

Product Information



FABP4 Blocking Peptide

Item No. 10006248

FABP4 is primarily expressed in adipocytes but is also expressed in activated macrophages, indicating that the protein plays a critical role in foam cell formation and thus atherosclerosis.^{1,2} FABP4 was first purified from murine 3T3-L1 adipocytes and then from human adipocytes and was named as ALBP.^{3,4} The expression of FABP4 is developmentally regulated by fatty acids, PPAR γ agonists, and insulin.^{5,6} In mice, deficiency of FABP4 prevents the development of hyperinsulinemia and insulin resistance in genetic and diet-induced obesity.^{7,8} Cayman's FABP4 polyclonal antibody can be used in western blot and immunocytochemistry with samples of human, rat, and murine origin. The blocking peptide can be used together with the antibody for a negative control in these procedures.

Laboratory Procedures

This vial contains 200 μ g peptide in 200 μ l TBBS, pH 7.4, containing 0.1% BSA and 0.02% sodium azide. The FABP4 blocking peptide (amino acids 103-118) can be used in conjunction with Cayman's FABP4 Polyclonal Antibody (Item No. 10004944) to block protein-antibody complex formation during immunochemical analysis of FABP4.

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

1. Mix the FABP4 Polyclonal Antibody (Item No. 10004944) and blocking peptide together in a 1:1 (v/v) ratio in a microfuge tube. For example, mix 20 μ l of antibody and 20 μ l of peptide.*
2. Incubate for one hour at room temperature with occasional mixing prior to further dilution and application of the mixture to the immunoblot.
3. Dilute the mixture 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.

References

1. Perrella, M.A., Pellacani, A., Layne, M.D., *et al.* Absence of adipocyte fatty acid binding protein prevents the development of accelerated atherosclerosis in hypercholesterolemic mice. *FASEB J.* **15**, 1774-1776 (2001).
2. Boord, J.B., Maeda, K., Makowski, L., *et al.* Adipocyte fatty acid-binding protein, aP2, alters late atherosclerotic lesion formation in severe hypercholesterolemia. *Arterioscler. Thromb. Vasc. Biol.* **22**, 1686-1691 (2002).
3. Matarese, V. and Bernlohr, D.A. Purification of murine adipocyte lipid-binding protein. Characterization as a fatty acid- and retinoic acid-binding protein. *J. Biol. Chem.* **263**(28), 14544-14551 (1988).
4. Baxa, C.A., Sha, R.S., Buelt, M.K., *et al.* Human adipocyte lipid-binding protein: Purification of the protein and cloning of its complementary DNA. *Biochemistry* **28**, 8683-8690 (1989).
5. Distel, R.J., Robinson, G.S., and Spiegelman, B.M. Fatty acid regulation of gene expression. Transcriptional and post-transcriptional mechanisms. *J. Biol. Chem.* **267**(9), 5937-5941 (1992).
6. Melki, S.A. and Abumrad, N.A. Expression of the adipocyte fatty acid-binding protein in streptozotocin-diabetes: Effects of insulin deficiency and supplementation. *J. Lipid Res.* **34**, 1527-1534 (1993).
7. Hotamisligil, G.S., Johnson, R.S., Distel, R.J., *et al.* Uncoupling of obesity from insulin resistance through a targeted mutation in aP2, the adipocyte fatty acid binding protein. *Science* **274**, 1377-1379 (1996).
8. Uysal, K.T., Scheja, L., Wiesbrock, S.M., *et al.* Improved glucose and lipid metabolism in genetically obese mice lacking aP2. *Endocrinology* **141**(9), 3388-3396 (2000).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10006248

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 via email to 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, 07/19/2012

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