

BAP3 (NT) antibody (pAb)

Rabbit Anti-Human BAP3 (NT)

Instruction Manual

Catalog Number	PK-AB718-4503
Synonyms	BAP3 Antibody: BAI1-associated protein 3, BAIAP3
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	BAP3 antibody was raised in rabbits against a 12 amino acid peptide from near the amino terminus of human BAP3.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
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. BAP3 antibody can be used for the detection of BAP3 by Western blot at 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	Shiratsuchi T, Oda K, Nishimori H, et al. Cloning and characterization of BAP (BAI-associated protein 3), a C2 domain-containing protein that interacts with BAI1. <i>Biochem. Biophys. Res. Comm.</i> 1998; 251:158-65. Palmer RE, Lee SB, Wong JC, et al. Induction of BAIAP3 by the EWS-WT1 chimeric fusion implicates regulated exocytosis in tumorigenesis. <i>Cancer Cell</i> 2002; 2:497-505.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4503P; BAP3 Peptide Cat.No. PK-AB718-1220; SK-N-SH Cell Lysate

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