

Nicastrin (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Nicastrin (Anterior Pharynx Defective 2, APH-2)

Instruction Manual

Catalog Number	PK-AB718-3983
Synonyms	Nicastrin Antibody: Anterior pharynx defective 2, APH-2
Description	Nicastrin, in addition to presenilin, PEN2, and APH-1 forms the γ -secretase protein complex, a membrane-bound aspartyl protease that can cleave certain proteins at peptide bonds buried within the hydrophobic environment of the lipid bilayer. This cleavage is responsible for a key step in signaling from several cell-surface receptors and is thought to be required for the generation of the neurotoxic amyloid peptides that are central to the pathogenesis of Alzheimer's disease. Like the tumor necrosis factor- α -converting enzyme (TACE) and the b-site cleavage enzyme (BACE) protease families, γ -secretase will cleave the amyloid precursor protein (APP), but within the intramembrane region of APP, resulting in either the non-toxic p3 (from the α and γ cleavage site) or the toxic A β amyloid peptide (from the β and γ cleavage site). It is thought that accumulation of the A β peptide is the precursor to Alzheimer's disease. Nicastrin is also thought to be involved in cell proliferation and signaling.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Nicastrin antibody was raised against a 17 amino acid peptide from near the carboxy terminus of human Nicastrin.
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 μ 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, IHC, IF Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Nicastrin antibody can be used for detection of Nicastrin by Western blot at 0.5 - 2 μ g/mL. Antibody can also be used for immunohistochemistry starting at 5 μ g/mL. For immunofluorescence start at 20 μ g/mL.
Images	Available upon request.
References	Weihofen A and Martoglio B. Intramembrane-cleaving proteases: controlled liberation of proteins and bioactive peptides. Trends Cell Biol. 2003; 13:71-8. Periz G and Fortini ME. Functional reconstitution of γ -secretase through coordinated expression of presenilin, Nicastrin, aph-1, and pen-2. J. Neurosci. Res. 2004; 77:309-22. Selkoe DJ. The cell biology of b-amyloid precursor protein and presenilin in Alzheimer's disease. Trends Cell Biol. 1998; 8:447-53. Nguyen V, Hawkins C, Bergeron C, et al. Loss of nicastrin elicits an apoptotic phenotype in mouse embryos. Brain Res. 2006; 1086:76-84.
Images	Available upon request.
Related Products	Nicastrin Peptide, Catalog No. PK-AB718-3983P; Brain Lysate, Cat. No. PK-AB718-1403; Nicastrin Antibody, Cat. No. PK-AB718-3985; PEN2 Antibody, Catalog No. PK-AB718-3979 APP Antibody, Cat. No. PK-AB718-2136; APP Antibody, Cat. No. PK-AB718-2133; APH1 Antibody, Cat. No. PK-AB718-4001; BACE Antibody, Cat. No. PK-AB718-2253; BACE2 Antibody, Cat. No. PK-AB718-2249

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