

ACVR2B Protein

Recombinant human protein expressed in Sf9 cells

Catalog # A36-35G

Lot # 1254-4

Product Description

Recombinant human ACVR2B (182-488) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_001106](#).

Gene Aliases

ActR-IIB; ACTRIIB; MGC116908

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

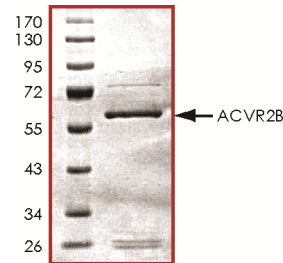
Scientific Background

ACVR2B or activin A receptor, type IIB is belong to the transforming growth factor-beta (TGF-beta) superfamily of structurally related signaling proteins that signal through a heteromeric complex of receptor serine kinases which include at least two type I (I and IB) and two type II (II and IIB) receptors (1). ACVR2B is help to regulate muscle growth (2). ACVR2B is a transmembrane protein that composed of a ligand-binding extracellular domain with cysteine-rich region, a transmembrane domain, and a cytoplasmic domain with predicted serine/threonine specificity.

References

- Bondestam, J. et.al: Assignment of ACVR2 and ACVR2B the human activin receptor type II and IIB genes to chromosome bands 2q22.2-q23.3 and 3p22 and the human follistatin gene (FST) to chromosome 5q11.2 by FISH. Cytogenet. Cell Genet. 87: 219-220, 1999.
- Lee, S.-J. et.al: Regulation of muscle growth by multiple ligands signaling through activin type II receptors. Proc. Nat. Acad. Sci. 102: 18117-18122, 2005.

Purity



The purity was determined to be **>85%** by densitometry. Approx. MW **61 kDa**.

ACVR2B Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number **A36-35G**

Specific Lot Number **1254-4**

Purity	>85%
Concentration	0.1 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.