

# AIFM3 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat AIFM3 (IN)

## Instruction Manual

Catalog Number	PK-AB718-4541
Synonyms	AIFM3 Antibody: Apoptosis-inducing factor mitochondrion-associated 3, AIFL
Description	Apoptosis, also known as programmed cell death, plays major roles in development and normal tissue turnover in addition to tumor formation. Recently a protein similar to the apoptosis-inducing factor (AIF) was cloned and designated AIFL (also known as AIFM3). AIFM3 is expressed ubiquitously and is predominantly localized to the inner membranes of mitochondria. Unlike AIF, AIFM3 does not translocate to the nucleus upon induction of apoptosis. However, overexpression of AIFM3, like AIF, induced cytochrome c release from the mitochondria, cleavage of caspase 3, and ultimately apoptosis, indicating AIFM3 induces apoptosis through caspase activation. Multiple isoforms of AIFM3 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	AIFM3 antibody was raised in rabbits against a 16 amino acid peptide from near the center of human AIFM3.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
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 <b>INote:</b> Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. AIFM3 antibody can be used for the detection of AIFM3 by Western blot at 1 - 2 µg/mL.
Images	Available upon request.
References	Jin Z and El Deiry WS. Overview of cell death signaling pathways. <i>Cancer Biol. Ther.</i> 2004; 4:139-63 Xie Q, Lin T, Zhang Y, et al. Molecular cloning and characterization of a human AIF-like gene with the ability to induce apoptosis. <i>J. Biol. Chem.</i> 2005; 280:19673-81.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4541P; AIFM3 Peptide Cat.No. PK-AB718-1303; Human Brain Tissue Lysate

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