[pSer²³⁹]VASP monoclonal antibody (16C2)

This antibody is covered by our Worry-Free Guarantee.

Citations: 12

View Online »

Ordering Information

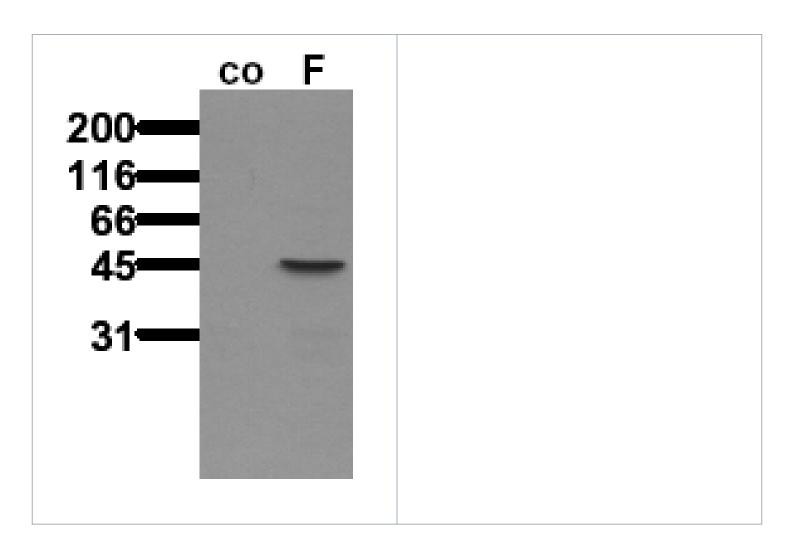
Order Online »

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.

lgG1

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

