

## Rabbit Anti-Human/Mouse/Rat IAP-like Protein 2 (Member in IAP Family)

### Instruction Manual

Catalog Number	PK-AB718-3017
Synonyms	ILP-2 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 prevented by the inhibitor of apoptosis (IAP) proteins. IAP proteins form a conserved gene family including IAP, XIAP/ILP-1/MIHA, and Livin/KIAP that bind to and inhibits specific proteases. A novel member in the IAP protein family was recently identified and designated ILP-2 for IAP-like protein-2. ILP-2 has high homology to ILP-1, but is encoded by a distinct gene that is solely expressed in testis of tested normal human tissues. ILP-2, unlike ILP-1, has no inhibitory effect on Fas and TNF induced apoptosis, but potently inhibits apoptosis induced by overexpression of Bax or by coexpression of caspase-9 with Apaf-1. ILP-2 interacts with the processed caspase-9. These results suggest that ILP-2 is a novel IAP family member with restricted specificity for caspase-9.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal ILP-2 antibody was raised with against a synthetic peptide corresponding to amino acids near the amino terminus of human ILP-2 (Genbank accession No. Q96P09).
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 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 <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>
Images	Available upon request.
References	Richter BW, Mir SS, Eiben LJ, et al. Molecular cloning of ILP-2, a novel member of the inhibitor of apoptosis protein family. Mol. Cell Biol. 2001;21(13):4292-301.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3017P HepG2 Cell Lysate, Cat. No. PK-AB718-1211 Apaf-1 Antibody (NT), Cat. No. PK-AB718-2013 Caspase 9 Antibody (IN1), Cat. No. PK-AB718-2071 Livin Antibody, Cat. No. PK-AB718-2505 cIAP Antibody (CT), Cat. No. PK-AB718-3325; XIAP Antibody (CT), Cat. No. PK-AB718-3331

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