



Capto™ MMC

Multimodal chromatography resin

Instructions for Use

Capto™ MMC is a multimodal salt-tolerant BioProcess™ resin for capture and intermediate purification of proteins from large feed volumes by packed bed chromatography.

Capto MMC increases productivity and reduces cost with:

- high dynamic binding capacity at high conductivity
- high volume throughput
- smaller unit operations

Table of Contents

1	BioProcess resins	3
2	Resin properties	3
3	Method design and optimization	6
4	Screening of binding and elution conditions	8
5	Optimization of throughput	8
6	Scaling up	13
7	Packing columns	14
8	Evaluation of column packing	28
9	Maintenance	31
10	Ordering information	34

Important

Read these instructions carefully before using the product.

Safety

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

1 BioProcess resins

BioProcess chromatography resins are developed and supported for production-scale chromatography. All BioProcess resins are produced with validated methods and are tested to meet manufacturing requirements. Secure ordering and delivery routines give a reliable supply of resins for production scale. Regulatory Support Files (RSF) are available for BioProcess resins to assist in process validation and submissions to regulatory authorities. BioProcess resins cover all purification steps from capture to polishing.

2 Resin properties

Capto MMC has a ligand with multimodal functionality (Fig 1). The multimodal functionality gives a different selectivity compared to traditional ion exchangers and also provides the possibility of binding proteins at high salt conditions.

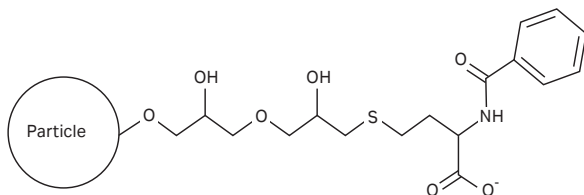


Fig 1. The Capto MMC ligand exhibits many functionalities for interaction with a target molecule. The most pronounced are ionic interaction, hydrogen bonding, and hydrophobic interaction.

Capto MMC is designed to increase speed and throughput in capture and intermediate purification. By offering high capacity at high salt concentrations, and high flow velocities with low back pressure, process cycle times can be reduced and productivity increased.

The resin is based on a high-flow agarose matrix. Typical flow velocities at large scale (1 m column diameter and 20 cm bed height) are 600 cm/h or over, with a back pressure below 3 bar (Fig 2). The highly cross-linked agarose base matrix gives the resin high chemical and physical stability. Characteristics such as capacity, elution behavior, and pressure-flow properties are unaffected by the solutions commonly used in process chromatography (Table 1).

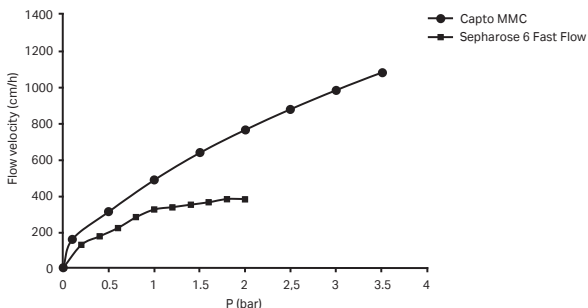


Fig 2. Pressure-flow properties for Capto MMC compared to Sepharose™ 6 Fast Flow. Running conditions: BPG 300 (30 cm i.d.), open bed at settled bed height equal to 20 cm, with water at 20°C.

Table 1. Characteristics of Capto MMC

Matrix	Highly cross-linked agarose, spherical
Functional group	Multimodal weak cation exchanger
Ionic capacity	0.07 to 0.09 mmol H ⁺ /mL resin
Particle size, d_{50v} ¹	~ 75 µm
Pressure-flow characteristics	≤ 600 cm/h at ≤ 0.3 MPa in a 1 m diameter column and 20 cm bed height (at 20°C using process buffers with the same viscosity as water) ²
Dynamic binding capacity, QB_{10%} ³	≥ 45 mg BSA/mL resin at 30 mS/cm
pH stability, operational ⁴	3 to 12
pH stability, CIP ⁵	3 to 14
Working temperature ⁶	4°C to 30°C
Chemical stability	Stable in commonly used aqueous buffers, 1 M acetic acid, 1.0 M NaOH ⁷ , 8 M urea, 6 M guanidine hydrochloride, 70% ethanol
Avoid	Oxidizing agents, cationic detergents
Autoclavability	17 min at 121°C in 0.05 M phosphate buffer, pH 7, 10 cycles

¹ Median particle size of the cumulative volume distribution.

² The pressure-flow characteristics describe the relationship between pressure and flow under the set circumstances. The pressure given shall not be taken as the maximum pressure of the resin.

³ Dynamic binding capacity at 10% breakthrough by frontal analysis at a mobile phase velocity of 300 cm/h in a Tricorn™ 5/100 column at 10 cm bed height (2 min residence time) for BSA in 50 mM sodium acetate, pH 4.75, 250 mM NaCl.

⁴ pH range where resin can be operated without significant change in function.

