

VASP (human)

polyclonal antibody

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: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-898-R100

100µl

Manuals, SDS & CofA

[View Online »](#)

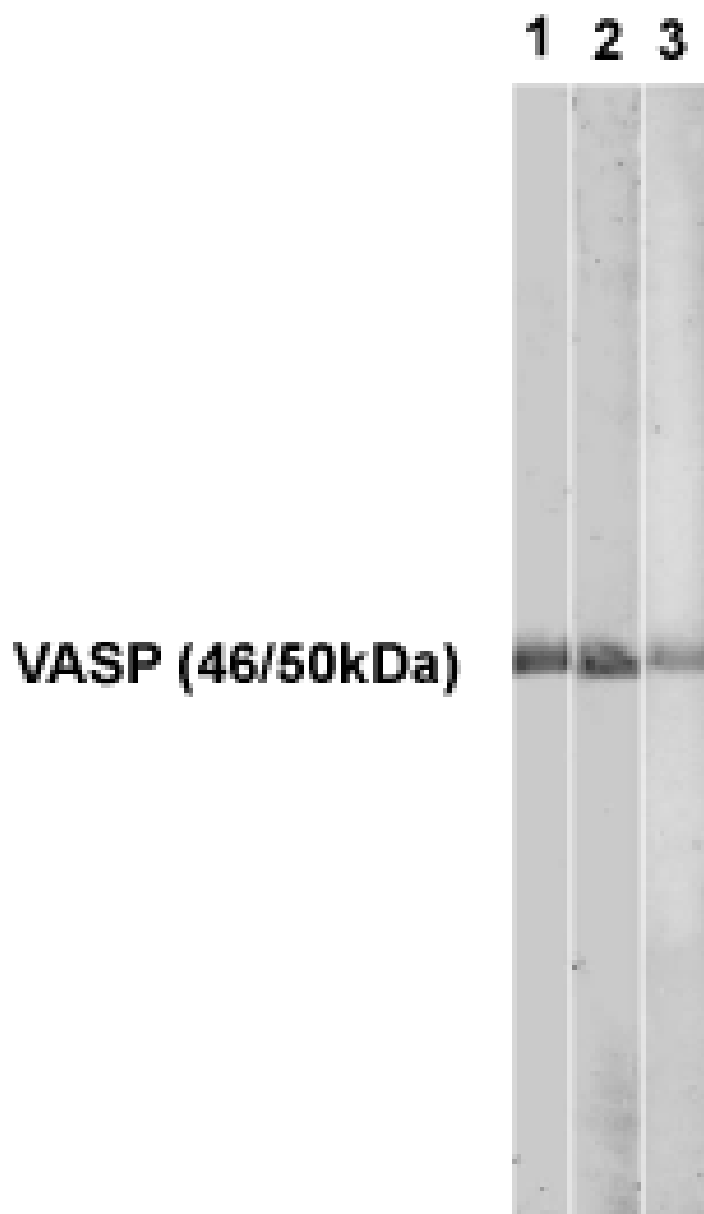


Figure: Western blot comparing three different PABs to VASP using 25µg human platelet protein. **Lane 1:** PAb to VASP (human) (Prod. N. ALX-210-898). **Lane 2:** PAb to VASP (M4) (Prod. N. ALX-210-725) (**Not available anymore!**). **Lane 3:** PAb to VASP (affinity purified) (Prod. N. ALX-210-880).

Handling & Storage

Handling	Avoid freeze/thaw cycles.
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
Application	IF, WB
Crossreactivity	Does not cross-react with mouse or rat VASP. For the detection of mouse VASP use Prod. No. ALX-210-880.
Formulation	Liquid. Contains 0.02% sodium azide.
Host	Rabbit
Immunogen	Recombinant human VASP (vasodilator stimulated phosphoprotein) fused to a His-tag.
Recommendation Dilutions/Conditions	Immunofluorescence of formaldehyde fixed cells (1:500-1:1000)Western Blot (1:1500-1:3000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Specificity	Recognizes unphosphorylated as well as phosphorylated (Ser ¹⁵⁷) 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-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

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

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

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