

Product Information



FABP4 Polyclonal Antibody

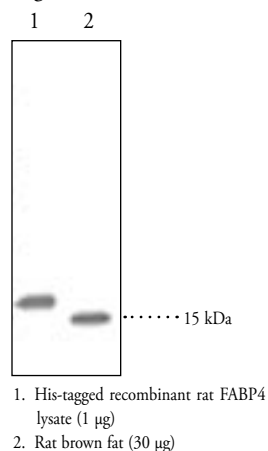
Item No. 10004944 • Lot No. XXXXXX

- Synonyms:** A-FABP, ALBP, aP2, Fatty Acid Binding Protein 4
- Contents:** This vial contains *lot specific* µg peptide affinity-purified IgG in *lot specific* µl TBS, pH 7.4, containing 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
- Host:** Rabbit
- Antigen:** This antibody was raised against human FABP4 amino acids 103-118 (TTIKRKREDDKLVVE). The antigen alignment with sequences from other species is as follows:
- | | |
|---------------|-------------------------------|
| Human FABP4 | T T I K R K R E D D K L V V E |
| Bovine FABP4 | T T I K R K l m D D K M V L E |
| Mouse FABP4 | T T I K R K R d g D K L V V E |
| Porcine FABP4 | T T I n R K i v D D K L V V E |
| Rabbit FABP4 | T T I K R K R E g D K L V V E |
| Rat FABP4 | T T I K R r R d g D K L V V E |
- Cross Reactivity:** (+) Human, mouse, and rat FABP4; other species not tested
- Stability:** ≥1 year at -20°C
- Applications:** Recommended dilutions for western blot (*lot specific* µg/ml) and immunocytochemistry (4 µg/ml). Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

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 mouse 3T3-L1 adipocytes and then from human adipocytes and was named 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}

Immunofluorescent staining procedure

1. Grow cells in 12 or 24 well plates until confluence.
2. Wash briefly with TBS, pH 7.4.
3. Fix the cells with 1% formaldehyde in TBS, pH 7.4, for 10 minutes.
4. Wash the cells with TBS containing 1% Triton-X 100 (TBSTX) 3 times, 10 minutes each.
5. Incubate the cells with 10% normal serum (from the same species in which the secondary antibody is raised) in TBSTX for 30 minutes.
6. Incubate the cells with the antibody (recommended starting concentration of 4 µg/ml. The optimal working condition should be determined by titration) for 1 hour.
7. Wash the cells with TBSTX 3 times, 10 minutes each.
8. Incubate the cells in the dark for 1 hour with a fluorochrome-conjugated secondary antibody at a concentration recommended by the provider.
9. Wash the cells with TBSTX 3 times, 10 minutes each.
10. Examine the staining under a fluorescent microscope with appropriate filter. Store the plate at 4°C in dark for later analysis.



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

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, 12/15/2011



References

1. Perrella, M.A., Pellacani, A., Layne, M.D., *et al.* *FASEB J.* **15**, 1774-1776 (2001).
2. Boord, J.B., Maeda, K., Makowski, L., *et al.* *Arterioscler. Thromb. Vasc. Biol.* **22**, 1686-1691 (2002).
3. Matarese, V. and Bernlohr, D.A. *J. Biol. Chem.* **263(28)**, 14544-14551 (1988).
4. Baxa, C.A., Sha, R.S., Buelt, M.K., *et al.* *Biochemistry* **28**, 8683-8690 (1989).
5. Distel, R.J., Robinson, G.S., and Spiegelman, B.M. *J. Biol. Chem.* **267(9)**, 5937-5941 (1992).
6. Melki, S.A. and Abumrad, N.A. *J. Lipid Res.* **34**, 1527-1534 (1993).
7. Hotamisligil, G.S., Johnson, R.S., Distel, R.J., *et al.* *Science* **274**, 1377-1379 (1996).
8. Uysal, K.T., Scheja, L., Wiesbrock, S.M., *et al.* *Endocrinology* **141(9)**, 3388-3396 (2000).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10004944

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