

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-2285
<b>Synonyms</b>	FAIM2 Antibody: Lifeguard, Fas apoptotic inhibitory molecule2, FAIM2, NMP35
<b>Description</b>	Programmed cell death regulates a number of biological processes such as normal organism development, tissue homeostasis, and removal of damaged cells. Disruption of this process has been implicated in a variety of diseases such as cancer. LFG is a recently identified protein that can inhibit the apoptotic signal transduced by the Fas receptor but not from the related tumor necrosis factor- $\alpha$ death signal. In this respect, LFG is functionally similar to the anti-apoptotic proteins FAIM, FLIP and Bcl-xL. LFG, a seven membrane spanning protein, can bind the Fas receptor but does not regulate Fas expression or inhibit binding of FADD to Fas. LFG is widely distributed, but highly expressed in the hippocampus and other neural tissues. LFG was also identified as the neural membrane protein 35 (NMP35) and its expression is known to be regulated by the Phosphatidylinositol 3-kinase-Akt/PKB pathway.
<b>Quantity</b>	100 $\mu$ g
<b>Source / Host</b>	Rabbit
<b>Subclass</b>	Rabbit IgG
<b>Immunogen</b>	Rabbit polyclonal LFG antibody was raised against a 16 amino acid peptide from near the amino terminus of human LFG (GenBank accession no. AAF06327).
<b>Specific Species React.</b>	Human, Mouse, Rat
<b>Formulation</b>	Antibody is supplied in PBS containing 0.02% sodium azide.
<b>Stabilizer</b>	sodium azide
<b>Concentration</b>	1 mg/ml
<b>Reconstitution</b>	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 $\mu$ l 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.
<b>Purification Method</b>	Affinity chromatography purified via peptide column.
<b>Applications</b>	FAIM2 antibody can be used for detection of FAIM2 by Western blot at 0.5 - 1 $\mu$ g/mL. Antibody can also be used for immunohistochemistry starting at 5 $\mu$ g/mL.
<b>Storage</b>	E, WB, IHC   Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. 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.
<b>References</b>	Lockshin RA, Osborne B, and Zakeri Z. Cell death in the third millennium. Cell Death Differ. 2000; 7:2-7 Somia NV, Schmitt MJ, Vetter DE, et al. LFG: an antiapoptotic gene that provides protection from Fas-mediated cell death. Proc. Natl. Acad. Sci. USA 1999; 96:12667-72 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. J. Exp. Med. 1999; 189:949-55.

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

# LFG antibody (pAb)

Rabbit Anti-Human/Mouse/Rat LFG (FAIM2)



Schweitzer B, Taylor V, Welcher AA, et al. Neural membrane protein 35 (NMP35): a novel member of a gene family which is highly expressed in the adult nervous system. Mol. Cell. Neurosci. 1998; 11:260-73

## Related Products

Blocking Peptide, Cat. No. PK-AB718-2285P; EL4 Cell Lysate, Cat. No. PK-AB718-1287

FLIP Antibody (NT), Cat. No. PK-AB718-1159; NGFR Antibody, Cat. No. PK-AB718-3593; FAIM Antibody, Cat. No. PK-AB718-2309

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