



430PR-01

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# Immobilized *p*-Aminophenyl Phosphoryl Choline

For Purifying C-Reactive Protein

(Cat. # 786-821)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

G-Biosciences' Immobilized *p*-Aminophenyl Phosphoryl Choline consists of a phosphoryl choline covalently linked to beaded agarose and is designed for the purification of C-reactive protein from plasma, ascites and other biological fluids.

CRP, C-reactive protein, is a pentameric protein found in the blood, the levels of which rise in response to inflammation, making CRP an acute-phase protein. Its physiological role is to bind to phosphocholine expressed on the surface of dead or dying cells (and some types of bacteria) in order to activate the complement system.

## ITEM(S) SUPPLIED (Cat. # 786-821)

Description	Size
Immobilized <i>p</i> -Aminophenyl Phosphoryl Choline	1ml resin

*Supplied as a 50% slurry in 0.02% sodium azide as a preservative*

## STORAGE CONDITIONS

Shipped as ambient temperature. Upon arrival store at 4°C. Do NOT freeze.

## ADDITIONAL MATERIALS REQUIRED

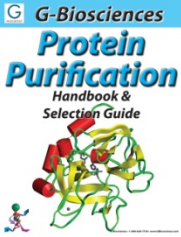
- Binding Buffer: 100mM Tris or Borate Buffer supplemented with 100-200mM NaCl, 1-2mM CaCl<sub>2</sub> at pH 8-8.5
- Elution Buffer: 100mM Tris or Borate Buffer supplemented with 100-200mM NaCl, 2mM EDTA at pH 8-8.5
- Storage Buffer: 10mM Sodium phosphate, 90mM NaCl, pH 8-8.5

## PROTOCOL

1. Prepare a suitable column by packing an appropriate volume of Immobilized *p*-Aminophenyl Phosphoryl Choline into a disposable column.
2. Apply 2 column volumes of Binding Buffer to equilibrate the resin.
3. Add approximately 0.5CV of purified sample to the resin.
4. Incubate for 1 hour at room temperature with constant mixing.
5. Wash the column with at least 5CV of Binding Buffer.
6. Elute bound C-reactive protein with Elution Buffer. Collect 0.5ml fractions and monitor elution with a suitable protein assay.
7. After elution, perform a buffer exchange by dialysis with the Storage Buffer.

## 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](http://www.GBiosciences.com) or contact us.

Last saved: 8/6/2012 CMH



[www.GBiosciences.com](http://www.GBiosciences.com)