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

⁶ Capto MMC can be used under cold-room conditions, but for some proteins the capacity can decrease.

⁷ 1.0 M NaOH must only be used for cleaning purposes.

3 Method design and optimization

Aim

The aim of designing and optimizing a method for the separation of biomolecules is to identify conditions that promote binding of the highest amount of target molecule, in the shortest possible time with the highest possible product recovery and purity. For optimization of binding conditions, pH, and conductivity (salt concentration) must be screened. Binding at relatively high conductivity can be expected with Capto MMC, but it depends on the pH. Binding can also be expected at higher pH values compared to traditional ion exchangers.

A pH gradient experiment in which the target molecule is loaded in analytical amounts on Capto MMC and eluted in an increasing pH gradient helps establishing the elution pH. A separate experiment with different salt concentrations in the pH gradient can also be included for better understanding. Knowing the elution pH is useful when setting up the screening experiments for binding and elution. Establishing the elution pH through a gradient run is more reliable than basing the pH selection on the isoelectric point of the target molecule.

Design of Experiments

The fact that Capto MMC allows capture of proteins at high conductivity in many cases limits the use of increasing salt concentrations as an efficient way of eluting proteins. Optimal elution is often achieved by a combination of changes in pH, buffer concentration, and eluting salt. Design of Experiments (DoE) is an effective tool for investigation of the effect of several parameters on protein recovery in order to establish the optimal elution protocol.

Examples of the DoE approach are described in the application note *High-throughput screening for elution conditions on Capto MMC using PreDicator plates (CY13156)*. Alternatively, a stepwise elution optimization protocol can be applied. For rapid screening of binding or elution conditions with low sample consumption, PreDicator™ plates (*Instructions for Use 28925834, Data file CY13663*) are preferably used.

For further reading, refer to the application note *High-throughput screening and process development for capture of recombinant pro-insulin from E. coli (CY13364)* where optimization of a chromatographic capture step for pro-insulin from *E. coli* on Capto MMC is described.

4 Screening of binding and elution conditions

Screening of binding and elution conditions can be done either with traditional column chromatography or in batch format (PreDicator plates). The procedure for the two formats is the same, but by using the latter format screening is quicker and much less sample is consumed, which allows a larger experimental space to be screened. Below, only the plate approach is described, but the procedure is essentially the same for both formats.

5 Optimization of throughput

The dynamic binding capacity (DBC) for the target protein must be determined by frontal analysis using real process feedstock. Because the DBC is a function of the flow velocity applied during sample application, the breakthrough capacity must be defined over a range of different residence times (flow velocities) to show the optimum level of throughput.

Loading conditions

The initial elution pH of the target molecule can be established by performing a pH gradient experiment as described above. PreDicator plates are used to screen for best loading conditions. Vary the salt concentration (typically 0 to 300 mM) and pH (0.5 to 2.5 pH units below the elution pH) and determine the static binding capacity under the different conditions. The screening gives the conditions, that is, the salt concentration and pH at which the capacity is the highest.

Before sample loading, adjust the pH and conductivity of the sample to the target loading conditions. This is done by buffer exchange or by direct adjustment of pH and salt concentration. Buffers normally used for ion exchange chromatography can also be used for Capto MMC (Table 2).

Table 2. Recommended buffers

pH interval	Buffer ¹	Concentrations ²
4 to 5.5	Acetate	20 to 100 mM
3.0 to 6.5	Citrate	20 to 200 mM
5.5 to 6.5	Bis-Tris	20 to 50 mM
6 to 7.5	Phosphate	20 to 200 mM
7.5 to 9.0	Tris	20 to 50 mM
8.8 to 10.6	Glycin-NaOH	20 to 100 mM

¹ The choice of buffer systems and salts can affect performance.

² Conductivity can be adjusted by adding salt or by varying the buffer concentration.

Loading density

The capacity obtained in plates is static binding capacity and not dynamic binding capacity (DBC). The loading density for the elution study must be based on the DBC under optimal binding conditions. How loading density (amount of loaded target molecules per volume resin) is specified varies. Two examples of loading density specifications are:

- 70% of the dynamic binding capacity at 10% breakthrough, and
- 80% of the dynamic binding capacity at 1% breakthrough.

Elution conditions

Screening of elution conditions is done in PreDicator plates in a similar manner as screening of binding conditions. The pH (typically 0.5 to 2.5 pH units above the initial elution pH established earlier) and salt concentration (typically 0 to 1 or 2 M) are screened. The loading density is set based on the DBC (above).

An example of an elution screening study is shown in Figure 3 (*Application note CY13156*). The figure compares the results obtained in PreDicator plates with results obtained in columns. For best elution profile, a rapid change in pH is desirable. Increasing the ionic strength of the buffer will facilitate this. The results with NH_4^+ salt illustrate that other salts than NaCl can be more effective.

