

VASP polyclonal antibody (affinity purified)

VASP (vasodilator stimulated phosphoprotein) is a proline-rich protein substrate of cAMP- and cGMP-dependent protein kinases. Phosphorylation of VASP at Ser-157 causes a mobility shift in SDS gel electrophoresis from 46 to 50 kDa, which has been used as a convenient marker to monitor cyclic nucleotide-dependent protein kinase activity. VASP is the founding member of the Ena-VASP protein family, comprising the *Drosophila* protein Enabled (Ena), its mouse homologue Mena (mammalian Enabled), and mouse EVL (Ena-VASP-like protein). With these proteins VASP shares a conserved overall domain organization:

- a) the conserved N-terminal Ena-VASP homology domain 1 (EVH1), which mediates binding to a proline-rich motif
- b) a more divergent proline-rich central domain (which is responsible for profilin binding)
- c) a conserved C-terminal EVH2 domain.

VASP is expressed in a variety of mammalian cell types and tissues. In cultured cells, VASP is associated with focal adhesions, cell-cell contacts, microfilaments, and highly dynamic membrane regions. From in vitro binding data VASP has been suggested to link profilin to zyxin, vinculin, and the *Listeria* spp. surface protein ActA, respectively. Functional evidence indicates that VASP is a crucial factor involved in the enhancement of actin filament formation and the actin-dependent motility of intracellular bacterial pathogens.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-8800

25µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Vasodilator stimulated phosphoprotein (affinity purified)
Application	ICC, IP, WB
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.02% sodium azide.
Host	Rabbit
Immunogen	Recombinant human VASP.
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (formaldehyde fixed cells, 1µg/ml, Immunoprecipitation (2µg/ml)Western Blot (0.1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Porcine
Specificity	Detects both the 46kDa (Ser ¹⁵⁷ dephospho) and 50kDa (Ser ¹⁵⁷ phospho) forms of VASP.
UniProt ID	P50552

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-](mailto:info-usa@enzolifesciences.com)
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-](mailto:info-eu@enzolifesciences.com)
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-](mailto:info-fr@enzolifesciences.com)
be@enzolifesciences.com

France
Phone: +33 472 440 655
[info-](mailto:info-fr@enzolifesciences.com)
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
[info-](mailto:info-de@enzolifesciences.com)
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-](mailto:info-uk@enzolifesciences.com)
uk@enzolifesciences.com

Last modified: May 29, 2024