

# PRODUCT INFORMATION



## FABP4 (human recombinant) FITC conjugated

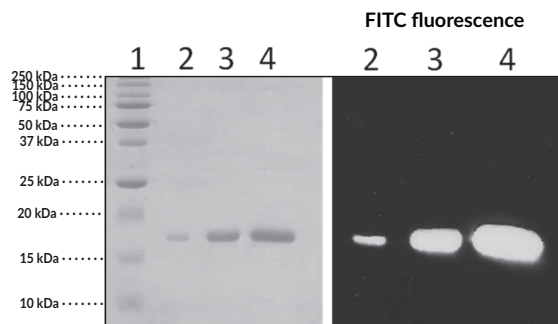
Item No. 9001068

### Overview and Properties

**Synonyms:** Adipocyte FABP, A-FABP, ALBP, aP2, Fatty Acid Binding Protein 4  
**Source:** Recombinant N-terminal His-tagged protein expressed in *E. coli*  
**Amino Acids:** 1-132 (full-length)  
**Molecular Weight:** 19 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥9 months  
**Purity:** ≥95% estimated by SDS-PAGE  
**Supplied in:** 50 mM sodium phosphate, pH 7.2, with 100 mM sodium chloride and 20% glycerol  
**Protein Concentration:** *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: FABP-FITC (1 µg)  
Lane 3: FABP-FITC (5 µg)  
Lane 4: FABP-FITC (10 µg)

Representative gel image shown; actual purity may

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, 02/07/2023

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

Fatty acid binding protein 4 (FABP4) is a member of the intracellular lipid-binding protein (iLBP) family.<sup>1</sup> It is composed of a  $\beta$ -barrel, which forms the lipid-binding site, and two  $\alpha$ -helices that form a lid on the  $\beta$ -barrel.<sup>2</sup> FABP4 is primarily expressed in adipocytes but has also been found in activated macrophages.<sup>3,4</sup> It is involved in lipid transport and metabolism, insulin resistance, and angiogenesis, and its expression is induced by long-chain fatty acids and peroxisome proliferator-activated receptor  $\gamma$  (PPAR $\gamma$ ) agonists.<sup>1,5-8</sup> Knockout of *Fabp4* prevents the development of atherosclerosis in *ApoE*<sup>-/-</sup> mice and increases body weight gain and insulin sensitivity in *ob/ob* mice.<sup>5,9</sup> Serum levels of FABP4 are increased in patients with colorectal cancer, and serum and synovial fluid levels of FABP4 are increased in patients with rheumatoid arthritis.<sup>7,10</sup> Cayman's FABP4 (human recombinant) FITC conjugated protein is conjugated to fluorescein isothiocyanate (FITC) and can be used for fluorescence applications.

## References

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(7)**, 1096-1116 (2002).
2. Reese-Wagoner, A., Thompson, J., and Banaszak, L. Structural properties of the adipocyte lipid binding protein. *Biochim. Biophys. Acta* **1441(2-3)**, 106-116 (1999).
3. 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(22)**, 8683-8690 (1989).
4. 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(10)**, 1774-1776 (2001).
5. 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).
6. 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(5291)**, 1377-1379 (1996).
7. Guo, D., Lin, C., Lu, Y., et al. FABP4 secreted by M1-polarized macrophages promotes synovitis and angiogenesis to exacerbate rheumatoid arthritis. *Bone Res.* **10(1)**, 45 (2022).
8. Bastie, C., Holst, D., Gaillard, D., et al. Expression of preoxisome proliferator-activated receptor PPAR $\delta$  promotes induction of PPAR $\gamma$  and adipocyte differentiation in 3T3C2 fibroblasts. *J. Biol. Chem.* **274(31)**, 21920-21925 (1999).
9. Lim, G.P., Yang, F., Chu, T., et al. Ibuprofen suppresses plaque pathology and inflammation in a mouse model for Alzheimer's disease. *J. Neurosci.* **20(15)**, 5709-5714 (2000).
10. Zhang, Y., Zhao, X., Deng, L., et al. High expression of FABP4 and FABP6 in patients with colorectal cancer. *World J. Surg. Oncol.* **17(1)**, 171 (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