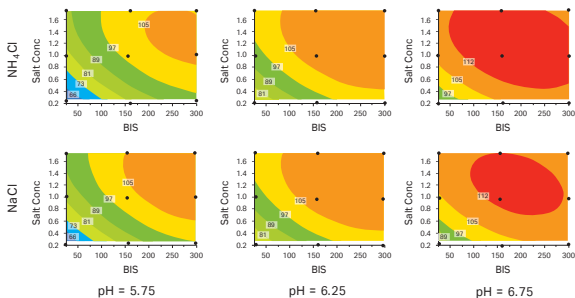
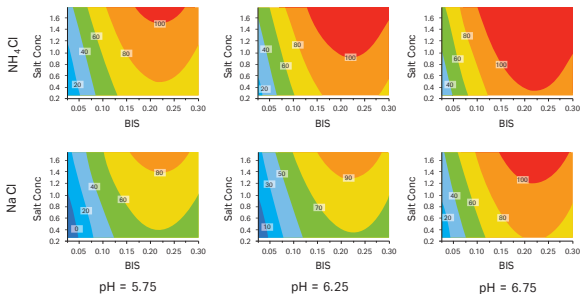
A**B**

Fig 3. Contour plots for the recovery of BSA in A) PreDictor plates and B) Tricorn column (data from Application Note *Optimizing elution conditions on Capto MMC using Design of Experiments (CY14386)*). Recovery is plotted as a function of salt concentration and buffer ionic strength (BIS) at three different pH values for the two salt types NaCl and NH_4Cl . Experimental data points are shown as black dots.

The effect of urea and organic modifiers ethanol (EtOH) and isopropanol (IPA) on dynamic binding capacity is shown in Figure 4. The decreased capacity in the presence of urea and organic modifiers suggests that these can be used to improve elution efficiency.

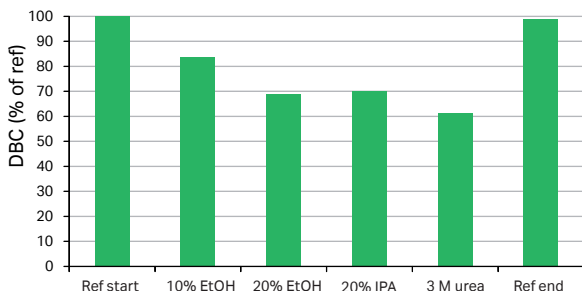


Fig 4. Dynamic binding capacity (DBC) of BSA on Capto MMC. Effect of organic modifiers and urea. DBC is plotted as the percentage of the initial capacity (Ref start).

General purification protocol

- Adjust pH and salt concentration to match optimal conditions found during screening.
- Equilibrate column with loading buffer of the same pH and salt concentration as the sample.
- Apply sample onto the column.
- Wash out unbound material with loading buffer.
- Elute sample under optimal elution conditions established during screening. Gradient (pH or salt) or step-elution can be used.
- Regenerate column to elute bound material.
- Clean-in-place.
- Re-equilibrate.

6 Scaling up

After optimizing the method at laboratory scale, the process can be scaled up, using the following general approach:

1. Select the bed volume according to required binding capacity.
2. Select a column diameter to obtain a bed height of 10 to 45 cm. To maximize the potential of Capto MMC at a large scale, a typical bed height range is 20 cm or higher.
3. Scale-up is typically done by keeping bed height and flow velocity constant, while increasing bed diameter and volumetric flow rate.

General recommendations for scale-up

However, as optimization is done with small column volumes, in order to save sample and buffer, some parameters like the dynamic binding capacity can be optimized using shorter bed heights than those being used at the final scale. As long as the residence time is constant, the binding capacity for the target molecule remains the same.

Other factors, such as clearance of critical impurities, can change when column bed height is changed and must be validated using the final bed height. The residence time is approximated as the bed height (cm) divided by the flow velocity (cm/h) applied during sample loading.

7 Packing columns

Recommended columns

Table 3. Recommended columns for Capto MMC

Column	Inner diameter (mm)	Bed volume ¹	Bed height (cm)
Laboratory scale			
Tricorn 5/100	5	2 mL	10
Tricorn 10/100	10	8 mL	10
HiScale™ 10/40 ²	10	8 to 20 mL	25
HiScale 16/20	16	20 to 40 mL	max 20
HiScale 16/40	16	20 to 70 mL	max 35
HiScale 26/20	26	53 to 106 mL	max 20
HiScale 26/40	26	53 to 186 mL	max 35
HiScale 50/20	50	196 to 393 mL	max 20
HiScale 50/40	50	196 to 687 mL	max 35
Production scale			
AxiChrom™ ³	50 to 200	0.2 to 12.5 L	max 40
AxiChrom ³	300 to 1 000	7 to 314 L	max 40
BPG ^{3,4}	100 to 300	1 to 28 L	max 40
ChromaFlow™ standard ^{3,5}	400 to 800	12 to 151 L	max 30

¹ Bed volume range calculated from 10 cm bed height to maximum bed height.

² Packing methods for bed heights up to 25 cm are provided.

