

PRODUCT INFORMATION



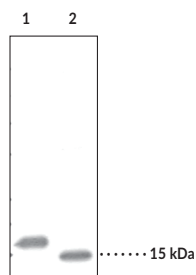
FABP4 Polyclonal Antibody

Item No. 10004944

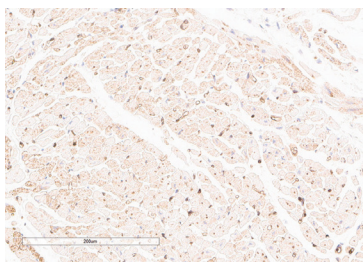
Overview and Properties

Contents:	This vial contains peptide affinity-purified IgG.
Synonyms:	Adipocyte FABP, A-FABP, ALBP, aP2, Fatty Acid Binding Protein 4
Immunogen:	Synthetic peptide from the C-terminal region of human FABP4
Cross Reactivity:	(-) FABP3 and FABP5
Species Reactivity:	(+) Human, Mouse, and Rat
Uniprot No.:	P15090
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Immunofluorescence (IF), Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution for IHC is 1:80 and the recommended starting dilution for IF and WB is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

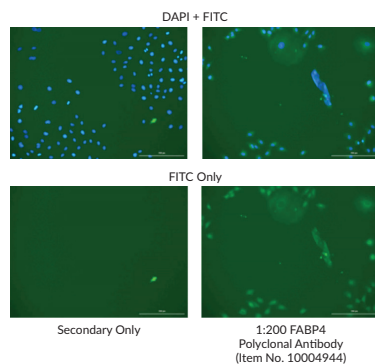
Images



Lane 1: His-tagged recombinant rat FABP4 lysate (1 µg)
Lane 2: Rat brown fat (30 µg)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human heart tissue after heat-induced antigen retrieval in pH 6.0 citrate buffer. After incubation with FABP4 Polyclonal Antibody (Item No. 10004944) at a 1:80 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).



Immunofluorescence analysis of paraformaldehyde-fixed A549 cells. After incubation with FABP4 Polyclonal Antibody (Item No. 10004944), at a 1:200 dilution (or negative control), cells were incubated with FITC-labeled anti-rabbit IgG (Item No. 10006588), followed by DAPI nuclear stain. Images show FITC alone or both fluorescence channels to highlight nuclear staining (where applicable).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/02/2019

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

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}

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).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM