

# Product Information



## FABP2 Polyclonal Antibody

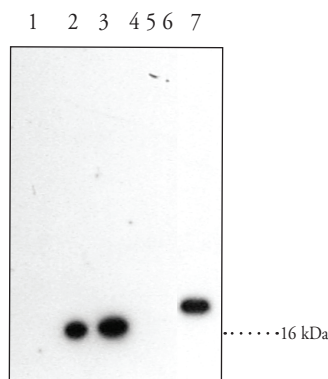
Catalog No. 10010019

- Contents:** This vial contains 300 µg peptide affinity-purified IgG in 500 µl TBS, pH 7.4, containing 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
- Synonyms:** Fatty Acid Binding Protein 2, Intestinal Fatty Acid Binding Protein, I-FABP
- Antigen:** Human FABP2 amino acids 33-40 (AHDNLKLT). This sequence is identical to that found in mouse and rat FABP2.
- Host:** Rabbit
- Cross Reactivity:** (+) human and rat FABP2; other species not tested but also expected to work with murine and bovine; (-) recombinant FABP1, 3, 4, and 5
- Stability:** ≥1 year at -20°C
- Applications:** The recommended starting dilution for western blot is 1:200 (3 µg/ml). Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Fatty acid binding protein 2 (FABP2) is one of nine known cytosolic FABPs ranging in size from 14-15 kDa containing 127-133 amino acids.<sup>1</sup> Members of this protein family exhibit high affinity for small lipophilic ligands and were named according to the tissue from which they were initially isolated. Studies suggest that FABPs are involved in the uptake and metabolism of fatty acids, in the maintenance of cellular membrane fatty acid levels, in intracellular trafficking of these substrates, in the regulation of specific enzymes of lipid metabolic pathways, and in the modulation of cell growth and differentiation.<sup>2</sup> FABP family members have highly conserved three dimensional structures and 22-73% amino acid sequence similarity. FABP2 is composed of ten antiparallel β strands that form a barrel that binds ligand in a bent conformation. FABP2 polymorphism has been suggested to be associated with gender specific obesity and increased risk of diabetes.<sup>1</sup>

### 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**, 1096-1116 (2002).
2. Massolini, G. and Calleri, E. Survey of binding properties of fatty acid-binding proteins chromatographic methods. *J. Chromatogr. B* **797**, 255-268 (2003).



1. Rat Recombinant FABP1 (0.4 µg)
2. Rat Recombinant FABP2 (0.025 µg)
3. Rat Recombinant FABP2 (0.050 µg)
4. Human Recombinant FABP3 (0.4 µg)
5. Murine Recombinant FABP4 (0.4 µg)
6. Murine Recombinant FABP5 (0.4 µg)
7. Human Duodenum Homogenate (30 µg)

### Related Products

FABP4 Polyclonal Antibody - Cat. No.10004944 • FABP4 (murine recombinant) - Cat. No.10005191 • FABP1 (rat recombinant) - Cat. No. 10005200 • FABP3 (human recombinant) - Cat. No. 10007432 • FABP5 (murine recombinant) - Cat. No. 10007433 • Arachidonic Acid-biotinamide - Cat. No. 10007466 • AFABP (human) EIA Kit - Cat. No. 10007614 • FABP2 (rat recombinant) - Cat. No. 10007938 • FABP4 (human recombinant) - Cat. No. 10009549 • FABP7 (human recombinant) - Cat. No. 10009551 • FABP4 (murine recombinant) Western Ready Control - Cat. No. 10009676 • FABP2 Blocking Peptide - Cat. No. 10010020 • FABP5 (human recombinant) - Cat. No. 10010364

### 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 NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG 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, 10/01/2008