

ADCK5 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # A13-11G

Lot # A1158-1

Product Description

Recombinant human ADCK5 (68-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [BC101659](#).

Gene Aliases

FLJ35454

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

ADCK5 belongs to the protein kinase superfamily, ADCK protein kinase family that contains a protein kinase domain (1). The function of this protein is not yet clear. It is not known if it has protein kinase activity and what type of substrates it would phosphorylate. An important paralog of this gene is ADCK1.

References

1. Manning G. et al: The protein kinase complement of the human genome. Science. 2002 Dec 6;298(5600): 1912-34.

Purity



The purity of ADCK5 protein was determined to be **>75%** by densitometry. Approx. MW **83 kDa**.

ADCK5 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	A13-11G
Lot #	A1158-1
Purity	>75%
Concentration	0.05 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.