

# ILPIP (IN) antibody (pAb)

## Rabbit Anti-Human/Mouse/Rat IAP-like Protein Interacting Protein

### Instruction Manual

Catalog Number	PK-AB718-3199
Synonyms	ILPIP Antibody:
Description	ILPIP (ILP-interacting protein) has recently been shown to interact with XIAP, a member of the IAP (Inhibitor of Apoptosis) protein family. Together, these two proteins synergistically activate the TAK1/JNK1 signal transduction pathway, which acts to protect against the interleukin-1 $\beta$ converting enzyme of Fas-induced apoptosis. This protein was independently isolated and characterized as polyploidy-associated protein kinase (PAPK), a member of the Ste20/germinal center kinase family that modulates cytoskeletal organization and cell survival, suggesting that its protective role may be a function of its kinase activity. Although two mRNA sequences have been isolated, only the shorter isoform has been observed.
Quantity	100 $\mu$ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal ILPIP antibody was raised against a synthetic peptide corresponding to amino acids near the center of human ILPIP (Genbank accession NP_061041).
Purification Method	Affinity chromatography purified via peptide column.
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 $\mu$ 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, 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. ILPIP antibody can be used for the detection of ILPIP by Western blot at 1 to 2 $\mu$ g/mL. Antibody can also be used for immunohistochemistry starting at 10 $\mu$ g/mL.
Images	Available upon request.
References	Sanna MG, Correia JS, Luo Y, et al. (2002) ILPIP, a novel anti-apoptotic protein that enhances XIAP-mediated activation of JNK and protection of apoptosis. J. Biol. Chem. 277:30454-62. Sanna MG, Duckett CS, Richter BW, et al. (1998) Selective activation of JNK1 is necessary for the anti-apoptotic activity of hILP. Proc. Natl. Acad. Sci USA 95:6015-20. Nishigaki K, Thompson D, Yugawa T, et al. (2003) Identification and characterization of a novel Ste20/germinal center kinase-related kinase, polyploidy-associated protein kinase. J. Biol. Chem. 278:13520-30.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3199P Human brain Lysate, Cat. No. PK-AB718-1303 ILPIP Antibody (CT), Cat. No. PK-AB718-3201 XIAP Antibody (CT), Cat. No. PK-AB718-3331 TAK1 Antibody (NT), Cat. No. PK-AB718-3385

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