

# [pSer239]VASP monoclonal antibody (16C2) (FITC conjugate)

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

Citations: 9

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-240F-C100	FITC - 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	Flow Cytometry, ICC
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	Immunocytochemistry (use at 1-10µg/ml; may tolerate 0.5% formaldehyde fixation)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 Ser<sup>239</sup>-phosphorylated VASP.

**Technical Info / Product Notes**

**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!**

**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)