

[pSer¹⁵⁷]VASP monoclonal antibody (5C6)

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

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-403-C100	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, FUNC, ICC, IP, WB
Application Notes	Functional Application: monitoring of PKA activity.
Clone	5C6
Crossreactivity	Does not cross-react with the non-phosphorylated form of VASP.
Formulation	Lyophilized from 1ml of PBS containing 0.09% sodium azide, PEG, and sucrose.
Host	Mouse
Immunogen	Synthetic peptide corresponding to a portion of human VASP (vasodilator-stimulated phosphoprotein) phosphorylated at Ser157 conjugated to KLH.
Isotype	IgG1
Purity Detail	Thiophilic adsorption and size exclusion chromatography purified.

Recommendation Dilutions/Conditions

Immunocytochemistry (1-10µg/ml; may tolerate 0.5% formaldehyde fixation)Western Blot> (1µg/ml for AP/BCIP/NBT (MTT) detection; 0.5µg/ml for HRPO/ECL detection)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

Technical Info / Product Notes

The use of MAb to VASP (activated) (Ser¹⁵⁷) (5C6) for diagnostic or therapeutic purposes is patented! The antibody is supplied for research use only!

Replacement for **ADI-905-768**

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)