³ Packing instructions for Capto MMC in process-scale columns are described in Application Note CY29524.

⁴ The pressure rating for BPG 450 is too low for Capto MMC resins.

⁵ Larger pack stations might be required at larger diameters.

All large-scale columns can be supplied as variable bed height columns. Do not choose large diameter columns if the bed height is low.

Packing Tricorn 5/100 and 10/100 columns

The following instructions are for packing Tricorn 5/100 and Tricorn 10/100 columns with a 10 cm bed height.

For more details about packing Tricorn columns, refer to the Instructions *Tricorn Empty High Performance Columns* (28409488).

Materials needed

- Capto MMC
- glass filter funnel
- plastic spoon or spatula
- filter flask
- measuring cylinder
- 10 mM NaCl in distilled water

Equipment needed

ÄKTA™ systems, or a stand-alone pump, depending on the flow rate required, can be used for packing. The pump filter unit and the flow restrictor must be removed due to the high flow velocity used in the column packing in order to decrease the system back pressure.

For packing Tricorn 5/100, an additional Tricorn 5/100 tube is used as packing tube which is connected by a Tricorn Packing Connector 5-5.

When packing Tricorn 10/100 column, use a Tricorn Packing Equipment 10/100 that includes the 10-mm packing connector, 100-mm glass tube (to be used as packing tube), and bottom unit with filter holder, cap, and stop plug.

When working with large volumes, real feed or repeated loading, Tricorn coarse filter kits are recommended to reduce the risk of clogging. Use Tricorn 5 Coarse Filter Kit (11001253) or Tricorn 10 Coarse Filter Kit (11001254).

Calculating amount of resin

The amount of resin needed can be calculated by: column cross-sectional area (cm^2) \times bed height (cm) \times compression factor (sedimented resin bed height/packed resin bed height). The compression factor is approximately 1.1 for Capto MMC.

Washing the resin

Equilibrate all materials to room temperature. Capto MMC is delivered in 20% ethanol and must first be washed with the recommended packing solution, 10 mM NaCl. Place the glass filter funnel on the filter flask. Pour the resin into the funnel and wash 5 times with 1 to 2 mL of 10 mM NaCl/mL resin. Gently stir with a spatula between additions.

Preparing the packing slurry

Transfer the washed resin from the funnel to a beaker. Add 10 mM NaCl to obtain a 60% slurry concentration. Measure the slurry concentration after the slurry has settled overnight in a measuring cylinder or use the Slurry Concentration Kit (29096100). Tricorn columns can also be packed with an excess of resin that can be removed during packing (step 13).

Packing procedure

To pack the column, use 10 mM NaCl in distilled water as packing solution and proceed as follows:

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

- | | |
|---|---|
| 1 | Rinse the column and packing tube in 10 mM NaCl. |
| 2 | Insert a bottom filter into the filter holder. |
| 3 | Wet the filter and the O-ring on the filter holder by dipping the filter holder into water, buffer, or 20% ethanol. |
| 4 | Insert the filter holder into the column tube. Make sure that the "keyed" part of the filter holder fits into the slot on the threaded section on the column tube. Screw the end cap onto the column tube. |
| 5 | Screw a suitable Tricorn packing connector onto the top of the column tube. The Tricorn packing connector must be fitted with suitable O-rings (included with the Tricorn packing connector). Screw the Tricorn packing tube into the upper fitting of the Tricorn packing connector. |
| 6 | Attach the column and the packing tube vertically to a lab stand. |
| 7 | Pour some packing solution into the tube and make sure that the liquid drips from the column bottom outlet. Insert a stop plug into the bottom unit when approximately 1 cm of packing solution remains. |

Step	Action
8	Prime the system and the column inlet capillary with packing solution.
9	Fill both column tube and packing tube with slurry. Prevent formation of air bubbles by pouring the slurry along a thin capillary.
10	Attach an extra bottom unit or an adapter unit to the top of the packing tube. Connect the pump to the top of the packing tube and remove the stop plug from the bottom of the column tube.
11	Increase the flow and pack the resin at 540 cm/h (Tricorn 5/100: 1.8 mL/min, Tricorn 10/100: 7.1 mL/min). When the liquid above the resin bed is clear, continue packing for 10 min.
12	Pack the resin for an additional 10 min at 3000 cm/h (Tricorn 5/ 100: 9.8 mL/min, Tricorn 10/100: 39.3 mL/min).
13	Switch off the pump and connect a stop plug into the bottom unit. Remove the packing tube and packing connector (the liquid is allowed to pour down outside the column tube). If necessary, remove excess resin with a Pasteur pipette or spatula by resuspending the top of the packed bed. Make sure that the resin surface is as even as possible.
14	Add packing solution up to the upper edge of the column tube.

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

- | | |
|----|--|
| 15 | Place a pre-wetted filter on top of the fluid in the column. |
|----|--|

Note:

The top coarse filter is inserted using another procedure. Refer to the separate instructions included in Tricorn coarse filter kits.

