



Product Information

MagSelect™ SA

Product Number **M 0444**

Storage Temperature 2–8 °C

TECHNICAL BULLETIN

Synonym: Streptavidin Magnetic Agarose Beads

Product Description

Streptavidin, a 66 kDa homotetrameric protein isolated from *Streptomyces avidinii*, is similar to avidin with respect to its high affinity for biotin ($K_a \sim 10^{15} \text{ M}^{-1}$).¹⁻⁴

Streptavidin is slightly anionic (pI ~5.6) and non-glycosylated - both properties contributing to the relatively low non-specific binding compared to egg white avidin (a glycoprotein with pI ~10.5).

The binding of biotinylated biomolecules to streptavidin is virtually irreversible. Very harsh conditions including low pH and denaturants are needed to completely remove the biotinylated molecules. This attribute makes MagSelect™ SA beads very useful in a variety of affinity capture applications. Examples include capture of biotinylated macromolecules and complexes such as proteins, antibodies, lectins, nucleic acids, receptors, and ligands.⁵⁻¹¹ MagSelect SA beads can also be used for direct purification of desthiobiotinylated proteins and peptides. This allows the molecule of interest to be eluted from the beads under mild conditions, using competitive displacement with biotin.¹²

MagSelect SA is composed of streptavidin attached to paramagnetic iron impregnated 6% agarose beads, with an average diameter of 50 μm . The resulting linkage has an effective spacer length of 14 atoms. The magnetic properties allow for very rapid separation of the beads from a suspension, significantly accelerating manipulations, such as repetitive washings, or processing of multiple samples performed in multiwell plates. That leads to faster experimentation, better reproducibility, and more accurate quantitation of the proteins of interest.

Binding capacity:

- Minimum 175 nmoles of biotin per ml of chromatographic medium
- minimum 2 mg of biotinylated BSA per ml of chromatographic medium

Specificity:

MagSelect SA has been tested for specificity of binding biotinylated proteins. Non-specific binding of 5% or less is observed when the reagent is used in a RIPA Buffer system (Product Code R 0278).

Reagents

MagSelect SA is supplied as a 50% suspension in 0.01 M sodium phosphate, pH 7.2, containing 50% glycerol, 0.15 M NaCl, and 15 ppm Kathon® CG/ICP II (as a preservative).

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

For ease of handling, the glycerol concentration may be reduced by dilution with a neutral buffer such as Phosphate Buffered Saline, pH 7.4, (PBS, Product Code P 3813) prior to use.

Equilibration for Batch Format - For the procedure described in this bulletin, 50 μl (100 μl of suspension) of MagSelect SA is utilized. Transfer the required amount of resin slurry to an appropriately sized microcentrifuge tube. When pipetting, use a wide orifice tip or cut ~1 mm off the end of a regular pipette tip to allow for unrestricted flow of the bead suspension. To equilibrate the medium, add 5 volumes of PBS, mix, and centrifuge for 30 seconds at 2,000 $\times g$ (e.g., 5,000 rpm in a Eppendorf® 5415 C microcentrifuge). Carefully remove the supernatant using a micropipette. Repeat the equilibration step two more times.

Storage/Stability

MagSelect SA ships on wet ice and storage at 2–8 °C is recommended. The product, as supplied in 50% glycerol with preservative, may be stored at –20 °C. However, freezing the magnetic beads in the absence of 50% glycerol will irreversibly damage the bead structure. For long term storage of the medium following equilibration of the beads and removal of glycerol, the medium should be stored at 2–8 °C with a preservative (e.g. 15 ppm of Kathon CG/ICP II, Product Code 48178U) and **not frozen**.

Procedures

A. Procedure for Immunoaffinity Batch Purification Using a Biotinylated Antibody as a Ligand

1. Transfer the required amount of equilibrated medium (see Preparation Instructions) to an appropriately sized microcentrifuge tube. For this procedure, 50 µl of MagSelect SA is utilized.
2. Add the biotinylated antibody dissolved in an appropriate neutral buffer such as PBS. To prevent non-specific binding of non-biotinylated molecules, detergents or blocking agents may also be included in the solution. Fifty µl of MagSelect SA will bind ~50 µg of a biotinylated antibody. For optimal binding, incubate with mixing for a minimum of 15 minutes at room temperature.
3. Wash away any unbound antibody with PBS. Specifically, add 500 µl of PBS, then mix and place on an appropriate magnetic device for 20 seconds. Remove the supernatant. Repeat this step five more times.
4. Apply the sample (i.e. mixture containing antigen) to the microcentrifuge tube. For optimal binding, incubate with mixing for a minimum of 15 minutes at room temperature.
5. Wash away any unbound biomolecules with PBS. Specifically, add 500 µl of PBS then mix and place on an appropriate magnetic device for 20 seconds. Remove the supernatant. Repeat this step five more times.
6. Elute the sample (i.e. antigen) with 0.1 M acetic acid or 0.1 M glycine HCl (pH 2.5) or other elution buffer to dissociate the antibody-antigen binding interaction. Immediately neutralize eluted samples, which are now ready for further analysis.

