

# PRODUCT INFORMATION



## FABP5 (mouse recombinant)

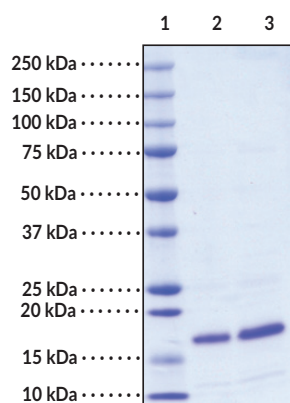
Item No. 10007433

### Overview and Properties

**Synonyms:** E-FABP, Fatty Acid Binding Protein 5, Epidermal-FABP  
**Source:** Recombinant N-terminal hexahistidine-tagged protein expressed in *E. coli*  
**Molecular Weight:** 19.3 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥2 years  
**Purity:** *batch specific* (≥80% estimated by SDS-PAGE)  
**Supplied in:** 50 mM sodium phosphate, pH 7.2, containing 20% glycerol and 150 mM sodium chloride  
**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  
Lane 2: 2 µg  
Lane 3: 4 µg

Representative gel image shown

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, 11/29/2021

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 5 (FABP5) is one of nine known cytosolic FABPs ranging in size from 14-15 kDa containing 127-132 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.<sup>1</sup> 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.<sup>2</sup> FABP family members have highly conserved three dimensional structures and 22-73% amino acid sequence similarity. FABP5 is composed of ten antiparallel  $\beta$  strands that form a barrel. FABP5 is expressed in skin, brain, lens, capillary, endothelium, and retina. Elevated level of FABP5 have been observed in prostate and bladder carcinoma.<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(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