[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.

lsotype 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.

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