

BRAL1 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat BRAL1

Instruction Manual

Catalog Number	PK-AB718-4501
Synonyms	BRAL1 Antibody: Brain link protein-1, hyaluronan and proteoglycan link protein 2, HAPLN2
Description	BRAL1 is a member a superfamily consisting of several highly homologous hyaluronan and proteoglycan binding link proteins. BRAL1 is predominantly expressed in brain tissue and spinal cord. Like other members in the link-module superfamily, BRAL1 contains an immunoglobulin-like fold and two proteoglycan tandem repeats (PTRs). Its mRNA expression pattern is similar to other lectican proteoglycans, suggesting that BRAL1 may act to stabilize the binding between the extracellular matrix molecule hyaluronan and brevican. Immunostaining of mouse brain showed BRAL1 expression at P20 in the white matter of the developing cerebellum and in myelinated fiber tracts in the adult brain, suggesting that expression starts when axonal myelination occurs.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	BRAL1 antibody was raised in rabbits against a 12 amino acid peptide from near the center of human BRAL1.
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. BRAL1 antibody can be used for the detection of BRAL1 by Western blot at 1 - 2 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	Spicer AP, Joo A, and Bowling Jr. RA. A hyaluronan binding link protein gene family whose members are physically linked adjacent to chondroitin sulfate proteoglycan core protein genes. <i>J. Biol. Chem.</i> 2003; 278:21083-91. Hirakawa S, Oohashi T, Su W-D, et al. The brain link protein-1 (BRAL1): cDNA cloning, genomic structure, and characterization as a novel link protein expressed in adult brain. <i>Biochem. Biophys. Res. Comm.</i> 2000; 276:982-9. Oohashi T, Hirakawa S, Bekku Y, et al. Bral1, a brain-specific link protein, colocalizing with the versican V2 isoform at the nodes of Ranvier in developing and adult mouse central nervous systems. <i>Mol. Cell Neurosci.</i> 2002; 19:43-57.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4501P; BRAL1 Peptide Cat.No. PK-AB718-1303; Human Brain Tissue Lysate

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