

VAV1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # V14-31G

Lot # O915-2

Product Description

Recombinant human VAV1 (73-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_005428](#).

Gene Aliases

VAV

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.

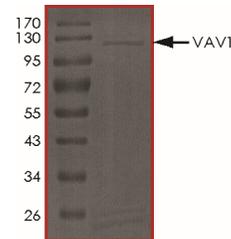
Scientific Background

VAV1 is a member of the VAV gene family which are guanine nucleotide exchange factors (GEFs) for Rho family GTPases. VAV1 activates pathways leading to actin cytoskeletal rearrangements and transcriptional alterations. VAV1 is important in hematopoiesis, playing a role in T-cell and B-cell development and activation (1). VAV1 has been identified as the specific binding partner of Nef proteins from HIV-1. Co-expression and binding of Nef with VAV1 initiates profound morphological changes, cytoskeletal rearrangements and the JNK/SAPK signaling cascade, leading to increased levels of viral transcription and replication. VAV1 proto-oncogene participates in the signaling processes that mediate the antigen-induced activation of B lymphocytes (2).

References

- Zhang, R. et.al: . Defective signalling through the T- and B-cell antigen receptors in lymphoid cells lacking the vav proto-oncogene. Nature 374: 470-473, 1995.
- Bustelo, X. R.et.al: Tyrosine phosphorylation of the VAV proto-oncogene product in activated B cells. Science 256: 1196-1199, 1992.

Purity



The purity of VAV1 was determined to be **>70%** by densitometry. Approx. MW **120kDa**.

VAV1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	V14-31G
Lot #	O915-2
Purity	>70%
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.