

# Phosphoserine monoclonal antibody (16B4)

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

Citations: 14

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable at -80°C up to 1 year, at 4°C up to 1 week.
<b>Handling</b>	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and freeze in liquid nitrogen.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Application</b>	ELISA, IP, WB
<b>Clone</b>	16B4
<b>Formulation</b>	Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.
<b>Host</b>	Mouse
<b>Immunogen</b>	Phosphoserine containing peptides conjugated to KLH.
<b>Isotype</b>	IgM
<b>Positive Control</b>	Included. (Prod. No. 840-012-1)
<b>Purity Detail</b>	Thiophilic adsorption and size exclusion chromatography purified.
<b>Recommendation Dilutions/Conditions</b>	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.
<b>Reconstitution</b>	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**

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)