- | | |
|----|--|
| 16 | Prepare the adapter unit by turning the guiding ring inside the adapter unit down to its end position so that the guiding ring is level with the bottom of the adapter unit. |
|----|--|

- | | |
|----|--|
| 17 | Wet the O-ring on the adapter unit by dipping it in water, packing solution, or 20% ethanol. |
|----|--|

- | | |
|----|--|
| 18 | Screw the guiding ring back 1.5 turns. |
|----|--|

- | | |
|----|--|
| 19 | Attach the top adapter unit onto the column tube, making sure the inner part of the guiding ring fits into the slot on the column tube threads. Make sure that there are no air bubbles. |
|----|--|

Note:

The top adapter must be connected but not fully screwed down.

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

Note:

Although it is possible to fit the adapter unit on the column tube without keying the inner locking device into the slot on the column tube, the adapter lock will not function. The consequence of this is that the adapter is not locked in position and accidental turning of the adapter is possible.

- | | |
|----|--|
| 20 | Connect the pump, remove the stop plug and start a flow (Tricorn 5/100: 1 mL/min, Tricorn 10/100: 5 mL/min). |
| 21 | Slowly screw the adapter unit down until the filter meets the bed surface. Make sure that the filter meets the bed horizontally. |
| 22 | Increase the flow to 3000 cm/h (Tricorn 5/100: 9.8 mL/min, Tricorn 10/100: 39.3 mL/min). |
| 23 | If the resin bed compresses, slowly screw the adapter unit down to the resin surface with maintained flow. |
| 24 | Pack the resin for 5 min. If the bed has compressed further, screw the adapter unit down to the resin surface. |
| 25 | Stop the flow and connect a stop plug to the bottom unit. |

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

- | | |
|----|--|
| 26 | Disconnect the pump. Screw the adapter unit down for a further 2/3 turns. Lock the adapter and attach a stop plug. If the adapter lock is attached correctly, it is not possible to turn the adapter unit. |
|----|--|

Note: *It is recommended to perform CIP after packing, as column packing involves open handling of the resin.*

The column is now ready to be tested.

Testing the packed column

See [Chapter 8 Evaluation of column packing, on page 28](#).

Packing HiScale columns

The following instructions are for packing HiScale 10/40, HiScale 16/20, 16/40, HiScale 26/20, 26/40, and HiScale 50/20, 50/40 columns with 10, 20, and 35 cm bed heights. For HiScale 10/40, packing methods are provided for 20 and 25 cm bed heights.

For more details about packing HiScale columns, refer to the Instructions *HiScale columns (10, 16, 26, 50) and accessories (28967470)*.

Materials needed

- Capto MMC
- HiScale column
- HiScale packing tube (depending on bed height)
- plastic spoon or spatula
- G3 glass filter funnel

- vacuum suction equipment
- filter flask
- measuring cylinder
- 20% ethanol with 0.4 M NaCl

Equipment

ÄKTA systems, or a stand-alone pump, depending on the flow rate required, can be used for packing.

Equilibrate all materials to room temperature.

Definitions

The height of a bed that has settled by gravity differs from the height of a bed that has settled at a given flow velocity (consolidated). Therefore, the compression factor (CF) must be separated from the packing factor (PF). The parameters for calculating the bed height are described below:

L_{settled}	Settled bed height (cm) Bed height measured after settling by gravity
L_{cons}	Consolidated bed height (cm) Bed height measured after settling at a given flow velocity
L_{packed}	Packed bed height (cm)
CF	Compression factor, $CF = L_{\text{settled}} / L_{\text{packed}}$
PF	Packing factor, $PF = L_{\text{cons}} / L_{\text{packed}}$
A_C	Cross-sectional area of the column (cm^2)
V_C	Column volume, $V_C = L_{\text{packed}} \times A_C$ (mL)
C_{slurry}	Slurry concentration (%)

Preparation of the slurry

To measure the slurry concentration, pour the resin in a measuring cylinder and let the resin settle in 20% ethanol at least overnight or use the Slurry Concentration Kit (29096100). This method can also be used for HiScale columns.

Washing the resin

Place a glass filter funnel onto a filter flask. Suspend the resin by shaking it, pour it into the filter funnel, and wash it as follows:

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

- | | |
|---|---|
| 1 | Wash the resin 5 times with 1 to 2 mL 20% ethanol with 0.4 M NaCl/mL resin. |
| 2 | Gently stir with a spatula between additions. |
| 3 | Transfer the washed resin from the funnel to a beaker and add 20% ethanol with 0.4 M NaCl to obtain the recommended slurry concentration. |

Main packing parameters

Table 4. Main packing parameters for HiScale 10/40

Parameter	HiScale 10/40	
Bed height (cm)	20	25
Slurry/packing solution	20% ethanol with 0.4 M NaCl	
Slurry concentration (%)	50	50
Packing factor (PF)	1.025	1.02
Settling flow velocity (cm/h)	76	76
Settling flow rate (mL/min)	1	1
Packing flow velocity (mL/min)	1910	1910
Packing flow rate (mL/min)	25	25
Conditioning flow velocity (cm/h)	1910	1910
Conditioning flow rate (mL/min)	25	25

