

Slitrk5 (IN2) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Slitrk5 (IN2)

Instruction Manual

Catalog Number	PK-AB718-4477
Synonyms	Slitrk5 Antibody: SLIT and NTRK-like family 5
Description	SLIT and NTRK-like family 5 (Slitrk5) is a member a protein family consisting of six homologous transmembrane proteins (Slitrk1-6) that share two conserved leucine-rich repeat domains in the extracellular domain and have significant homology to Slit, a secreted axonal growth-controlling protein. These proteins are also homologous to trk neurotrophin receptors in their intracellular domains. Expression of Slitrk proteins is highly restricted to neural and brain tumor tissues, but varies within the protein family. Like every other Slitrk protein except Slitrk1, overexpression of Slitrk5 inhibited neurite outgrowth in cultured neurons, suggesting that these proteins are involved in the control of neurite outgrowth. This antibody is predicted to have no cross-reactivity to other Slitrk proteins.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Slitrk5 antibody was raised in rabbits against a 15 amino acid peptide from near the center of human Slitrk5.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	This antibody is predicted to have no cross-reactivity to other Slitrk proteins.
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. Slitrk5 antibody can be used for detection of Slitrk5 by Western blot at 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	Aruga J and Mikoshiba K. Identification and characterization of Slitrk, a novel transmembrane protein family controlling neurite outgrowth. Mol. Cell Neurosci. 2003; 24:117-29. Aruga J, Yokota N, and Mikoshiba K. Human SLITRK family genes: genomic organization and expression profiling in normal and brain tumor tissue. Gene 2003; 315:87-94.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4477P; Slitrk5 Peptide Cat.No. PK-AB718-1212; 3T3 Cell Lysate

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