

[pSer239]VASP monoclonal antibody (16C2)

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

Citations: 12

[View Online »](#)

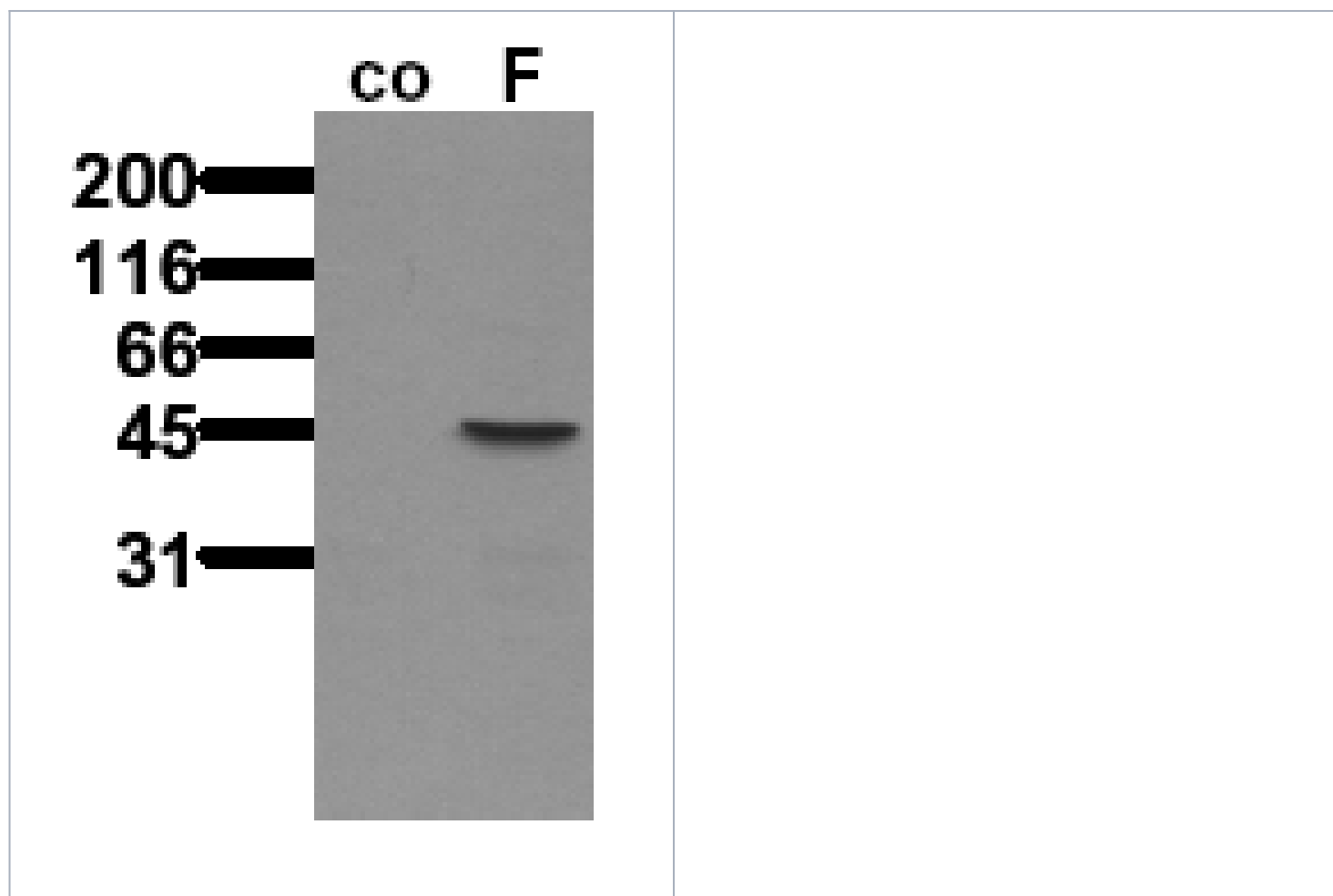
Ordering Information

[Order Online »](#)

ALX-804-240-C100	Purified - 100µg
------------------	------------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable at -80°C up to 1 year, at 4°C up to 3 months.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and freeze in liquid nitrogen.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Vasodilator stimulated phosphoprotein
Application	ELISA, Flow Cytometry, ICC, IP, WB
Clone	16C2
Crossreactivity	Does not cross-react with the non-phosphorylated form of VASP.
Formulation	Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.
Host	Mouse
Immunogen	Synthetic peptide corresponding to a portion of VASP (vasodilator stimulated phosphoprotein) phosphorylated at Ser239 (RKVpSKQE) conjugated to KLH.
Isotype	IgG1
Purity Detail	Thiophilic adsorption and size exclusion chromatography purified.

Recommendation Dilutions/Conditions

ELISA (use at 0.05µg/ml)Immunocytochemistry (use at 1-10µg/ml; may tolerate 0.5% formaldehyde fixation) , Immunoprecipitation (use 1-10 µg per 106 pervanadate-treated A431 cells)Western Blot (1µg/ml for AP/BCIP/NBT (MTT) detection; 0.5µg/ml for HRPO/ECL detection; recommended blocking buffer BPPT: 2x PBS 3% (w/v) BSA, 1% (w/v) PEG 4,000, 1% (w/v) polyvinylpyrrolidone, 0.1% (v/v) Tween 20)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Reconstitution

Reconstitute with 1ml water (15 minutes at room temperature).

Species Reactivity

Human, Mouse, Rabbit, Rat

Specificity

Recognizes Ser²³⁹-phosphorylated VASP.

Technical Info / Product Notes

Important! The use of MAb to VASP (activated) (Ser²³⁹) (16C2) for diagnostic or therapeutic purposes is patented! The antibody is supplied for research use only!

UniProt ID

P50552

Worry-free Guarantee

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

Last modified: May 29, 2024



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)