



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 O- Phosphorylethanolamine

For the Purification of C-Reactive Protein

(Cat. # 786-1310)



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

## INTRODUCTION

G-Biosciences Immobilized O-Phosphorylethanolamine is used in purification of calcium binding C-reactive proteins such as canine CRP<sup>1</sup> and human CRP<sup>2</sup>.

Immobilized O-Phosphorylethanolamine can also be used as substitute for p-Amino Phosphoryl Choline Agarose which is used in purification of C-reactive protein (CRP).

Immobilized O-Phosphorylethanolamine is more economical when compared to p-Amino Phosphoryl Choline Agarose

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-1310	Immobilized O-Phosphorylethanolamine	5ml

## STORAGE CONDITIONS

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

## SPECIFICATIONS

- Ligand: O-Phosphorylethanolamine
- Matrix: 6% cross-linked Agarose
- Ligand density: 3-5mg/ml resin
- Storage buffer: 20 % isopropanol
- Storage temperature: 4°C

## ADDITIONAL ITEMS REQUIRED

- Sample containing CRP
- Binding Buffer: 50 mM Tris, 150 mM NaCl, 2 mM CaCl<sub>2</sub>, pH 8.0
- Elution Buffer: 10 mM Tris, 150 mM NaCl, 2mM EDTA, pH 8.0
- CRP storage buffer: 10mM Sodium Phosphate, 90 mM NaCl

## PROTOCOL

1. Pour the required amount of resin into a disposable column (G-Trap™ 5ml FPLC Column, Cat. # 786-1291)
2. Equilibrate the resin with 2-3 column volumes (CV) of Binding Buffer.
3. Plug the bottom of the column with stop plug and then add the sample.
4. Incubate the sample with resin for 15-30 minutes at room temperature.
5. Remove the stop plug and collect the flow through
6. Wash the column with 5 CV of Binding Buffer
7. Elute the bound protein with Elution Buffer and collect fractions as 0.5 ml aliquot per vial. Check the OD<sub>280nm</sub> of the fractions to identify the fraction that has CRP.

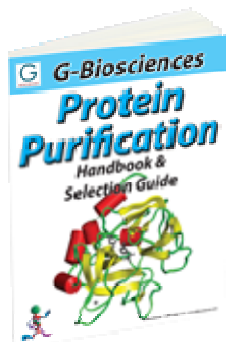
8. Collect the eluted fractions containing the CRP and dialyze the collected fractions against CRP storage buffer at 4°C overnight.

## REFERENCES

1. Onishi, T. et al. (1993). J. Vet. Med Sci. 56(2):417-419.
2. Pontet, M. et al. (1978). FEBS Lett. 88.172-175.
3. Hokama, Y. and Nakamura, R. M. (1987) J. Clin. Lab. Anal. 1: 15-27.

## 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.



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