

NIPSNAP (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat NIPSNAP (CT)

Instruction Manual

Catalog Number	PK-AB718-4535
Synonyms	NIPSNAP Antibody: Non-neuronal SNAP25-like protein, NIPSNAP1
Description	NIPSNAP is a member of an evolutionarily well conserved gene family and has a strong sequence similarity to the central portion of a protein encoded by <i>C. elegans</i> chromosome III between a 4-nitrophenylphosphatase (NIP) domain and non-neuronal SNAP25-like protein. Recent studies have indicated that NIPSNAP is involved in the regulation of the Ca ²⁺ -selective transient receptor potential vanilloid channel 6 (TRPV6). NIPSNAP1 associates with TRPV6 at the plasma membrane and inhibits TRPV6 currents. Other studies show that NIPSNAP's expression is reduced in the phenylketonuria (PKU) mouse brain, suggesting that NIPSNAP may play a role in memory.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	NIPSNAP antibody was raised in rabbits against an 18 amino acid peptide from near the carboxy terminus of human NIPSNAP.
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, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. NIPSNAP antibody can be used for the detection of NIPSNAP by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Seroussi E, Pan H-Q, Kedra D, et al. Characterization of the human NIPSNAP1 gene from 22q12: a member of a novel gene family. <i>Gene</i> 1998; 212:13-20. Schoeber JP, Topala CN, Lee KP, et al. Identification of Nipsnap1 as a novel auxiliary protein inhibiting TRPV6 activity. <i>Pflugers Arch.</i> 2008; epub. Surendran S, Tyring SK and Matalon R. Expression of calpastatin, minopontin, NIPSNAP1, rabaptin-5 and neuronatin in the phenylketonuria (PKU) mouse brain: possible role on cognitive defect seen in PKU. <i>Neurochem. Int.</i> 2005; 46:595-9.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4535P; NIPSNAP Peptide Cat.No. PK-AB718-1303; Human Brain Tissue Lysate

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