

# BAP3 (NT2) antibody (pAb)

## Rabbit Anti-Human BAP3 (NT2)

### Instruction Manual

Catalog Number	PK-AB718-4505
Synonyms	BAP3 Antibody: BAI1-associated protein 3, BAIAP3
Description	BAP3 was initially identified through interaction in a yeast two-hybrid system with the brain-specific angiogenesis inhibitor 1, a p53-target gene that encodes a seven-span transmembrane protein member of the secretin receptor family. BAP3 is predominantly expressed in the brain and possess high homology with Munc13 and synaptotagmin, suggesting that BAP3 may play a role in regulating neurotransmitter release. Recent experiments have shown that BAP3 is induced in certain tumors such as desmoplastic small round cell tumor. Ectopic expression of BAP3 in tumor cells dramatically enhances growth in low serum conditions and colony formation in soft agar, suggesting that the regulated exocytotic pathway may play a role in cancer cell proliferation. BAP3 is known to have two isoforms; this BAP3 antibody will recognize only isoform 2. Lower molecular weight bands may represent cleavage or degradation products.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	BAP3 antibody was raised in rabbits against a 13 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	BAP3 is known to have two isoforms; this BAP3 antibody will recognize only isoform 2. Lower molecular weight bands may represent cleavage or degradation products.
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 Note: 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 1 and 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 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-AB718-4505P; BAP3 Peptide Cat.No. PK-AB718-1220; SK-N-SH Cell Lysate

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