

Rabbit Anti-Human Survivin (Inhibitor of Apoptosis)

Instruction Manual

Catalog Number	PK-AB718-2235
Synonyms	Survivin Antibody: TIAP
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 prevented by the inhibitor of apoptosis (IAP) proteins. IAP proteins form a conserved gene family that binds to and inhibits cell death proteases. A novel IAP protein was recently identified and designated survivin, apoptosis inhibitor 4 (API4), and TIAP. Survivin/TIAP interacted with the processed form of caspase-3 and inhibited its proteolytic activity. Survivin/TIAP is predominantly expressed in tissues of embryos, transformed cell lines, and many human cancers and lymphomas.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal survivin antibody was raised against a peptide corresponding to amino acids near the carboxy-terminus of human survivin.
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 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.
Images	Survivin antibody can be used for detection of survivin by Western blot at 1 µg/mL. An approximate 17 kDa band should be detected. Antibody can also be used for immunocytochemistry starting at 10 µg/mL. For immunofluorescence start at 10 µg/mL.
References	Ambrosini G, Adida C, Altieri DC. A novel anti-apoptosis gene, survivin, expressed in cancer and lymphoma. <i>Nat Med</i> 1997;3:917-21 Li F, Ambrosini G, Chu EY, Plescia J, Tognin S, Marchisio PC, Altieri DC . Control of apoptosis and mitotic spindle checkpoint by survivin. <i>Nature</i> 1998;396:580-4 Kobayashi K, Hatano M, Otaki M, Ogasawara T, Tokuhisa T. Expression of a murine homologue of the inhibitor of apoptosis protein is related to cell proliferation. <i>Proc Natl Acad Sci USA</i> 1999;96:1457-62 (WD0600)
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2235P Survivin Antibody (NT), Cat. No. PK-AB718-2233 MOLT4 Cell Lysate, Cat. No. PK-AB718-1205 mSurvivin Antibody (CT), Cat. No. PK-AB718-2237

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