Table 5. Main packing parameters for HiScale 16/20 and HiScale 16/40

Parameter	HiScale 16/20	HiScale 16/40	
Bed height (cm)	10	20	35
Slurry/packing solution	20% ethanol with 0.4 M NaCl		
Slurry concentration (%)	50	50	50
Packing factor (PF)	1.10	1.10	1.02
Packing flow velocity (cm/h)	750	750	750
Packing flow rate (mL/min)	20.1	16.7	11.7
Conditioning flow velocity (cm/h)	750	750	420
Conditioning flow rate (mL/min)	25	25	14

Table 6. Main packing parameters for HiScale 26/20 and HiScale 26/40

Parameter	HiScale 26/20	HiScale 26/40	
Bed height (cm)	10	20	35
Slurry/packing solution	20% ethanol with 0.4 M NaCl		
Slurry concentration (%)	50	50	50
Packing factor (PF)	1.15	1.10	1.03
Packing flow velocity (cm/h)	750	750	750
Packing flow rate (mL/min)	66	66	66
Conditioning flow velocity (cm/h)	750	750	420
Conditioning flow rate (mL/min)	66	66	37

Table 7. Main packing parameters for HiScale 50/20 and HiScale 50/40

Parameter	HiScale 50/20	HiScale 50/40	
Bed height (cm)	10	20	35
Slurry/packing solution	20% ethanol with 0.4 M NaCl		
Slurry concentration (%)	50	50	50
Packing factor (PF)	1.15	1.15	1.03
Packing flow velocity (cm/h)	750	750	750
Packing flow rate (mL/min)	250	250	250
Conditioning flow velocity (cm/h)	750	750	420
Conditioning flow rate (mL/min)	250	250	140

Packing procedure

Step Action

- 1 Assemble the column according to the column instructions *HiScale columns (10, 16, 26, 50) and accessories (28967470)*.

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

- | | |
|---|---|
| 2 | Attach the column tube to a laboratory stand. |
| 3 | Connect the bottom adapter unit to the pump or a syringe and prime the bottom net with a slow flow of packing solution. This is easiest done if the nets are dry but if air is trapped under the net it can be removed by a light suction with a syringe. |
| 4 | Attach the bottom adapter unit in the bottom of the column tube and tighten the O-ring. |
| 5 | Fill the column with approximately 1 cm packing liquid using the pump/syringe. Disconnect the pump/syringe and put a stop plug on the outlet. |
| 6 | Place the packing tube on top of the column tube. |
| 7 | Connect the top adapter to the pump and prime it with a slow downward flow. The net needs to be facing the roof as this is done. If air is trapped under the net it can be removed by a light suction with a syringe. |
| 8 | Fill the column with slurry suspended in packing solution. If needed, top up the slurry with extra packing solution so the top adapter dips into the slurry to avoid air under the net. |
| 9 | Attach the top adapter unit on top of the packing tube. Tighten the O-ring firmly and remove the bottom stop plug. |

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

10	Start a downward flow with settling (HiScale 10/40) or packing flow velocity according to Tables 4, 5, 6, and 7 .
----	---

11	Let the flow run until the bed has consolidated.
----	--

Note:

For HiScale 10/40 columns, increase the flow to packing velocity according to [Table 4, on page 24](#). Let the flow run for 4 minutes.

12	Use the graduation marks on the column to measure the bed height. There might be a buildup of resin at the column wall after the bed is consolidated. To easier see where the top of the bed is, a light source can be used.
----	--

13	Calculate the final bed height by dividing the consolidated bed height with the desired packing factor.
----	---

$$L_{\text{packed}} = L_{\text{cons}} / \text{PF}$$

14	Turn off the flow and attach a stop plug in the bottom.
----	---

15	Dismount the top adapter from the packing tube.
----	---

16	Over a beaker or a sink, detach the packing tube from the column.
----	---

17	Remount the top adapter in the column tube. Make sure that no air is trapped under the net and lower the adapter down to 1 to 2 cm above the bed, making sure that the surface is not disturbed.
----	--

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

- | | |
|----|--|
| 18 | Tighten the O-ring on the adapter. Remove the bottom stop plug and carefully start turning the end cap down. While spilling out liquid through the bottom, proceed turning until the calculated final bed height is reached. |
| 19 | Make sure that the pressure peaks that occur during turning the end knob down do not exceed the pressure specifications of the resin. |
| 20 | Start a downward flow to flow condition the bed. The flow rate is shown in Tables 4, 5, 6, and 7 . |
| 21 | Let the flow run for about 10 column volumes. |

Note: *It is recommended to perform CIP after packing, as column packing involves open handling of the resin.*

The column is ready to be tested.

