

Rabbit Anti-Human NALP1 (Caspase Interacting Protein)

Instruction Manual

Catalog Number	PK-AB718-3037
Synonyms	NALP1 Antibody: CARD7, DEFCAP, NACHT domain- leucine-rich-repeat and PYD-containing protein 1
Description	NALP1 belongs to a family of cytoplasmic proteins that have been implicated in cell responses to apoptotic and inflammatory stimuli. It contains a caspase recruitment domain (CARD) in addition to a NACHT domain and leucine rich repeat (LRR). This protein interacts strongly with caspase 2 and weakly with caspase 9, as well as with Apaf-1. Transient overexpression of this gene in cultured cells was sufficient to induce apoptosis. NALP1 mRNA and protein expression levels in myeloid leukemia cells are increased following CREB (cAMP-response element binding protein) activation, suggesting that NALP1 may contribute to modulate the response of these cells to pro-inflammatory stimuli. Multiple alternatively spliced transcript variants encoding distinct isoforms have been found for NALP1.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal NALP1 antibody was raised against a peptide corresponding to 13 amino acids near the carboxy-terminus of human NALP1 (GenBank accession no. AAG30288).
Purification Method	Immunoaffinity chromatography purified IgG.
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. NALP1 antibody can be used for detection of NALP1 by Western blot at 1 to 4 µg/mL. Antibody can also be used for immunocytochemistry starting at 10 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	NALP1 antibody can be stored at 4°C, stable for one year. Tschopp J, Martinon F, and Burns K. NALPs: a novel protein family involved in inflammation. <i>Nat. Rev. Mol. Cell Biol.</i> 2003; 4:95-104. Chu ZL, Pio F, Xie Z, et al. A novel enhancer of the Apaf1 apoptosome involved in cytochrome c-dependent caspase activation and apoptosis. <i>J. Biol. Chem.</i> 2001; 276:9239-45. Hlaing T, Guo RF, Dille KA, et al. Molecular cloning and characterization of DEFCAP-L and -S, two isoforms of a novel member of the Ced-4 family of apoptosis proteins. <i>J. Biol. Chem.</i> 2001; 276:9230-8. Sanz C, Calasanz MJ, Andreu E, et al. NALP1 is a transcriptional target for cAMP-response-element-binding protein (CREB) in myeloid leukemia cells. <i>Biochem. J.</i> 2004; 384:281-6.
Images	Available upon request.
Related Products	U937 Lysate, Cat. No. PK-AB718-1215 K562 Lysate, Cat. No. PK-AB718-1204 NALP1 Blocking Peptide, Cat. No. PK-AB718-3037P
FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.	