

Instruction Manual

Catalog Number	PK-AB718-4659
Synonyms	CLIP170 Antibody: CAP-Gly domain-containing linker protein 1, CLIP, CYLN1, Restin, RSN
Description	CLIP170 was initially identified as a new type of intermediate filament associated protein that is highly expressed in Reed-Sternberg cells, the tumoral cells diagnostic for Hodgkin's disease. Later experiments showed that it is located at microtubule plus ends and is required for the binding of endocytic carrier vesicles. CLIP170 has also been suggested to act with LIS1, a protein implicated in brain development, to regulate dynein/dynactin binding microtubules. Other studies suggest that CLIP170 can influence the formation of lamellipodia and cell invasion by invasive breast cancer cells by regulating the release of kinesin and IQGAP1 from a complex of those proteins, CLIP170 and Rac1. At least two isoforms of CLIP170 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	CLIP170 antibody was raised in rabbits against a 17 amino acid peptide from near the carboxy terminus of human CLIP170.
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, IF Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. CLIP170 antibody can be used for detection of CLIP170 by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunofluorescence starting at 20 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Bilbe G, Delabie J, Bruggen J, et al. Restin: a novel intermediate filament-associated protein highly expressed in the Reed-Sternberg cells of Hodgkin's disease. <i>EMBO J.</i> 1992; 11:2103-13. Diamantopoulos GS, Perez F, Goodson HV, et al. Dynamic localization of CLIP-170 to microtubule plus ends is coupled to microtubule assembly. <i>J. Cell Biol.</i> 1999; 144:99-112. Pierre P, Scheel J, Richard JE, et al. CLIP-170 links endocytic vesicles to microtubules. <i>Cell</i> 70:887-900. Suzuki K and Takahashi K. Regulation of lamellipodia formation and cell invasion by CLIP-170 in invasive human breast cancer cells. <i>Biochem. Biophys. Res. Commun.</i> 2008; 368:199-204.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4659P; CLIP170 Peptide Cat. No. PK-AB718-1463; Rat Brain Tissue Lysate

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