

Instruction Manual

Catalog Number	PK-AB718-2309
Synonyms	FAIM Antibody: Fas apoptotic inhibitory molecule
Description	The susceptibility of primary splenic B cells to Fas-mediated apoptosis is regulated in a receptor-specific fashion. Terminal effectors of B cell Fas-resistance include the known anti-apoptotic proteins Bcl-xL, FLIP, and a recently identified protein termed FAIM. This molecule is broadly expressed in various tissues and exists in at least three isoforms. It is thought that resistance to Fas killing via increased expression of FAIM protects foreign antigen-specific B cells during interactions with FasL-bearing T cells whereas autoreactive B cells are deleted via Fas-dependent cytotoxicity. More recent results have indicated that FAIM interacts with both Trk and p75 neurotrophin receptor and may play a role in promoting neurite outgrowth in different neuronal systems by a mechanism involving the activation of NF-κB and the Ras-ERK pathway.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal FAIM antibody was raised against a 14 amino acid peptide from near the carboxy terminus of human FAIM (GenBank accession no. NP_060617).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually.
	FAIM antibody can be used for detection of FAIM by Western blot at 5 - 10 µg/mL.
Images	Available upon request.
References	Rothstein TL. Inducible resistance to Fas-mediated apoptosis in B cells. <i>Cell Res.</i> 2000; 10:245-66. Schneider TJ, Fischer GM, Donohoe TJ, et al. A novel gene coding for a Fas apoptosis inhibitory molecule (FAIM) isolated from inducibly Fas-resistant B lymphocytes. <i>J. Exp. Med.</i> 1999; 189:949-55. Sole C, Dolcet X, Segura MF, et al. The death receptor antagonist FAIM promotes neurite outgrowth by a mechanism that depends on ERK and NF-kappa B signaling. <i>J. Cell Biol.</i> 2004; 167:479-92. (06-01D)
Images	Available upon request.
Related Products	Human Spleen Tissue Lysate, Cat. No. PK-AB718-1306 FAIM2 Antibody, Cat. No. PK-AB718-2285 Blocking Peptide, Cat. No. PK-AB718-2309P FLIP Antibody (NT), Cat. No. PK-AB718-1159 NGFR Antibody, Cat. No. PK-AB718-3593
FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.	