

op18 (CT) antibody (pAb)

Rabbit Anti-Human/Mouse op18 (CT)

Instruction Manual

Catalog Number	PK-AB718-4235
Synonyms	op18 Antibody: oncoprotein 18, STMN1, Stathmin 1, Lag, LAP18
Description	Op18 belongs to the stathmin family of genes and encodes a ubiquitous cytosolic phosphoprotein that may function as an intracellular relay integrating several signaling pathways such as those involved in cell proliferation and differentiation. Op18 has also been shown to be involved in the regulation of the microtubule filament system by destabilizing microtubules, thereby preventing assembly and promoting the disassembly of microtubules. More recently, op18 has been implicated as a potential target of the ASK1-p38 MAP kinase cascade, suggesting that the ASK1-p38 cascade may regulate microtubule dynamics through op18. Op18 is highly expressed in a wide variety of human malignancies, including leukemia, prostate cancer, ovarian carcinoma, and breast carcinoma, suggesting that op18 may be an ideal target for anti-cancer therapeutics.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Op18 antibody was raised in rabbits against a 20 amino acid peptide from near the carboxy terminus of human op18.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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. Op18 antibody can be used for detection of op18 by Western blot at 0.5 - 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	Curmi PA, Gavet O, Charbaut E, et al. Stathmin and its phosphoprotein family: general properties, biochemical and functional interaction with tubulin. Cell Structure and Function 1999; 24:345-57. Belmont LD and Mitchison TJ. Identification of a protein that interacts with tubulin dimers and increases catastrophe rate of microtubules. Cell 1996; 84:623-31. Mizumura K, Takeda K, Hashimoto S, et al. Identification of op18/stathmin as a potential target of ASK1-p38 MAP kinase cascade. J. Cell Physiol. 2006; 206:363-70. Mistry SJ and Atweh GF. Role of stathmin in the regulation of the mitotic spindle. Mount Sinai J. Med. 2002; 69:299-304.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4235P; op18 Peptide Cat.No. PK-AB718-1287; EL4 Lysate

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