

AIF (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Apoptosis Inducing Factor

Instruction Manual

| | |
|---------------------|---|
| Catalog Number | PK-AB718-2267 |
| Synonyms | AIF Antibody: AIF |
| 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 large scale DNA fragmentation, which are the hallmarks of apoptosis, of the isolated nucleus and the nucleus in live cells by microinjection and apoptosis stimuli. 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 517 to 531 of human AIF . This sequence is identical to those of mouse and rat AIF . |
| Purification Method | Ion exchange chromatography purified. |
| 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, IHC 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 immunohistochemistry starting at 10 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 Daugas E, Susin SA, Zamzami N, et al. Mitochondrio-nuclear translocation of AIF in apoptosis and necrosis. FASEB J 2000;14:729-39 (WD0500) |
| Images | Available upon request. |
| Related Products | Blocking Peptide, Cat. No. PK-AB718-2267P K562 Lysate, Cat. No. PK-AB718-1204 |

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