



548PR-02

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

A Geno Technology, Inc. (USA) brand name

Immobilized Protein A

0.2ml Spin Column Kit

(Cat. # 786-828)



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

INTRODUCTION 3

ITEM(S) SUPPLIED 3

STORAGE CONDITIONS 3

SPECIFICATIONS 3

ADDITIONAL ITEMS REQUIRED 5

PREPARATION BEFORE USE 5

PROTOCOL 5

TROUBLESHOOTING 6

RELATED PRODUCTS 6

INTRODUCTION

Immobilized Protein A consists of recombinant protein A ligand covalently immobilized onto 4% highly cross-linked agarose. The dynamic binding capacity will vary depending on several factors such as target antibody, flow rate etc.

The recombinant protein A, produced in *E. coli*, has been designed to retain solely its binding properties for immunoglobulins. Immobilized protein A is widely used for the isolation and purification of a wide variety of immunoglobulins from a variety of species (see table below).

The Immobilized Protein A 0.2ml spin column kit is ideal for the small scale affinity purification of antibodies from a variety of samples. Each 0.2ml column enables purification of up to 8mg IgG from 25-500µl of serum or other sample.

ITEM(S) SUPPLIED

Part. #	Description	Size
442P-B	IgG Elution Buffer	2 x 30ml
001J	JAW Phosphate Buffered Saline Pack	For 1L
228C-B	Collection Tube, 2ml	100
440P-D	Immobilized Protein A Resin*	10 x 0.2ml columns
411S-B	Stoppers, Rubber (Small)	10/bag

**Immobilized Protein A is supplied as a 50% slurry in 20% ethanol/PBS solution*

STORAGE CONDITIONS

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

SPECIFICATIONS

- High binding capacity: >40mg human IgG/ml resin
- Ligand: Recombinant Staphylococcal Protein A lacking the albumin-binding domain produced in *E. coli*
- Bead size: 45-165µm
- Bead Structure: 4% highly cross-linked agarose

Species	Antibody Class	Protein A	Protein G	Protein A/G
Mouse	Total IgG	++++	++++	++++
	IgG ₁	+	+++	+++
	IgG _{2a}	++++	++++	++++
	IgG _{2b}	++++	++++	++++
	IgG ₃	+++	+++	++++
Human	Total IgG	++++	++++	++++
	IgG ₁	++++	++++	++++
	IgG ₂	++++	++++	++++
	IgG ₃	+	++++	++++
	IgG ₄	++++	++++	++++
Rat	Total IgG	+	++	+++
	IgG ₁	-	+	+++
	IgG _{2a}	-	++++	++++
	IgG _{2b}	-	++	+
	IgG _{2c}	++	+++	++++
Hamster	Total IgG	++	++	++
Guinea Pig	Total IgG	++++	++	++++
Rabbit	Total IgG	++++	+++	++++
Horse	Total IgG	++	++++	++++
Cow	Total IgG	++	++++	++++
Pig	Total IgG	+++	++	++++
Sheep	Total IgG	+	++	++++
Goat	Total IgG	+	++	++++
Chicken	Total IgG	-	-	-

Table 1: Relative affinity of Protein A, Protein G & Protein A/G for Immunoglobulins

ADDITIONAL ITEMS REQUIRED

- Storage Buffer: 10mM NaH₂PO₄, 150mM NaCl, 2.7mM KCl, pH 7.4, 20% ethanol
- Neutralization Buffer: 1M Tris-HCl, pH8.5
- Microcentrifuge

PREPARATION BEFORE USE

Binding Buffer: Add the entire contents of the JAW Phosphate Buffered Saline pack to 1 liter of deionized water. Stir until completely dissolved.

PROTOCOL

NOTE: *The Immobilized Protein A columns can be used up to 10 times without significant loss of binding capacity.*

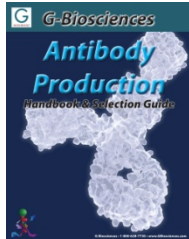
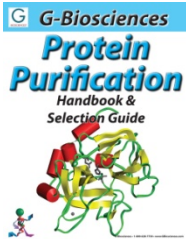
1. Allow the buffers and columns to equilibrate to room temperature.
2. Place the columns in the collection tubes and briefly centrifuge at 5,000xg for 0.5-1 minute to pack the resin.
3. Snap off the bottom of the column and remove the top cap. Return the columns to the collection tubes and centrifuge at 5,000xg for 1 minute to remove the storage buffer. Discard the flow through.
4. Equilibrate the resin by adding 400µl Binding Buffer to the column in a collection tube. Centrifuge the column for 1 minute at 5,000xg and discard the flow through. Repeat this step once.
5. Cap the bottom of the column and gently apply 25-500µl sample to the column by adding to the top of the resin. Do not disturb the gel bed. Cap the column.
6. Incubate the column for 10 minutes at room temperature with end-over-end mixing, when volumes allow mixing to occur.
7. Remove the top cap and then the bottom cap. Place the spin column in a new collection tubes and centrifuge at 5,000xg for 1 minute. The flow through contains the non-bound sample and can be analyzed to determine binding efficiency.
8. Transfer the column to a new collection tube and apply 400µl Binding Buffer to the resin. Mix briefly to suspend the resin and then centrifuge for 1 minute. Repeat this wash step two additional times.
9. Add 40µl Neutralization Buffer to three clean collection tubes. Place the spin column in one of the tubes.
10. Add 400µl IgG Elution Buffer to the spin column, mix gently and then centrifuge for 1 minute at 5,000xg. Transfer the spin column to another collection tube containing Neutralization Buffer. Repeat this step two more times.
11. Identify the immunoglobulin-containing fractions by measuring absorbances at 280nm or using a suitable protein assay. (NI-Protein Assay Cat. # 786-005)
12. If using for further downstream applications, exchange the buffer with our SpinOUT™ desalting columns or Tube-O-DIALYZER™ dialysis devices.
13. To regenerate the column, wash the column four times with Elution Buffer, followed by four washes with Storage Buffer. Do not allow the resin to dry out.

TROUBLESHOOTING

Issue	Possible Reason	Solution
No protein detected in eluted fractions	Initial sample is devoid of any antibody species or isotype that bind protein A.	Check Table 1 for relative affinity. Ensure by ELISA or isotyping that the sample contains the correct IgG type.
Large amount of antibody purified, but specific antibody of interest not detected	Antibody of interest has low affinity for the resin or is at low concentration	Use serum free media for cell supernatant samples
		Use affinity purification using a specific antigen coupled to a support (see Protein Purification Handbook for selection of supports)
Antibody of interest, but is denatured and inactive	Antibody is sensitive to low pH	Elute in a high salt buffer (5M LiCl, 10mM phosphate, pH7.2)
	Downstream application is sensitive to neutralized Elution Buffer	Desalt or dialyze sample

RELATED PRODUCTS

Download our Protein Purification and Antibody Production Handbooks.



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

<http://info.gbiosciences.com/complete-Antibody-Production-handbook/>

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

Last saved: 6/16/2017 CMH

This page is intentionally left blank



www.GBiosciences.com