B. Purification of Desthiobiotinylated Biomolecules in a Centrifuge Tube Batch Format

1. Transfer the required amount of equilibrated medium (see Preparation Instructions) to an appropriately sized microcentrifuge tube. For this procedure, 50 µl of MagSelect SA is used.
2. Add up to 12 nmoles of desthiobiotinylated sample, dissolved in an appropriate neutral buffer such as PBS. To facilitate optimal binding, incubate with mixing for a period of at least 15 minutes at room temperature.
3. Wash away any unbound biomolecules with PBS. Specifically, add 500 µl of PBS then mix and place on an appropriate magnetic device for 20 seconds. Remove the supernatant. Repeat this step five more times.
4. To elute the desthiobiotinylated biomolecules, add 200 µl of 2 mM biotin in PBS. Allow the solution to incubate at room temperature with mixing for at least 15 minutes.
5. Place on an appropriate magnetic device, then remove and save the supernatant containing the desthiobiotinylated biomolecules of interest for further analysis.

Related Products	Product Code
Anti-FLAG® BioM2 Antibody	F 9291
Protein A – Biotin	P 2165
Protein G – Biotin	P 8045
Desthiobiotin PEO Iodoacetamide	D 2192
Magnetic separators for microcentrifuge tubes	M 1167
Magnetic separators for tissue culture flasks	M 1292
Magnetic separators for centrifuge tubes	M 1542

References

1. Chaiet, L., and Wolf, F.J., The properties of streptavidin, a biotin-binding protein produced by *Streptomyces*. *Arch. Biochem. Biophys.*, **106**, 1-5 (1964).
2. Bayer, E.A., *et al.*, Isolation and properties of streptavidin. *Methods Enzymol.*, **184**, 80-89 (1990).
3. Bayer, E.A., and Wilchek, M., Avidin-Biotin Technology. *Methods in Molecular Biology*, Vol. 10, *Immunochemical Protocols*, p. 137-162 (1992).
4. Green, M.N., Avidin and Streptavidin. *Methods Enzymol.*, **184**, 51-67 (1990).
5. Von Boxberg, Y., *et al.*, Use of the biotin-avidin system for labeling, isolation and characterization of neural cell-surface proteins. *Eur. J. Biochem.*, **190**, 249-256 (1990).
6. Updyke, T.V., and Nicolson, G.L., Immunoaffinity isolation of membrane antigens with biotinylated monoclonal antibodies and immobilized streptavidin matrices. *J. Immunol. Methods*, **73**, 83-95 (1984).
7. Updyke, T.V., and Nicolson, G.L., Immunoaffinity isolation of membrane antigens with biotinylated monoclonal antibodies and streptavidin-agarose. *Methods Enzymol.*, **121**, 717-25 (1986).
8. Gretch, D.R. *et al.*, The use of biotinylated monoclonal antibodies and streptavidin affinity chromatography to isolate herpesvirus hydrophobic proteins or glycoproteins. *Anal. Biochem.*, **163**, 270-277 (1987).
9. Wilchek, M., and Bayer, E.A., Applications of Avidin-Biotin Technology: Literature Survey. *Methods Enzymol.*, **184**, 14-45 (1990).
10. Buckie, J.W., and Cook, G.M.W., Specific isolation of surface glycoproteins from intact cells by biotinylated concanavalin A and immobilized streptavidin. *Anal. Biochem.*, **156**, 463-472 (1986).
11. Diamandis, E.P., and Christopoulos, T.K., The biotin-(strept)avidin system: Principles and applications in biotechnology. *Clin. Chem.*, **37**, 625-636 (1991).
12. Hirsch, J.D. *et al.*, Easily Reversible Desthiobiotin Binding to Streptavidin, Avidin, and Other Biotin-Binding Proteins: Uses for Protein Labeling, Detection, and Isolation. *Anal. Biochem.*, **308**, 343-357 (2002).

Kathon is a registered trademark of Rohm and Haas Company.

Eppendorf is a registered trademark of Eppendorf-Netheler-Hinz GmbH.

CDM,KLF,MAM 12/04-2

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.