

Anti-VAV

Rabbit Polyclonal Antibody

Catalog # V14-63R

Lot # R157-4

Cited Applications

- Western blot (1:1000)
- Immunoprecipitation (2µg)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the VAV protein

Cross Reactivity

- Western blot VAV from human cells
- VAV from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 172-176 of human VAV

Formulation

PBS, pH 7.4, 0.02% NaN₃, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

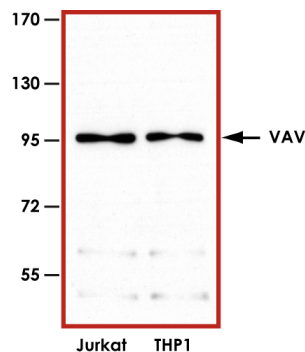
VAV is a proto-oncogene that is a member of the Dbl family of guanine nucleotide exchange factors (GEF) for the Rho family of GTP binding proteins. VAV plays a vital role in hematopoiesis and in T-cell and B-cell development as well as in their activation. VAV is the specific binding partner of Nef proteins from HIV-1 (1). Co-expression and binding of these VAV partners initiates profound morphological changes, cytoskeletal rearrangements and JNK/SAPK pathway activation leading to increased levels of viral transcription and replication. VAV is also the guanine nucleotide dissociation inhibitor for Ras.

References

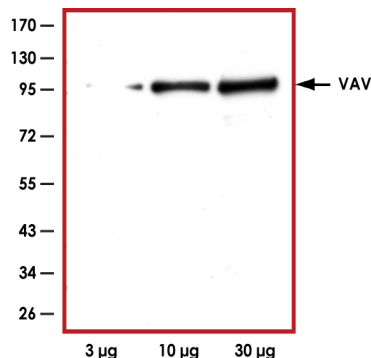
1. Fackler, O T. et al: Activation of Vav by Nef induces cytoskeletal rearrangements and downstream effector functions. *Molec. Cell* 3729-739, 1999.

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

Sample Data



Representative Immunoprecipitation with 2µg of Anti-VAV using 500µg of Jurkat and THP1 cell lysates.



Representative Western blot with Anti-VAV (1:1000) using 3µg, 10µg and 30µg of Jurkat cell lysate.

Anti-VAV

Rabbit Polyclonal Antibody

Catalog Number

V14-63R

Specific Lot Number

R157-4

Purification

Affinity chromatography

Concentration

1.0 µg/µL

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.