

FABP7 (IN) antibody (pAb)

Rabbit Anti-Human FABP7 (IN)

Instruction Manual

Catalog Number	PK-AB718-4259
Synonyms	FABP7 Antibody: Fatty acid binding protein 7, B-FABP, MRG, BLBP
Description	FABP7 was initially isolated from a human fetal brain cDNA library and whose mRNA was expressed in adult brain and muscle tissues at low levels. The protein encoded by this gene is a member of the fatty acid binding protein (FABPs) family, a group of small, highly conserved, cytoplasmic proteins that bind long-chain fatty acids and other hydrophobic ligands. FABPs are thought to play roles in fatty acid uptake, transport, and metabolism. FABP7 is a downstream gene of the Pax6 transcription factor and has been suggested to be essential for the maintenance of neuroepithelial cells during early cortical development. More recently, FABP7 was found to be frequently expressed in melanomas. Down-regulation of FABP7 through RNAi expression could reduce in vitro cell proliferation and Matrigel invasion, suggesting that FABP7 may be a potential target for the development of diagnostic and therapeutic tools.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	FABP7 antibody was raised in rabbits against a 17 amino acid peptide from near the center of human FABP7.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. FABP7 antibody can be used for detection of FABP7 by Western blot at 0.5 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL.
Images	Available upon request.
References	Shimizu F, Watanabe TK, Shinomiya H, et al. Isolation and expression of a cDNA for human brain fatty acid-binding protein (B-FABP). <i>Biochim. Biophys. Acta</i> 1997; 1354:24-8. Chmurzynska A. The multigene family of fatty acid-binding proteins (FABPs): function, structure and polymorphism. <i>J. Appl. Genet.</i> 2006; 47:39-48. Arai Y, Funatsu N, Numayama-Tsuruta K, et al. Role of FABP7, a downstream gene of pax6, in the maintenance of neuroepithelial cells during early embryonic development of the rat cortex. <i>J. Neurosci.</i> 2005; 25:9752-61. Goto Y, Matsuzaki Y, Kurihara S, et al. A new melanoma antigen fatty acid-binding protein 7, involved in proliferation and invasion, is a potential target for immunotherapy and molecular target therapy. <i>Cancer Res.</i> 2006; 66:4443-9.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4259P; FABP7 Peptide Cat.No. PK-AB718-1311; Human Breast Tissue Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.