

PRODUCT INFORMATION



FABP5 (human, recombinant)

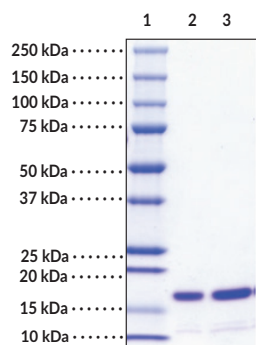
Item No. 10010364

Overview and Properties

Synonyms:	DA11 FABP, E-FABP, Epidermal-FABP, Fatty Acid Binding Protein 5, Keratinocyte FABP, Psoriasis-Associated FABP
Source:	Recombinant N-terminal hexahistidine-tagged protein expressed in <i>E. coli</i>
Amino Acids:	2-135 (Full length)
Uniprot No.:	Q01469
Molecular Weight:	18 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	<i>batch specific</i> (≥90% estimated by SDS-PAGE)
Supplied in:	50 mM sodium phosphate, pH 7.2, with 20% glycerol and 100 mM sodium chloride
Protein Concentration:	<i>batch specific</i> 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: FABP5 (1 μg)
Lane 3: FABP5 (2 μg)

Representative gel image shown; actual purity may vary between each batch.

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, 06/04/2024

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.¹ Members of this protein family exhibit high affinity for lipophilic ligands and were named according to the tissue from which they were initially isolated.¹ FABP5 is found in skin, brain, lens, capillary epithelium, and retina. Studies suggest that FABPs are involved in the uptake and metabolism of fatty acids, in the maintenance of cellular 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. The presence of one or more disulphide bridges is a unique feature of FABP5 that may be physiologically relevant.² Elevated levels of FABP5 have been observed in prostate and bladder carcinoma.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**, 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).

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