

XAF-1 (IN) antibody (pAb)

Rabbit Anti-Human XIAP Associated Factor 1 (Inhibitor of XIAP)

Instruction Manual

Catalog Number	PK-AB718-3409
Synonyms	XAF-1 Antibody: XAF-1
Description	XAF-1 binds to XIAP, an inhibitor of caspases-3, -7, and -9, and triggers its relocation from the cytosol to the nucleus. Overexpression of XAF-1 results in the neutralization of XIAP's ability to inhibit cell death. XAF-1 is normally expressed in all adult and fetal tissues but was found to be present in very low levels in a variety of cancer cell lines. In contrast, XIAP levels have been shown to be high in a majority of cell lines. Low XAF-1 and high basal expression of XIAP may therefore play a critical role in maintaining survival of cancer cell lines. Both IFN- α 2 and IFN- β can induce XAF-1 mRNA in all cells examined but induction of XAF-1 protein (as observed by immunoblot analysis) was seen only in cell lines sensitive to the apoptotic effects of IFNs.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal XAF-1 antibody was raised against a synthetic peptide corresponding to 16 amino acids near the middle of human XAF-1 (Genbank accession CAA68030).
Purification Method	Affinity chromatography purified via peptide column.
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 μ 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.
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, 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.
	XAF-1 antibody can be used for the detection of XAF-1 by Western blot at 1 - 4 μ g/mL. Multiple isoforms of XAF-1 exist and a doublet at ~30 kDa can be seen using anti-XAF-1 (IN). For immunofluorescence start at 20 μ g/mL.
Images	Available upon request.
References	Liston P, Fong W, Kelly NL, et al. Identification of XAF1 as an antagonist of XIAP anticaspase activity. <i>Nature Cell Biol.</i> 2001;3:128-33. Deveraux QL, Takahashi R, Savesan GS, and Reed JC. X-linked IAP is a direct inhibitor of cell-death proteases. <i>Nature</i> 1997;388:300-4. Fong WG, Liston P, Rajcan-Separovic E, et al. Expression and genetic analysis of XIAP-associated factor 1 (XAF1) in cancer cell lines. <i>Genomics</i> 2000;70:113-122. Leaman DW, Chawla-Sarkar M, Vyas K, et al. Identification of X-linked inhibitor of apoptosis associated factor-1 as an interferon-stimulated gene that augments TRAIL/Apo2L-induced apoptosis. <i>J. Biol. Chem.</i> 2002;277:28504-11.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3409P Human spleen Lysate, Cat. No. PK-AB718-1306 XAF-1 Antibody (CT), Cat. No. PK-AB718-3207 XIAP Antibody (CT), Cat. No. PK-AB718-3331

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