

NAIP (CT) antibody (pAb)

Rabbit Anti-Human Neuronal Apoptosis Inhibitor Protein

Instruction Manual

Catalog Number	PK-AB718-3315
Synonyms	NAIP Antibody:
Description	Neuronal apoptosis inhibitor protein (NAIP) was the first human inhibitor of apoptosis protein (IAP) identified and was discovered by its association with the neurodegenerative disorder spinal muscular atrophy. Members of the IAP family contain one to three copies of an approximately 70 amino acid motif termed baculovirus IAP repeat (BIR); these BIRs promote protein-protein interactions with various caspases such as caspase-3, -7, and -9 as well as members of the TRAF family of signal molecules. Unlike other IAPs however, NAIP requires ATP to bind caspase-9 and is not inhibited by the IAP-inhibiting molecule Smac/DIABLO, suggesting that NAIP is unique among the IAPs in its regulation of its activity. Finally, although only one human NAIP protein has been identified, other shorter NAIP mRNA transcripts have been reported.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal NAIP antibody was raised against a synthetic peptide corresponding to 13 amino acids at the C-terminus of human NAIP (Genbank accession No. AAC52047).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	Minor, lower molecular weight bands may represent alternately spliced forms.
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. NAIP antibody can be used for the detection of NAIP by Western blot at 0.5 to 2 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	Roy N, Mahadevan MS, McLean M, et al. The gene for neuronal apoptosis inhibitory protein is partially deleted in individuals with spinal muscular atrophy. Cell 1995; 80:167-78. Liston P, Fong WG, Korneluk RG. The inhibitors of apoptosis: there is more to life than Bcl2. Oncogene 2003; 22:8568-80. Verhagen AM, Coulson EJ, and Vaux DL. Inhibitor of apoptosis proteins and their relatives: IAPs and other BIRPs. Genome Biol. 2001; 2:reviews3009.1–reviews3009.10. Davoodi J, Lin L, Kelly J, et al. Neuronal apoptosis-inhibitory protein does not interact with Smac and requires ATP to bind caspase-9. J. Biol. Chem. 2004; 279:40622-8.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3315P Smac/DIABLO Antibody, Cat. No. PK-AB718-2409 Casp-9 Antibody (IN1), Cat. No. PK-AB718-2071 PC-3 Cell Lysate, Cat. No. PK-AB718-1216

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