

PIST / GOPC (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat PIST (Autophagy Related Protein)

Instruction Manual

Catalog Number	PK-AB718-3631
Synonyms	PIST Antibody: Golgi associated PDZ and coiled-coil motif containing, GOPC
Description	Autophagy, the process of bulk degradation of cellular proteins through an autophagosomal-lysosomal pathway is important for normal growth control and may be defective in tumor cells. It is involved in the preservation of cellular nutrients under starvation conditions as well as the normal turnover of cytosolic components and is negatively regulated by TOR (Target of rapamycin). PIST, a PDZ-containing protein, was discovered in a yeast two-hybrid system as a binding partner to Beclin-1, a Bcl-2-interacting protein homologous to the yeast autophagy gene <i>apg6</i> . Experiments with mutant PIST proteins lacking the PDZ domain showed that PIST interaction with Beclin-1 through its coiled-coil domain can modulate Beclin-1 activity and suggest that PIST interactions with other proteins through its PDZ domain may regulate the activity of PIST and Beclin-1.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal PIST antibody was raised against a 16 amino acid peptide from near the carboxy terminus of human PIST (Genbank accession No. AAG00572).
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.h application have to be determined individually. PIST antibody can be used for the detection of PIST by Western blot at 1 - 4 µg/mL. Antibody can also be used for immunohistochemistry starting at 1 µg/mL. For immunofluorescence start at 10 µg/mL.
Images	Available upon request.
References	Gozuacik D and Kimchi A. Autophagy as a cell death and tumor suppressor mechanism. <i>Oncogene</i> . 2004; 23:2891-906. Kisen GO, Tessitore L, Costelli P, et al. Reduced autophagic activity in primary rat hepatocellular carcinoma and ascites hepatoma cells. <i>Carcinogenesis</i> 1993; 14:2501-5. Kamada Y, Funakoshi T, Shintani T, et al. Tor-mediated induction of autophagy via Apg1 protein kinase complex. <i>J. Cell. Biol.</i> 2000; 150:1507-13. Yu Z, Horton A, Bravin M, et al. A novel protein complex linking the ?2 glutamate receptor and autophagy: implications for neurodegeneration in Lurcher mice. <i>Neuron</i> 2002; 35:921-33.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-3631P Rat colon Cell Lysate, Cat. No. PK-AB718-1472 PIST Antibody (IN), Cat. No. PK-AB718-3633 TOR Antibody, Cat. No. PK-AB718-3485 Beclin-1 Antibody (CT), Cat. No. PK-AB718-3611

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