



274PR-01

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Immobilized Iminobiotin Resin

(Cat. #786-599)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

G-Biosciences Immobilized Iminobiotin Resin consists of iminobiotin, a cyclic guanido analog of biotin, covalently coupled to 6% crosslinked agarose. The resin allows for the purification of avidin, streptavidin and Neutraavidin and their subsequent gentle elution using non-denaturing elution buffers.

The normal biotin-avidin complex requires strong denaturing buffers, i.e. 8M guanidine•HCl, to denature the avidin and release the biotin, which obviously destroys the native and functional aspects of the avidin. The iminobiotin-avidin complex will form at >pH9.5 and can be dissociated at pH4.0 with gentle elution buffers, including 50mM ammonium acetate, pH4.0 with 0.5M NaCl.

ITEM(S) SUPPLIED (Cat. #786-599)

Description	Size*
Iminobiotin, Immobilized	5ml resin

** Immobilized Iminobiotin resin is supplied as a 50% slurry with 0.02% sodium azide as a preservative.*

STORAGE CONDITIONS

It is shipped at ambient temperature. Upon arrival, store refrigerated at 4°C, **DO NOT FREEZE**. This product is stable for 1 year at 4°C.

SPECIFICATIONS

Binding Capacity: >2mg avidin/ml resin
Bead Structure: 6% cross-linked agarose

ADDITIONAL ITEMS

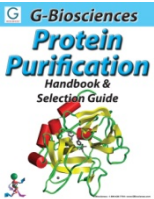
- Avidin, streptavidin or Neutraavidin labeled antibody or protein
- Binding/Wash Buffer: 50mM ammonium carbonate, pH11.0 with 0.5M NaCl
- Elution Buffer: 50mM ammonium acetate, pH4.0 with 0.5M NaCl
- Columns (optional): G-Biosciences offers columns for a large range of resin volumes (Cat. # 786-718 to 786-724)

PROCEDURE

1. Allow the resin and reagents to equilibrate to room temperature.
2. Add an appropriate volume of homogenous iminobiotin resin to a suitable column.
3. Wash the resin with 4-5 resin volumes of Binding/Wash Buffer.
4. Apply the sample to the column and add the bottom then top cap. Incubate at room temperature for 30 minutes.
5. Elute the protein with 1 resin volume of Elution Buffer. Repeat the elution 3-6 times and monitor elutions by absorbance at 280nm.
6. The samples can be used directly for SDS PAGE, or alternatively, can be dialyzed for specific downstream applications.
7. Wash the resin with 4-5 resin volumes of Binding/Wash Buffer. For long term storage supplement the Binding/Wash Buffer with 0.02% sodium azide. Store upright at 4°C. The resin can be used at least 10 times.

RELATED PRODUCTS

Download our Protein Purification Handbook.



<http://info2.gbiosciences.com/complete-protein-purification-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 3/18/2014 TNN



www.GBiosciences.com