

Anti-VASP

Rabbit Polyclonal Antibody

Catalog # V12-63R

Lot # R133-6

Cited Applications

- Western blot (1:1000)

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

Specificity

Recognizes the VASP protein

Cross Reactivity

- Western blot human VASP

VASP from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 155-159 of human VASP

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

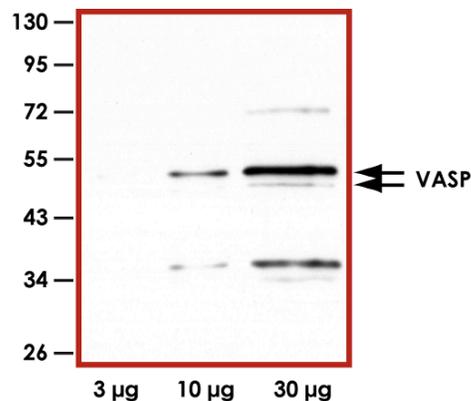
VASP or Vasodilator-stimulated phosphoprotein is a member of the Ena-VASP protein family which contains an EHV1 N-terminal domain that binds proteins containing E/DFPPPPXD/E motifs and targets Ena-VASP proteins to focal adhesions (1). VASP also contains proline-rich domain that bind SH3 and WW domain-containing proteins. VASP is associated with filamentous actin formation and likely plays a key role in cell adhesion and motility. VASP is also involved in the intracellular signaling pathways that regulate integrin-extracellular matrix interactions. VASP is regulated by the cyclic nucleotide-dependent kinases PKA and PKG.

References

1. Brindle, N P J. et al: The focal-adhesion vasodilator-stimulated phosphoprotein (VASP) binds to the proline-rich domain in vinculin. *Biochem. J.* 318:753-757, 1996.

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 Western blot with Anti-VASP (1:1000) using 3µg, 10µg and 30µg of THP1 cell lysate.

Anti-VASP

Rabbit Polyclonal Antibody

Catalog Number

V12-63R

Specific Lot Number

R133-6

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.