

# Immobilized *p*-Aminophenyl Phosphoryl Choline Gel

20307

1156.4

Number	Description
20307	<b>Immobilized <i>p</i>-Aminophenyl Phosphoryl Choline Gel, 5mL</b> Support: Crosslinked 6% beaded agarose Binding Capacity: $\geq$ 3mg of human C-reactive protein per milliliter of gel  <b>Storage:</b> Upon receipt store at 4°C. Product is shipped at ambient temperature.

## Introduction

Within the ascites fluid of humans and rabbits exist several different types of immunological proteins that provide aid during an inflammatory response. One of these proteins is the C-reactive protein that has been linked to several biological functions including activation of the classical complement pathway, enhancement of phagocytosis and interaction with certain subpopulations of T-lymphocytes. In the late 1970s, Volanakis and colleagues discovered a method of isolating and studying C-reactive protein using the protein's affinity for phosphoryl choline.<sup>1</sup> Thermo Scientific Immobilized *p*-Aminophenyl Phosphoryl Choline Gel enables purification of C-reactive protein in a convenient column format that is quick and easy to use.

## Additional Materials Required

- Sample purified from ascites fluid by centrifugation and filtration
- Binding Buffer: 0.1M Tris or Borate Buffer, 0.1-0.2M NaCl, 1-2mM CaCl<sub>2</sub>; pH 8-8.5
- Elution Buffer: 0.1M Tris or Borate Buffer, 0.1-0.2M NaCl, 2mM EDTA; pH 8-8.5
- Storage Buffer: 0.01M sodium phosphate, 0.09M NaCl

## Procedure for C-reactive Protein Purification

1. Prepare a 5mL column by packing gel into a disposable column.
2. Equilibrate column by applying two column volumes of Binding Buffer.
3. Add 2-3mL of purified sample to column.
4. Incubate for 1 hour at room temperature.
5. Wash with 5 column volumes of Binding Buffer.
6. Elute bound protein with Elution Buffer. Collect fractions in 0.5mL aliquots.
7. After elution, micro-concentrate the protein and dialyze into the Storage Buffer.

## Cited Reference

1. Volanakis, J.E. *et al.* (1978). C-reactive protein: Purification by affinity chromatography and physicochemical characterization. *J Immunol Meth* 23:285-95.

---

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at [www.thermoscientific.com/pierce](http://www.thermoscientific.com/pierce). For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2012 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.