

# XIAP (CT) antibody (pAb)

## Rabbit Anti-Human/Mouse X-linked Inhibitor of Apoptosis Protein

### Instruction Manual

Catalog Number	PK-AB718-3331
Synonyms	XIAP Antibody:
Description	Apoptosis, or programmed cell death, is related to many diseases, such as cancer. Apoptosis is triggered by a variety of stimuli including members in the TNF family and can be prevented by the inhibitor of apoptosis (IAP) proteins. IAP proteins form a conserved gene family that binds to and inhibits cell death proteases. The X-chromosome linked inhibitor of apoptosis (XIAP) contains 3 baculoviral IAP repeat (BIR) motifs that are essential and sufficient for the binding and inhibition of caspases-3, -7, and -9. Upregulation of XIAP expression can protect cells from apoptosis induced by low level radiation; conversely, decreased XIAP expression by antisense targeting resulted in increased cell death following low level radiation. Two negative regulators, termed XAF-1 and Smac, can bind and inhibit XIAP activity.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal XIAP antibody was raised against a synthetic peptide corresponding to 13 amino acids at the C-terminus of human XIAP (Genbank accession No. NP_001158).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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, IF <span style="float: right;">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.</span> XIAP antibody can be used for the detection of XIAP by Western blot at 0.5 to 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2 µg/mL. For immunofluorescence start at 10 µg/mL.
Images	Available upon request.
References	Schimmer AD. Inhibitor of apoptosis proteins: translating basic knowledge into clinical practice. <i>Cancer Res.</i> 2004; 64:7183-90. Deveraux QL, Takahashi R, Savesan GS, et al. X-linked IAP is a direct inhibitor of cell-death proteases. <i>Nature</i> 1997; 388:300-4. Deveraux QL, Leo E, Stennicke HR, et al. Cleavage of human inhibitor of apoptosis protein XIAP results in fragments with distinct specificities for caspases. <i>EMBO J.</i> 1999; 18:5242-51. Holcik M, Yeh C, Korneluk RG, et al. Translational upregulation of X-linked inhibitor of apoptosis (XIAP) increases resistance to radiation induced cell death. <i>Oncogene</i> 2000; 19:4174-7.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3331P Smac/DIABLO Antibody (CT), Cat. No. PK-AB718-2409 XAF-1 Antibody (CT), Cat. No. PK-AB718-3207 Human kidney Lysate, Cat. No. PK-AB718-1305

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