

AIF (NT) antibody (pAb)

Rabbit Anti-Human Apoptosis Inducing Factor

Instruction Manual

Catalog Number	PK-AB718-2239
Synonyms	AIF Antibody:
Description	Apoptosis is characterized by several morphological nuclear changes including chromatin condensation and nuclear fragmentation. These changes are triggered by the activation of members of caspase family, caspase activated DNase, and several novel proteins. A novel gene, the product of which causes chromatin condensation and DNA fragmentation, was recently identified, cloned, and designated apoptosis inducing factor (AIF). Like the critical molecules, cytochrome c and caspase-9, in apoptosis, AIF localizes in mitochondria. AIF translocates to the nucleus when apoptosis is induced and induces mitochondria to release the apoptogenic proteins cytochrome c and caspase-9. AIF induces chromatin condensation and DNA fragmentation, which are the hallmarks of apoptosis, of the isolated nucleus and the nucleus in live cells by microinjection. AIF is highly conserved between human and mouse and widely expressed.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal AIF antibody was raised against a peptide corresponding to amino acids near the amino terminus of mature human AIF .
Purification Method	Ion exchange chromatography purified.
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, ICC, IF 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. AIF antibody can be used for detection of AIF by Western blot at 1 µg/mL. A 67 kDa band should be detected. Antibody can also be used for immunocytochemistry starting at 2 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Zamzami N, Kroemer G. Condensed matter in cell death. Nature 1999;401:127-8 Susin SA, Lorenzo HK, Zamzami N, et al. Molecular characterization of mitochondrial apoptosis-inducing factor. Nature 1999;397:441-6 (WD0500)
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2239P K562 Cell Lysate, Cat. No. PK-AB718-1204 Caspase-9 Antibody (IN1), Cat. No. PK-AB718-2071

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