrap columns equilibrated with 5 to 10 CV 20% ethanol at the following temperatures:

Product	Storage temperature
HiScreen Capto Core 400	2°C to 8°C
HiScreen Capto Core 700	
HiTrap Capto Core 400	4°C to 30°C
HiTrap Capto Core 700	

Do not freeze.

Make sure that the columns are sealed tightly, to prevent them from drying out.

## 8 Troubleshooting

Problem	Possible cause	Corrective action
High back pressure during the run	Solutions with high viscosity are used.	Decrease the flow rate.
	The column is clogged.	Clean the column, see <a href="#">Chapter 4 Cleaning-In-Place (CIP)</a> , on page 15.
Unstable pressure curve during sample loading	Air bubbles trapped in the sample pump.	If possible, de-gas the sample using a vacuum de-gasser.
Gradual broadening of the eluate peak	Insufficient elution and CIP, caused by contaminant buildup in the column.	Optimize the elution conditions, the CIP protocol, or perform CIP more frequently.
Gradual decrease in yield	Insufficient elution and CIP.	Optimize the elution conditions, the CIP protocol, or perform CIP more frequently.
Gradual increase in CIP peaks	Insufficient elution and CIP.	Optimize the elution conditions, the CIP protocol, or perform CIP more frequently.
High back pressure during CIP	Protein precipitation in the column.	Make sure that all components are eluted from the column.  Optimize elution conditions, or perform a CIP at a lower flow rate.
Decreased column performance despite optimized elution and CIP	Column longevity, which mainly depends on the sample type and sample preparation.	Use a new column.

## 9 Ordering information

Product	Quantity	Product code
HiScreen Capto Core 400	1 × 4.7 mL	17372410
HiScreen Capto Core 700	1 × 4.7 mL	17548115
HiTrap Capto Core 400	5 × 1 mL	17372411
HiTrap Capto Core 700	5 × 1 mL	17548151

Related product	Quantity	Product code
Capto Core 400	25 mL	17372401
	100 mL	17372402
	1 L	17372403
Capto Core 700	25 mL	17548101
	100 mL	17548102
	1 L	17548103
	5 L <sup>1</sup>	17548104

<sup>1</sup> Process-scale quantities are available. Visit [cytiva.com/bioprocess](https://www.cytiva.com/bioprocess), or contact your local Cytiva representative.

Accessories HiScreen	Quantity	Product code
Fingertight connector 1/16" male, narrow	8	28401081
<i>(For connecting HiTrap and HiPrep columns with 1/16" fittings to ÄKTA systems)</i>		
Union 1/16" male - male i.d. 0.5 mm	2	28954326
<i>(For connecting two columns with 1/16" fittings in series)</i>		
Fingertight stop plug, 1/16" <sup>1</sup>	5	11000355
<i>(For sealing a HiScreen column)</i>		

<sup>1</sup> One fingertight stop plug is connected to the inlet and the outlet of each HiScreen column at delivery.

Accessories HiTrap	Quantity	Product code
Connector 1/16" male/luer female <i>(For connecting syringe to top of HiTrap column)</i>	2	18111251
Tubing connector flangeless/M6 female <i>(For connecting tubing to bottom of HiTrap column)</i>	2	18100368
Tubing connector flangeless/M6 male <i>(For connecting tubing to top of HiTrap column)</i>	2	18101798
Union 1/16" female/M6 male <i>(For connecting to original FPLC System through bottom of HiTrap column)</i>	6	18111257

## 10 Related literature on the web

User documentation and other literature related to the products, such as application notes, selection guides, handbooks, and data files can be downloaded from the web. Follow the steps below to access the documentation.

Step	Action
------	--------

- |   |   |
|---|---|
| 1 | Go to the product web page on <a href="https://www.cytiva.com">cytiva.com</a> . |
| 2 | Navigate to <b>Documents</b> .  |
| 3 | Select the type of document and download the chosen literature.                 |

**Give feedback on this document**

Visit [cytiva.com/techdocfeedback](https://cytiva.com/techdocfeedback) or scan the QR code.



## cytiva.com

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corporation or an affiliate doing business as Cytiva.

ÄKTA, ÄKTA pure, Capto, HiPrep, HiScreen, HiTrap, MidiTrap, MiniTrap, and Sephadex are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Any other third-party trademarks are the property of their respective owners.

© 2020–2025 Cytiva

For local office contact information, visit [cytiva.com/contact](https://cytiva.com/contact)

28995881 AK V:8 03/2025