

ACVR2A Protein

Recombinant human protein expressed in Sf9 cells

Catalog # A35-35H

Lot # 1025-1

Product Description

Recombinant human ACVR2A (161-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [NM_001616](#).

Gene Aliases

ACTRII; ACVR2

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 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.

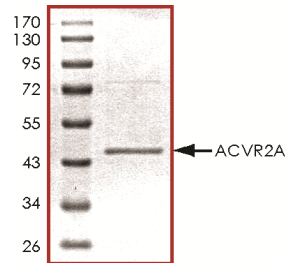
Scientific Background

ACVR2A encodes activin A type II receptor that are dimeric growth and differentiation factors which belong to the transforming growth factor-beta (TGF-beta) superfamily of structurally related signaling proteins. Activins is an autocrine/paracrine regulator in the human placenta (1). ACVR2A is highly expressed in human brain and ovary (2). Activins 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. These receptors are all transmembrane proteins, composed of a ligand-binding extracellular domain with cysteine-rich region, a transmembrane domain, and a cytoplasmic domain with predicted serine/threonine specificity.

References

1. Donaldson, C. J. et.al: Molecular cloning and binding properties of the human type II activin receptor. *Biochem. Biophys. Res. Commun.* 184: 310-316, 1992.
2. Peng, C. et.al: Expression of the type II activin receptor gene in the human placenta. *Endocrinology* 133: 3046-3049, 1993.

Purity



The purity was determined to be **>90%** by densitometry. Approx. MW **47 kDa**.

ACVR2A Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number **A35-35H**

Specific Lot Number **1025-1**

Purity	>90%
Concentration	0.1µg/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.