

Instruction Manual

Catalog Number	C-60068
Synonyms	ACVR2A, ACTRIIA, ACTR-IIA
Description	<p>ACVR2A takes part in various biological processes including mesoderm induction, neural cell differentiation, bone remodeling, hematopoiesis, carcinogenesis, and inflammation. ACVR2A which is a receptor for Actv A, Actv B and inhibin A mediates induction of adipogenesis by GDF-6.</p> <p>Human Recombinant ACVR2A produced in Sf9 cells is a single, glycosylated polypeptide chain of 124 amino acids (20-135a.a.), having a molecular mass of 14.4 kDa (molecular size on SDS-PAGE will appear at approximately 40-57 kDa). ACVR2A is fused to 8 amino acid His-tag at C-terminus and has been purified using proprietary chromatographic techniques.</p>
Quantity	10 µg
Molecular Mass	14.4 kDa
Source	Sf9 insect cells
Biological-Activity	NA
Specific Activity	NA
Formulation	Sterile-filtered, clear protein solution (0.5 mg/ml) containing phosphate-buffered saline (pH 7.4) and 10% glycerol.
Reconstitution	Please Note: Always centrifuge product briefly before opening vial. The dissolved protein can be diluted into other aqueous buffers and stored at -20°C for future use.
Purity	Greater than 95.0% as determined by SDS-PAGE.
Endotoxin Level	< 0.1 ng per µg protein (1EU/µg)
Storage	<p>Store at 4°C if entire vial will be used within 2-4 weeks.</p> <p>Store frozen at -20°C for longer periods of time. For long-term storage it is recommended to add a carrier protein (0.1% endotoxin-free HSA or BSA; e.g. Cat.No. C-69500A).</p> <p>Avoid multiple freeze-thaw cycles.</p>

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