Phosphoserine monoclonal antibody (16B4)

This antibody is covered by our Worry-Free Guarantee.

Citations: 14

View Online »

Ordering Information

Order Online »

ALX-804-167-C100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable at -80°C up to 1 year, at 4°C up to 1 week.

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

Application ELISA, IP, WB

Clone 16B4

Formulation Lyophilized from 1ml of 2x PBS containing 0.09% sodium

azide, PEG, and sucrose.

Host Mouse

Immunogen Phosphoserine containing peptides conjugated to KLH.

lsotype lgM

Positive Control Included. (Prod. No. 840-012-1)

Purity Detail Thiophilic adsorption and size exclusion chromatography

purified.

Recommendation Dilutions/Conditions ELISA (use at 0.05µg/ml)Western Blot (1µg/ml 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. Do not use

milk/casein for blocking and dilution! Replace by

BSA!)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

Species independent

Specificity

Recognizes a broad range of serine-phosphorylated proteins in crude cell extracts, strongly interacting with pSK and pSP motifs, thus specifically recognizing substrates of MAP/SAP kinases and CDC kinases. Please note that phosphoserine detection by MAbs always depends on the surrounding aa sequence.

Technical Info / Product Notes

Replacement for ADI-905-805

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-