

NPTX2 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat NPTX2 (Neuronal pentraxin 2, NP2, NARP)

Instruction Manual

Catalog Number	PK-AB718-4573
Synonyms	NPTX2 Antibody: Neuronal pentraxin 2, NP2, NP-II, neuronal activity-regulated protein, NARP
Description	Neuronal pentraxin (NPTX) 2 was initially identified as a member of the pentraxin family having high homology to NPTX1. Unlike NPTX1 however, NPTX2 is expressed in testes, pancreas, liver, heart, skeletal muscle as well as brain. NPTX2 possesses the lectin properties common to the pentraxin family, promotes neurite outgrowth, and is rapidly regulated by neuronal activity. NPTX1 and NPTX2 form heterocomplexes that contribute to both activity-independent and -dependent excitatory synaptogenesis. Recently, NPTX2 has been found to be highly upregulated in Parkinsonian substantia nigra and localizes to Lewy bodies and Lewy neurites in sporadic Parkinson's disease (PD), suggesting that it is involved in the pathway dysregulation that underlies PD. This NPTX2 antibody is predicted to be specific to NPTX2 and not recognize NPTX1.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	NPTX2 antibody was raised in rabbits against a 16 amino acid peptide near the center of the human NPTX2.
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. NPTX2 antibody can be used for detection of NPTX2 by Western blot at 0.5 and 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Hsu YC and Perin MS. Human neuronal pentraxin II (NPTX2): conservation, genomic structure, and chromosomal localization. <i>Genomics</i> 1995; 28:220-7. Tsu CC, Copeland GG, Gilbert DJ, et al. Narp, a novel member of the pentraxin family, promotes neurite outgrowth and is dynamically regulated by neuronal activity. <i>J. Neurosci.</i> 1996; 16:2463-78. Xu D, Hopf C, Reddy R, et al. Narp and NP1 form heterocomplexes that function in developmental and activity-dependent synaptic plasticity. <i>Neuron</i> 2003; 39:513-28. Moran LB, Hickey L, Michael GJ, et al. Neuronal pentraxin II is highly upregulated in Parkinson's disease and a novel component of Lewy bodies. <i>Acta Neuropathol.</i> 2008; 115:471-8.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4573P; NPTX2 Peptide Cat. No. PK-AB718-1403; Mouse Brain Tissue Lysate

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