

## Rabbit Anti-Human Ipaf (ICE Protease Activating Factor)

### Instruction Manual

<b>Catalog Number</b>	PK-AB718-3107
<b>Synonyms</b>	Ipaf Antibody: Ipaf, CARD12, CLANA
<b>Description</b>	Apoptosis is related to many diseases and induced by a family of cell death receptors and their ligands. Cell death signals are transduced by death domain containing adaptor molecules and proteases including several members of the caspase family. Another family of proteins that functions as a critical regulator of apoptosis and NF-κB signaling pathways is the CED-4/Apaf-1 (apoptosis protein activating factor-1) protein family. Ipaf (ICE protease activating factor) is a CED-4/Apaf-1 family member that activates caspase-1/ICE and can induce apoptosis in human cells in a caspase-1 dependent manner. Ipaf and caspase-1 are thought to interact with each other through the association of the Ipaf amino-terminal CARD (caspase recruitment domain) and amino-terminal CARD of caspase-1.
<b>Quantity</b>	100 µg
<b>Source / Host</b>	Rabbit
<b>Immunogen</b>	Rabbit polyclonal Ipaf antibody was raised against a synthetic peptide corresponding to amino acids near the C-terminus of human Ipaf (Genbank accession NP_067032).
<b>Purification Method</b>	Affinity chromatography purified via peptide column.
<b>Clone / IgG Subtype</b>	Polyclonal antibody
<b>Species Reactivity</b>	Human
<b>Specificity</b>	
<b>Formulation</b>	Antibody is supplied in PBS containing 0.02% sodium azide.
<b>Reconstitution</b>	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.
<b>Storage &amp; Stability</b>	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.
<b>Applications</b>	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> Ipaf polyclonal antibody can be used for the detection of Ipaf by Western blot at 1 and 2 µg/mL. A 110 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 10 µg/mL. For immunofluorescence start at 10 µg/mL.
<b>Images</b>	Available upon request.
<b>References</b>	Li P, Nijhawan D, Budihardjo I, et al. Cytochrome c and dATP-dependent formation of Apaf-1/caspase-9 complex initiates an apoptotic protease cascade. Cell 1997; 91:479-89. Poyet J-L, Srinivaula SM, Tnani M, et al. Identification of Ipaf, a human caspase-1-activating protein related to Apaf-1. J. Biol. Chem. 2001; 276:28309-13. Geddes BJ, Wang L, Huang WJ, et al. Human CARD12 is a novel CED4/Apaf-1 family member that induces apoptosis. Biochem. Biophys. Res. Commun. 2001; 284:77-82.
<b>Images</b>	Available upon request.
<b>Related Products</b>	Blocking Peptide, Cat. No. PK-AB718-3107P Human PBL Lysate, Cat. No. PK-AB718-1371 Apaf Antibody (CT), Cat. No. PK-AB718-2015 Apaf Antibody (NT), Cat. No. PK-AB718-2013.

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