

Product Information



FABP3 Blocking Peptide

Catalog No. 10273

Fatty acid binding protein 3 (FABP3) is among the nine known cytosolic fatty acid binding proteins ranging from 14-15 kDa in molecular mass containing 127-132 amino acids and was named after the first tissue of isolation.¹ FABP3 is the most widely distributed FABP. It is found in heart, skeletal and smooth muscle, mammary epithelial cells, aorta, distal tubules of the kidney, lung, brain, placenta, and ovary. Studies have shown the role of this protein as a potential biomarker for myocardial injury especially for early detection of acute myocardial infarction (AMI).¹

Laboratory Procedures

This vial contains 200 µg of peptide in 200 µl of TBS, pH 7.4, containing 0.1% BSA and 0.02% sodium azide. The FABP3 blocking peptide (human FABP3 amino acids 44-55) can be used in conjunction with Cayman's FABP3 Monoclonal Antibody (Catalog No. 10233) to block protein-antibody complex formation during the immunochemical analysis of FABP3.

Store this peptide solution at -20°C. It will be stable for at least one year. To block antibody/protein complex formation, the following procedure is recommended:

1. Mix the monoclonal antibody and blocking peptide together in a 1:5 (v/v) ratio in a microfuge tube. For example, mix 10µl of antibody and 50µl of peptide.*
2. Incubate for 1 hour at room temperature with occasional mixing.
3. Dilute the reagents to the final working antibody concentration and apply to the slide or membrane as usual.

*This is a recommended mixture. The minimum amount of peptide needed for complete blocking has not been precisely determined and may vary depending on the sample being analyzed. The amount of peptide required may need to be increased if sufficient blocking does not occur.

Reference

1. Zimmerman, A.W. and Veerkamp, J.H. New insights into the structure and function of fatty acid-binding proteins. *Cell. Mol. Life Sci.* **59**, 1096-1116 (2002).

Related Products

FABP3 Monoclonal Antibody (Clone CC68) - Cat. No. 10233 • FABP3 (human recombinant) - Cat. No. 10007432 • 1,8-ANS - Cat. No. 10010219

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 07/06/2009

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com