



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

A Geno Technology, Inc. (USA) brand name

Immobilized Jacalin

For the purification of Immunoglobulin A (IgA)

(Cat. #786-167)



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

INTRODUCTION

Jacalin, or *Artocarpus integrifolia* lectin, is a tetrameric two-chain lectin with a molecular weight of 66kDa. Jacalin is a α -D-galactose binding lectin purified from jack-fruit (*Artocarpus integrifolia*) seeds. Applications include isolating IgA from human serum and colostrums, isolating human plasma glycoproteins and histochemistry. Jacalin also binds IgD.

ITEMS SUPPLIED Cat. # 786-167

Description	Size
Immobilized Jacalin	2ml Resin

*SUPPLIED AS 50% SLURRY CONTAINING 0.02% SODIUM AZIDE AS A PRESERVATIVE.

ADDITIONAL MATERIALS

1X Phosphate Buffered Saline (PBS)

Elution Buffer: 0.1M Melibiose (6-O- α -D-Galactopyranosyl-D-glucose) or 0.1M α -D-galactose in PBS

- *Melibiose is a disaccharide formed by an alpha linkage between galactose and glucose (D-Gal- α (1 \rightarrow 6)-D-Glc)*

Disposable columns

STORAGE CONDITIONS

The kit is shipped at ambient temperature. Upon arrival, store at 4°C (DO NOT FREEZE).

SPECIFICATIONS

Binding Capacity: 1-3mg human IgA/ml resin

Loading: \approx 4.5mg jacalin/ml of resin

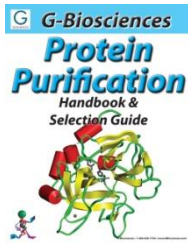
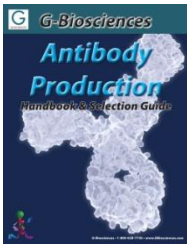
Support: 6% cross-linked agarose

PROTOCOL FOR HUMAN IGA PURIFICATION

1. Equilibrate the Immobilized Jacalin to room temperature and pack a suitable amount into a column.
2. Wash the column with 5 column volumes of 1X PBS
3. Dilute the human serum 1:1 with PBS and add to the column.
4. Wash the column with a further 5 column volumes of PBS. To ensure adequate washing, monitor the flow through using an absorbance of 280nm and continue washing until no further change is seen.
5. To recover the bound IgA, add repeating single column volumes of Elution Buffer and monitor elution with absorbance readings at 280nm.
6. To remove the melibiose or galactose sugar, perform a buffer exchange into PBS using desalting columns. We recommend our SpinOUT™ GT600 columns. Dialysis can also be used and we recommend our Tube-O-DIALYZER™ dialysis devices.
7. Regenerate the columns by washing with >20 column volumes of PBS and then store the columns in deionized water supplemented with 0.02% sodium azide.

RELATED PRODUCTS

Download our Antibody Production and Protein Purification Handbook.



<http://info2.gbiosciences.com/complete-antibody-production-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: 2/17/2014 TN



www.GBiosciences.com