Testing the packed column

See [Chapter 8 Evaluation of column packing, on page 28](#).

8 Evaluation of column packing

Intervals

Test the column efficiency to check the quality of packing. Testing should be done after packing, at regular intervals during the working life of the column, and when separation performance is seen to deteriorate.

Column efficiency testing

The best method of expressing the efficiency of a packed column is in terms of the height equivalent to a theoretical plate (HETP) and the asymmetry factor (A_s). These values are easily determined by applying a sample such as 1% acetone solution to the column. Sodium chloride can also be used as a test substance. Use a concentration of 0.8 M NaCl in water as sample and 0.4 M NaCl in water as eluent.

For more information about column efficiency testing, consult the application note *Column efficiency testing (CY13149)*.

Note: *The calculated plate number varies according to the test conditions and must only be used as a reference value. It is important that test conditions and equipment are kept constant so that results are comparable. Changes of solute, solvent, eluent, sample volume, flow velocity, liquid pathway, temperature, etc., affect the results.*

Sample volume and flow velocity

For optimal results, the sample volume must be at maximum 2.5% of the column volume and the liquid velocity 30 cm/h. If an acceptance limit is defined in relation to column performance, the column plate number can be used as one of the acceptance criteria for column use.

Method for measuring HETP and A_s

Calculate HETP and A_s from the UV curve (or conductivity curve) as follows:

$$\text{HETP} = \frac{L}{N}$$

L = bed height (cm)

N = number of theoretical plates

$$N = 5.54 \times \left(\frac{V_R}{W_h} \right)^2$$

V_R = volume eluted from the start of sample application to the peak maximum

W_h = peak width measured as the width of the recorded peak at half of the peak height

V_R and W_h are in the same units

The concept of reduced plate height is often used for comparing column performance.

The reduced plate height (h) is calculated as follows:

$$h = \frac{HETP}{d_{50v}}$$

d_{50v} = median particle size of the cumulative volume distribution

As a guideline, a value of < 3 for (h) is very good.

The peak should be symmetrical, and the asymmetry factor as close to 1 as possible (a typical acceptable range could be $0.8 < A_s < 1.8$).

A change in the shape of the peak is usually the first indication of bed deterioration due to excessive use.

Peak asymmetry factor calculation:

$$A_s = \frac{b}{a}$$

a = ascending part of the peak width at 10% of peak height

b = descending part of the peak width at 10% of peak height

Figure 5 shows a UV trace for acetone in a typical test chromatogram from which the HETP and A_s values are calculated.

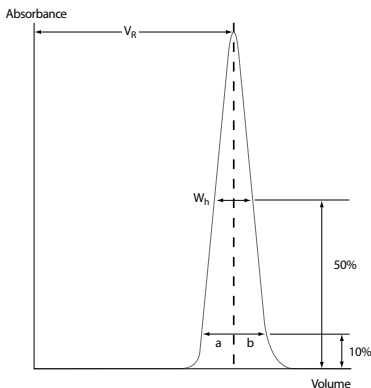


Fig 5. A typical test chromatogram showing the parameters used for HETP and A_s calculations.

9 Maintenance

For the best performance from Capto MMC over a long working life time, follow the procedures described below.

Equilibration

After packing, and before a chromatographic run, equilibrate with start buffer by washing with at least 5 bed volumes, or until the column effluent shows stable conductivity and pH values.

Regeneration

After each separation, elute any reversibly bound material with a high ionic strength solution (e.g., 2 M NaCl in buffer) and at the same time increase pH to about 10 to 11. Regenerate the resin by washing with reversed flow direction with at least 5 bed volumes of buffer, or until the column effluent shows stable conductivity and pH values.

Cleaning-In-Place (CIP)

CIP is a procedure that removes impurities or contaminants such as lipids, endotoxins, and precipitated or denatured proteins that remain in the packed column after regeneration. These types of impurity or contamination occur frequently when working with crude sample. Regular CIP prevents the buildup of impurities or contaminants in the resin bed and helps to maintain the capacity, flow properties, and general performance of the resin. A specific CIP protocol must be designed for each process according to the type of impurities or contaminants present.

It is recommended to perform a CIP:

- before first-time use – especially after packing the column – or after long-term storage
- between cycles
- when an increase in back pressure or a reduction in column performance is observed
- to prevent potential cross-contamination or carry-over, when the same column is used for purification of different proteins or protein lots and batches.

Recommended protocols for removing specific contaminants or impurities are described below.

CIP protocols

Precipitated, hydrophobically bound proteins or lipoproteins	Wash with at least 3 column volumes of 1.0 M NaOH at 40 cm/h with reversed flow direction. Contact time 1 to 2 hours, depending on feed.
Ionically bound proteins	Wash with 0.5 to 2 column volumes of 2 M NaCl with reversed flow direction. Contact time 10 to 15 min.
Lipids and very hydrophobic proteins	Wash with 2 to 4 column volumes of up to 70% ethanol, 30% isopropanol ¹ , or 1% to 5% 1-propanol with reversed flow direction. Contact time 1 to 2 hours, depending on feed. 1-propanol has a higher flash point and might be preferred in an industrial environment. Alternatively, wash with 2 to 4 column volumes of 0.1% nonionic detergent with reversed flow direction. Contact time 1 to 2 hours, depending on feed.

