

CaptureSelect™ hCG Affinity Matrix

Catalog Numbers 1943410250, 1943410500, 19434101L, 19434105L

Pub. No. MAN0016318 **Rev.** A.0

Product description

The CaptureSelect™ hCG Affinity Matrix purifies human Chorionic Gonadotropin (hCG) from complex source materials (recombinant sources) in a single step.

The matrix has selectivity for the alpha chain of human Chorionic Gonadotropin. It also cross-binds with LH, FSH, and TSH that contain this alpha chain. The 14 kDa single domain [VHH] antibody fragment provides high selectivity and is immobilized on high-quality chromatography matrix.

Product advantages

The CaptureSelect™ hCG Affinity Matrix offers:

- High recovery and purity in a single step
- Selective binding of the alpha chain of hCG
- Compatibility with FPLC systems

Product specifications

Ligand	Human Chorionic Gonadotropin affinity ligand
Binding specificity	Alpha chain of human Chorionic Gonadotropin
Matrix and particle size	Aldehyde-activated agarose, 65 µm
Dynamic binding capacity	~4 g of hCG/L of matrix
Shipping solution	20% (v/v) ethanol

Conditions for use

Parameter	Conditions for use
Equilibration buffer	20 mM Tris or PBS, pH 7.0–7.5
Elution buffer	50 mM citric acid, pH 3.0
Strip buffer	Any of the following: <ul style="list-style-type: none"> • 0.1 M glycine, pH 2.0 • 0.5–1.0 M acetic acid • Citric acid
Flow rate	50–150 cm/h
Pressure limit	≤ 2 bar
Cleaning solution	Any of the following: <ul style="list-style-type: none"> • 0.5–1.0 M acetic or citric acid • 10 mM NaOH, pH 12 (Higher concentrations affect the functionality of the affinity ligand on the matrix.) • PAB (120 mM phosphoric acid, 167 mM acetic acid, and 2.2% (v/v) benzyl alcohol) (Rogers <i>et al.</i>, 2009) Freshly prepare PAB every 4–5 days and store protected from light to minimize radicals that affect the functionality of the matrix.
Storage solution	20% (v/v) ethanol
Operating temperature	2–25°C

Contents and storage

Product	Cat. No.	Amount	Storage
CaptureSelect™ hCG Affinity Matrix	1943410250	250 mL	Room temperature for < 2 weeks
	1943410500	500 mL	
	19434101L	1 L	
	19434105L	5 L	

Caution: For manufacturing, processing or repacking.

Flow characteristics

You can use agarose-based CaptureSelect™ affinity matrices at flow rates of 50–250 cm/h (Figure 1).

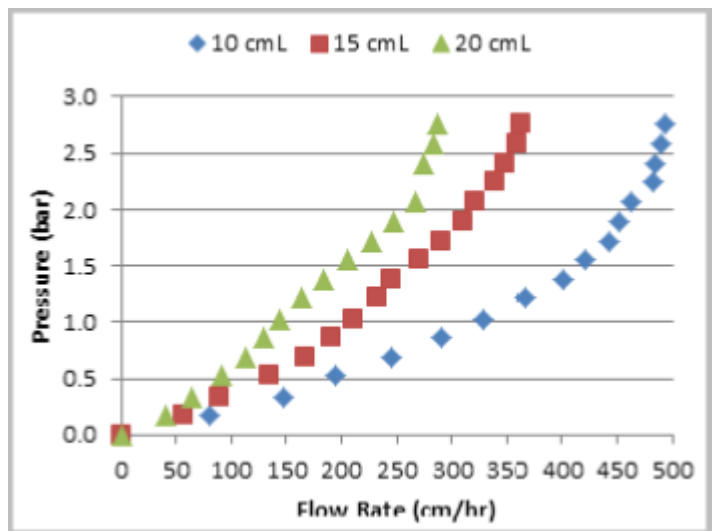


Figure 1 Typical pressure-flow curve for CaptureSelect™ resin at increasing bed heights: 10 cm (blue), 15 cm (red), and 20 cm (green). (10-cm diameter column packed at 3 bar; mobile phase=0.1 M NaCl).

