

Instruction Manual

Catalog Number	PK-AB718-4663P
Quantity	50 µg
Source	17 amino acids near the carboxy terminus of human BICD1
Formulation	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH ₂ PO ₄ , 10 mM Na ₂ HPO ₄ , 130 mM NaCl) containing 0.1% bovine serum albumin and 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	Store BICD1 peptide at -20°C, stable for one year.
Application	BICD1 peptide is used for blocking the activity of BICD1 antibody.
References	Baens M and Marynen P. A human homologue (BICD1) of the drosophila bicaudal-D gene. Genomics 1997; 45:601-6.

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