

PRODUCT INFORMATION



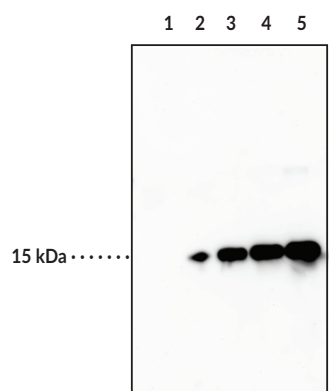
FABP3 Monoclonal Antibody (Clone CC68)

Item No. 10233

Overview and Properties

Contents:	This vial contains 200 µg of protein A-purified IgG2b
Synonyms:	H-FABP, Heart Fatty Acid Binding Protein, Fatty Acid Binding Protein 3, Cardiac FABP
Immunogen:	Synthetic peptide from human FABP3 amino acids 44-55
Cross Reactivity:	(-) Mouse FABP1, Mouse FABP2, and Mouse FABP4
Species Reactivity:	(+) Human and Rat FABP3; (-) Mouse FABP3
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2 with 50% glycerol and 0.02% sodium azide
Clone:	CC68
Host:	Mouse
Isotype:	IgG _{2b}
Applications:	Western blot; the recommended starting dilution 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: FABP4 WRC (4 µl)
Lane 2: FABP3 (10 ng)
Lane 3: FABP3 (100ng)
Lane 5: FABP3 (200 ng)

Description

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, 09/20/2018

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

Fatty acid binding protein 3 (FABP3) is one of nine known cytosolic FABPs ranging in size from 14-15 kDa containing 127-132 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.¹ 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 modulation of cell growth and differentiation.² FABP family members have highly conserved three dimensional structures and 22-73% amino acid sequence similarity. FABP3 is composed of ten antiparallel β strands that form a barrel and is the most widely distributed FABP. It is found in heart, skeletal and smooth muscle, mammary epithelial cells, aorta, distal tubules of the kidney, lung, brain, placenta, and ovary. FABP3 is a potential biomarker for myocardial injury, especially for early detection of acute myocardial infarction (AMI).¹

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* **797(1-2)**, 255-268 (2003).