The resin can be operated at flow rates up to 250 cm/h, with a pressure drop that allows use in conventional low-pressure chromatography columns and systems. However, for optimal binding capacity and elution efficiency, we recommend flow rates of 50–150 cm/h. A low flow rate results in longer contact time of the load with the affinity matrix and drives the binding capacity. In addition, the elution fraction is more concentrated at a lower flow rate.

We recommend that you optimize each of your specific processes to achieve the best conditions for process time, binding capacity, and elution efficiency.

Guidelines for use - FPLC

For optimal matrix performance, optimize the conditions in the following procedure for your application.

1. Pack the column as described in *CaptureSelect™ Affinity Matrices: Guidelines for Packing* (Pub. no. MAN0009645).
2. Attach the packed column to the FPLC system.
3. Equilibrate the matrix with 10 column volumes (CVs) of equilibration buffer.
4. Determine the volume of sample to load, based on the dynamic binding capacity, concentration of the target molecule, and the column size. Optimum loading is at physiological pH. Avoid acidic conditions, which decrease binding efficiency.
5. Load the sample on the column.
6. Wash the sample with 5–10 CVs of equilibration buffer. To optimize washing efficiency, you can add NaCl to the equilibration buffer (up to 1.0 M).
7. Elute with 3–5 CVs of elution buffer.
8. Re-equilibrate the column in equilibration buffer.

9. Strip the column with 0.1 M glycine (pH 2.0), citric acid, or acetic acid (0.5–1.0 M).
10. Re-equilibrate the column in equilibration buffer to prepare the column for another purification run.
11. If the column will not be used immediately, store the matrix according to the storage parameters provided in “Conditions for use” on page 1.

Cleaning guidelines

Resin lifetime depends on how the resin is used and cleaned. Therefore, we recommend that you specifically evaluate each purification process.

Typical cleaning procedures for CaptureSelect™ resins include combinations of acidic cleaning followed by low concentrations of NaOH, before storing in 20% (v/v) ethanol at neutral pH (Eifler *et al.*, 2014). Figure 2 shows the results of our cleaning solution testing.

To optimize column cleaning, consider these guidelines:

- Pump the cleaning solution through the column for 15–20 minutes in upflow.
- Incorporate a static hold to increase the time that the cleaning solution is in the column while minimizing the volume of cleaning solution required.
- When a combination of acidic and mildly caustic cleaning agents is used, apply the NaOH solution as a final cleaning agent to minimize the risk of irreversibly binding impurities on the column.
- In some purification processes, 20% (v/v) isopropanol (with or without acid) and 6.0 M guanidine-HCl can help remove discoloration.
- The CaptureSelect™ hCG Affinity Matrix was exposed to several cleaning agents for up to 96 hours at ambient temperature. The functionality of the resin was measured every 24 hours to test compatibility of the matrix with these cleaning agents (Figure 2).

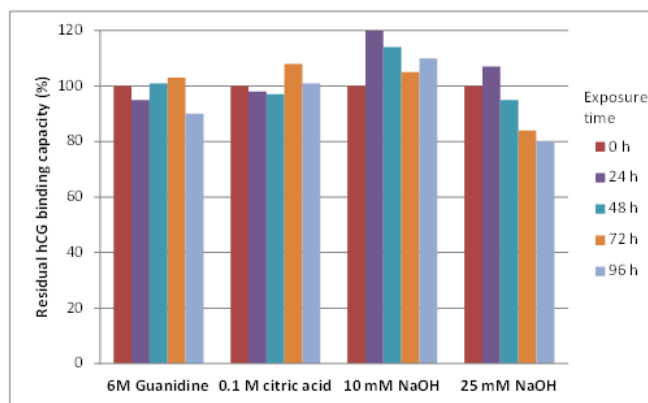


Figure 2 The CaptureSelect™ hCG Affinity Matrix was exposed to several cleaning agents for up to 96 hours at ambient temperature. The functionality of the resin was measured every 24 hours to test compatibility of the matrix with the cleaning agents. The results show that the matrix is compatible with acidic and mild caustic cleaning agents for up to 96 hours at ambient temperature. In addition, chaotropic agents like 6.0 M Guanidine- HCl are compatible with the resin.