¹ Consult local safety regulations to verify if safety precautions are required for the use of 70% ethanol or 30% isopropanol.

Sanitization

To reduce microbial contamination of the packed column, sanitization using 0.5 to 1.0 M NaOH with a contact time of 1 hour is recommended. The CIP protocol for precipitated, hydrophobically bound proteins or lipoproteins removes bound contaminants and sanitizes the resin effectively. For more information about the use of NaOH for sanitization, refer to *Application note (CY13951)*.

Storage

Store unused resin in its container at a temperature of 4°C to 30°C. Make sure that the screw top is fully tightened. Packed columns must be equilibrated in 20% ethanol to prevent microbial growth. After storage, equilibrate with at least 5 bed volumes of starting buffer before use.

10 Ordering information

Product	Pack size	Product code
Capto MMC	25 mL	17531710
	100 mL	17531702
	1 L	17531703
	5 L ¹	17531704
	10 L ¹	17531705
	60 L ¹	17531760

¹ Pack sizes available upon request.

All bulk resins products are supplied in suspension in 20% ethanol. For additional information, including data file, contact your local Cytiva representative.

Related product	Pack size	Product code
HiTrap™ Capto MMC	1 × 1 mL	29400465
	5 × 1 mL	11003273
	5 × 5 mL	11003275
HiTrap Capto IEX Selection kit	5 × 1 mL	28934388
PreDictor Capto MMC, 6 µL	4 x 96-well filter plates	28925814
PreDictor Capto MMC, 20 µL	4 x 96-well filter plates	28925815

Related product	Pack size	Product code
PreDictor Capto MMC, 50 µL	4 x 96-well filter plates	28925816
PreDictor Capto MMC, 2, 4, 6, 8, 20, and 50 µL	4 x 96-well filter plates	28943281
PreDictor RoboColumn Capto MMC, 200 µL	one row of eight columns	28986084
PreDictor RoboColumn Capto MMC, 600 µL	one row of eight columns	28986178
HiScreen™ Capto MMC	1 × 4.7 mL	28926980
ReadyToProcess™ Capto MMC	1 L (80/200)	28951118
	2.5 L (126/200) ¹	28929120
	5 L (178/200) ¹	29146145
	10 L (251/200) ¹	28929121
	12.4 L (251/200) ¹	29589044
	15 L (359/200) ¹	29372600
	20 L (359/200) ¹	28929122
	57 L (600/200) ¹	29428374
Tricorn 5/20 column	1	28406408
Tricorn 5/50 column	1	28406409
Tricorn 5/100 column	1	28406410
Tricorn 5/150 column	1	28406411
Tricorn 5/200 column	1	28406412
Tricorn 10/20 column	1	28406413
Tricorn 10/50 column	1	28406414
Tricorn 10/100 column	1	28406415
Tricorn 10/150 column	1	28406416
Tricorn 10/200 column	1	28406417
Tricorn 10/300 column	1	28406418
Tricorn 10/600 column	1	28406419
HiScale 10/40	1	29360550
HiScale 16/20	1	28964441

Related product	Pack size	Product code
HiScale 16/40	1	28964424
HiScale 26/20	1	28964514
HiScale 26/40	1	28964513
HiScale 50/20	1	28964445
HiScale 50/40	1	28964444

¹ Pack sizes on request.

Related literature		Reference
Handbook	Multimodal chromatography	CY14738
Data files	Capto MMC	CY13468
	HiScreen preppacked columns	CY13473
	PreDicator 96-well filter plates and Assist Software	CY13663
Application notes	High-throughput screening and process development for capture of recombinant pro-insulin from <i>E. coli</i>	CY13364
	Efficient purification of the pertussis antigens toxin, filamentous haemagglutinin, and pertactin in chromatography workflows	CY13819
	High-throughput screening for elution conditions on Capto MMC using PreDicator plates	CY13156
	Optimizing elution conditions on Capto MMC using Design of Experiments	CY14386
	Use of sodium hydroxide for cleaning and sanitization of chromatography resins and systems	CY13951
Instructions	Tricorn Empty High Performance Columns	28409488
	HiScale columns (10, 16, 26, 50) and accessories	28967470
	Packing Capto adhere and Capto MMC using verified methods	CY29524

Page intentionally left blank



Give feedback on this document

Visit 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, AxiChrom, BioProcess, Capto, ChromaFlow, HiScale, HiScreen, HiTrap, PreDictor, ReadyToProcess, Sepharose, and Tricorn are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

RoboColumn is a trademark of Repligen GmbH.

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

© 2020–2024 Cytiva

For local office contact information, visit cytiva.com/contact

11003505 AH V:2 03/2024