



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

A Geno Technology, Inc. (USA) brand name

N-acetyl-D-Glucosamine Agarose

(Cat. #786-270)



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

INTRODUCTION

N-Acetyl-D-Glucosamine Agarose is used in carbohydrate affinity chromatography for purification of wheat germ (*Triticum vulgaris*) lectin, other N-acetyl-glucosamine specific lectins such as lectin from *Dolichos biflorus* and carbohydrate binding proteins/enzymes.

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-270	N-Acetyl-D-Glucosamine Agarose	5 ml resin

Supplied as a 50% suspension in 50% glycerol containing 0.01% sodium azide

STORAGE CONDITION

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

SPECIFICATIONS

- Ligand: N-acetyl-D-glucosamine
- Matrix bead structure: 6 % cross-linked agarose
- Matrix bead size: 45-160 µm
- Matrix activation: Epoxy
- Matrix binding to ligand: via –OH groups
- Spacer arm: 12 atoms
- Binding capacity: ≥ 25 mg wheat germ lectin/ml of resin

ADDITIONAL ITEMS REQUIRED

- Columns
- Binding buffer: 0.1M sodium phosphate, 0.15M sodium chloride, pH7.2 or other suitable buffer.
- Elution buffer: 0.1-0.2 M N-acetyl-D-glucosamine in binding buffer

PROTOCOL

1. Aliquot the desired volume of N-Acetyl-D-Glucosamine Agarose in the column.
2. Wash the column with 10 column volumes (CV) of working or binding buffer.
3. Equilibrate the column with 10 CV of working or binding buffer.
4. Load the protein solution and wash the unbound protein.

NOTE: Wash to remove all unbound protein. The unbound protein in the eluent can be monitored by checking the absorbance at 280 nm.

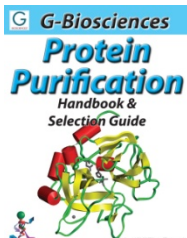
5. Elute the specifically bound lectin or carbohydrate binding protein by competitive elution with 0.1 to 0.3 M elution sugar added to the working buffer or perform elution at different pH and/or ionic strength.

RESIN CLEANING AND REGENERATION

1. Wash the resin with 10 CV of binding buffer or working buffer containing 0.1 to 0.3M elution sugar.
2. Wash the resin with 10 CV of 2 M NaCl solution.
3. Wash the resin with 10 CV of working buffer and store the resin in working buffer or storage solution at 4°C.

RELATED PRODUCTS

Download our Protein Purification Handbook.



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

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

Last saved: 8/2/2016 CMH



www.GBiosciences.com