Example application—FPLC

In this example, hCG was bound and eluted from the column. After the resin was loaded, the column was equilibrated, then eluted. Conditions were as follows:

- **Column** – CaptureSelect™ hCG Affinity Matrix
- **Equilibration buffer** – PBS
- **Load** – Partially purified hCG
- **Elution buffer** – 0.1M glycine, pH 3.0
- **Flow** – 200 cm/h

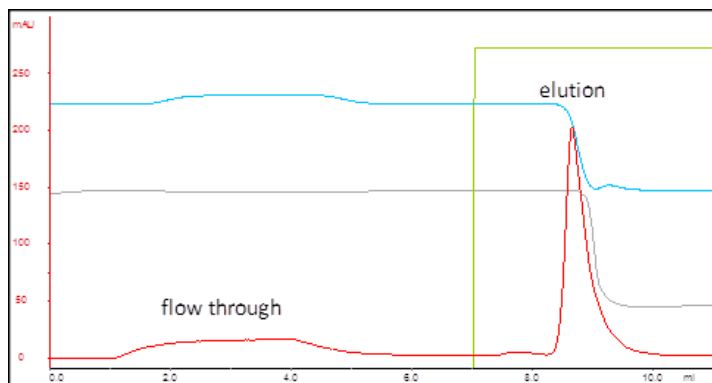


Figure 3 Chromatogram of partially purified hCG using the CaptureSelect™ hCG Affinity Matrix.

Red line: OD280 signal; Blue line: conductivity; Grey line: pH value

Supporting products

A biotinylated anti-gonadotropin conjugate is available. Applications for the conjugate include:

- ELISA
- Western blot
- Gyros™ Gyrolab™-based immunoassays
- Label-free detection platforms, such as those based on surface plasmon resonance (Biacore™ and IBIX-MX96 systems) and bio-layer interferometry (ForteBio™ Octet™ systems)

In addition, a ligand leakage ELISA is available for detecting possible leached ligand in the elution fractions of the CaptureSelect™ hCG Affinity Matrix.

Product	Size	Cat. no.
Biotin Anti-gonadotropin Conjugate	100 µg	7103412100
	500 µg	7103412500
CaptureSelect™ hCG Ligand Leakage ELISA	1 assay	810341001
	10 assays	810341010

Regulatory Support File

A Regulatory Support File (RSF) is available that contains detailed information about the resin and the manufacturing process. For more information about the RSF, contact your local sales representative.

For more information

For more information on CaptureSelect™ products and ligand leakage ELISA products, go to www.thermofisher.com/captureselect.

Customer and technical support

Visit thermofisher.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
 - Product support, including:
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
 - Order and web support
 - Product documentation, including:
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)
- Note:** For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

References

- Rogers, M. *et al.* 2009. Development of a rapid sanitization solution for silica-based protein A affinity adsorbents. *Journal of Chromatography A*. 1216:4589–4596.
- Eifler, N. *et al.* 2014. Development of a novel affinity chromatography resin for platform purification of lambda fabs. *Biotechnology Progress* DOI:10.1002/btpr.1958.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0016318

Revision	Date	Description
A.0	30 November 2016	New document.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Biacore is a trademark of GE Healthcare Bio-Sciences AB. ForteBio and Octet are trademarks of Pall Fortebio Corp. Gyros and Gyrolab are trademarks of Gyros Patent AB.