

# [pSer<sup>239</sup>]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

<b>Use/Stability</b>	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.
<b>Handling</b>	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and freeze in liquid nitrogen.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Vasodilator stimulated phosphoprotein
<b>Application</b>	ELISA, Flow Cytometry, ICC, IP, WB
<b>Clone</b>	16C2
<b>Crossreactivity</b>	Does not cross-react with the non-phosphorylated form of VASP.
<b>Formulation</b>	Liquid. 0.5mg/ml in 2x PBS containing 0.09% sodium azide, PEG and sucrose.
<b>Host</b>	Mouse
<b>Immunogen</b>	Synthetic peptide corresponding to a portion of VASP (vasodilator stimulated phosphoprotein) phosphorylated at Ser239 (RKVpSKQE) conjugated to KLH.
<b>Isotype</b>	IgG1
<b>Purity Detail</b>	Thiophilic adsorption and size exclusion chromatography purified.
<b>Recommendation Dilutions/Conditions</b>	ELISA (use at 0.05µg/ml)Immunocytochemistry (use at 1-10µg/ml; may tolerate 0.5% formaldehyde fixation)Immunoprecipitation (use 1-10 µg per 10 <sup>6</sup> 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.
<b>Species Reactivity</b>	Human, Mouse, Rabbit, Rat

<b>Specificity</b>	Recognizes Ser <sup>239</sup> -phosphorylated VASP.
<b>Technical Info / Product Notes</b>	<b>Important! The use of MAb to VASP (activated) (Ser<sup>239</sup>) (16C2) for diagnostic or therapeutic purposes is patented! The antibody is supplied for research use only!</b>
<b>UniProt ID</b>	P50552
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)