

FEZ2 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat FEZ2 (IN)

Instruction Manual

Catalog Number	PK-AB718-4405
Synonyms	FEZ2 Antibody: Fasciculation and elongation protein zeta-2, zygin 2
Description	Fasciculation and elongation protein zeta-2 (FEZ2) is a homolog to the mammalian FEZ1, itself an ortholog of the <i>C. elegans</i> UNC-76. In contrast to FEZ1, FEZ2 mRNA is widely expressed in mouse tissues. FEZ2 interacts with protein kinase C (PKC)-z substrate and induces neurite extension of PC12 cells when co-expressed with a constitutively active form of PKC-z, suggesting FEZ2 may play an important role in the morphological changes of various cells by associating with PKC-z in a tissue non-specific manner. FEZ2 can interact with FEZ1 through its c-terminal regions and especially its coiled-coil region. At least two isoforms of FEZ2 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	FEZ2 antibody was raised in rabbits against a 15 amino acid peptide from near the center of human FEZ2.
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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. FEZ2 antibody can be used for the detection of FEZ2 by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunofluorescence starting at 5 µg/mL. For immunofluorescence start at 5 µg/mL.
Images	Available upon request.
References	Fujita T, Ikuta J, Okajima T, et al. Identification of a tissue-non-specific homologue of axonal fasciculation and elongation protein zeta-1. <i>Biochem. Biophys. Res. Commun.</i> 2004; 313:738-44. Kuroda S, Nakagawa N, Tokunaga C, et al. Mammalian homologue of the <i>Caenorhabditis elegans</i> UNC-76 protein involved in axonal outgrowth is a protein kinase C z-interacting protein. <i>J. Cell Biol.</i> 1999; 144:403-11. Assmann EM, Alborghetti MR, Camargo MER, et al. FEZ1 dimerization and interaction with transcription regulatory proteins involves its coiled-coil region. <i>J. Biol. Chem.</i> 2006; 281:9869-81.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4405P; FEZ2 Peptide Cat.No. PK-AB718-1212; 3T3 Cell Lysate

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