

# NALP2 / PYPAF2 antibody (pAb)

## Rabbit Anti-Human/Mouse/Rat NALP2 (Caspase Interacting Protein)

### Instruction Manual

Catalog Number	PK-AB718-3027
Synonyms	NALP2 Antibody: NACHT domain-, leucine-rich-repeat-, and PYD-containing protein 2, PYRIN-containing APAF1-like protein2, PAN1, PYPAF2
Description	NALP2 belongs to a family of cytoplasmic proteins that have been implicated in cell responses to apoptotic and inflammatory stimuli. Unlike the prototypical NALP protein NALP1, NALP2 only contains a NACHT domain, leucine rich repeat (LRR), and pyrin-containing domain (PYD). This protein interacts with the adapter protein ASC in addition to CARD8 and caspase-1 to form an inflammasome with high proIL-1 $\beta$ -processing activity and is thought to function as a modulator of NF- $\kappa$ B and procaspase-1 activation in macrophages. It has also been suggested that NALP2, in addition to other NALP family members, can act as innate sensors of pathogens in a manner similar to the toll-like receptors (TLRs). At least two alternatively spliced transcript variants encoding distinct isoforms have been found for NALP2.
Quantity	100 $\mu$ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal NALP2 antibody was raised against a peptide corresponding to 12 amino acids near the carboxyterminus of human NALP2 (GenBank accession no. NP_060322).
Purification Method	Immunoaffinity chromatography purified IgG
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 I Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. NALP2 antibody can be used for detection of NALP2 by Western blot at 1 to 2 $\mu$ g/mL.
Images	Available upon request.
References	Tschopp J, Martinon F, and Burns K. NALPs: a novel protein family involved in inflammation. Nat. Rev. Mol. Cell Biol. 2003; 4:95-104. Bruey JM, Bruey-Sedano N, Newman R, et al. PAN1/NALP2/PYPAF2, an inducible inflammatory mediator that regulates NF- $\kappa$ B and caspase-1 activation in macrophages. J. Biol. Chem. 2004; 279:51897-907. Agostini L, Martinon F, Burns K, et al. NALP3 forms an IL-1 $\beta$ -processing inflammasome with increased activity in Muckle-Wells autoinflammatory disorder. Immunity 2004; 20:319-25. Martinon F and Tschopp J. NLRs join TLRs as innate sensors of pathogens. TRENDS Imm. 2005; 26:447-54.
Images	Available upon request.
Related Products	PC-3 Cell Lysate, Cat. No. PK-AB718-1216; Blocking Peptide, Cat. No. PK-AB718-3027P NALP1 Antibody, Cat. No. PK-AB718-3037; ASC Antibody (CT), Cat. No. PK-AB718-2287 CARD8 Antibody, Cat. No. PK-AB718-3187; Caspase 1 Antibody (CT), Cat. No. PK-AB718-3459

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