

[pSer239]VASP monoclonal antibody (16C2) (biotin conjugate)

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-240B-C100	Biotin - 100µg
-------------------	----------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Reconstituted antibody is stable for 1 year when stored at -80°C. Thaw aliquots at +37°C. Thawed aliquots may be stored at +4°C for 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	Liquid. 0.5mg/ml in 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.

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)