

PRODUCT INFORMATION



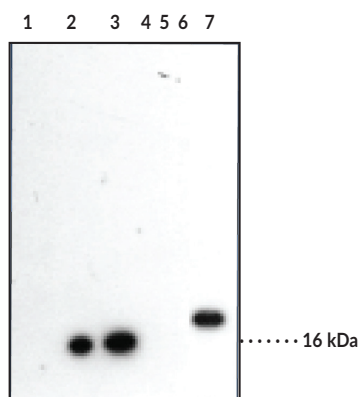
FABP2 Polyclonal Antibody

Item No. 10010019

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
Synonyms:	Fatty Acid Binding Protein 2, Intestinal-Fatty Acid Binding Protein, I-FABP
Immunogen:	Synthetic peptide from an internal region of human FABP2
Cross Reactivity:	(+) FABP2; (-) FABP1, FABP3, FABP4, FABP5, FABP7
Species Reactivity:	(+) Human, rat; other species not tested
Uniprot No.:	P12104
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	TBS, pH 7.4, with 50% glycerol, 0.1% BSA, and 0.02% sodium azide
Applications:	Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



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

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, 10/18/2023

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 2 (FABP2) is one of nine known cytosolic FABPs ranging in size from 14-15 kDa containing 127-133 amino acids.¹ 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 modulation of specific enzymes of lipid metabolic pathways, and in the regulation of cell growth and differentiation.² 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.¹ Cayman's FABP2 Polyclonal Antibody can be used for western blot applications. The antibody recognizes FABP2 at 15.207 kDa from human and rat samples.

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. Massolini, G. and Calleri, E. Survey of binding properties of fatty acid-binding proteins chromatographic methods. *J. Chromatogr. B. Analyt. Technol. Biomed. Life Sci.* **797(1-2)**, 255-268 (2003).

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