

BICD1 (CT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat BICD1 (Bicaudal D homolog 1)

Instruction Manual

Catalog Number	PK-AB718-4663
Synonyms	BICD1 Antibody: Bicaudal D homolog 1
Description	BICD1 was initially identified as the human homolog to Drosophila Bicaudal-D protein that forms part of the cytoskeleton and mediates the correct sorting of mRNAs for oocyte- and axis-determining factors during oogenesis. BICD1 can bind to dynein-dynactin and co-localizes with Rab6A on the trans-Golgi network and cytoplasmic vesicles, suggesting BICD1 regulates COPI-independent Golgi-ER transport. BICD1 will also interact with Rab6B, a Rab6 isoform that is expressed primarily in brain. Live cell imaging showed bi-directional movement of Rab6B structures in SK-N-SH neurites, indicating that BICD1 is involved in the Rab6B regulation of retrograde cargo transport in neuronal cells. At least two isoforms of BICD1 are known to exist; this antibody will only recognize the larger isoform.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	BICD1 antibody was raised in rabbits against a 15 amino acid peptide from near the carboxy terminus of human BICD1.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	At least two isoforms of BICD1 are known to exist; this antibody will only recognize the larger isoform.
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, ICC, IF I Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. BICD1 antibody can be used for detection of BICD1 by Western blot at 1 µg/mL. Antibody can also be used for immunocytochemistry starting at 2.5 µg/mL. For immunofluorescence start at 10 µg/mL.
Images	Available upon request.
References	Baens M and Marynen P. A human homologue (BICD1) of the drosophila bicaudal-D gene. <i>Genomics</i> 1997; 45:601-6. Matanis T, Akhmanova A, Wulf P, et al. Bicaudal-D regulates COPI-independent Golgi-ER transport by recruiting the dynein-dynactin motor complex. <i>Nat. Cell Biol.</i> 2002; 986-92. Wanshers BF, van de Vorstenbosch R, Schlager MA, et al. A role for the Rab6B Bicaudal-D1 interaction in retrograde transport in neuronal cells. <i>Exp. Cell Res.</i> 2007; 313:3408-20.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4663P; BICD1 Peptide Cat. No. PK-AB718-1224; Daudi Cell Lysate

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