

# Phosphoserine polyclonal antibody

Protein phosphorylation is a complex biochemical process that is regulated by protein kinases that catalyze the transfer of the gamma-P of ATP to a recipient protein that acts as a substrate. This post-translational modification is primarily directed onto serine, threonine, and tyrosine amino acid residues within a protein. Changes in the phosphorylation status of phosphoproteins can alter their biological function, and aberrations in signal transduction cascades have been linked to a number of diseases, including cancer, diabetes, heart disease, inflammation and neurological disorders.

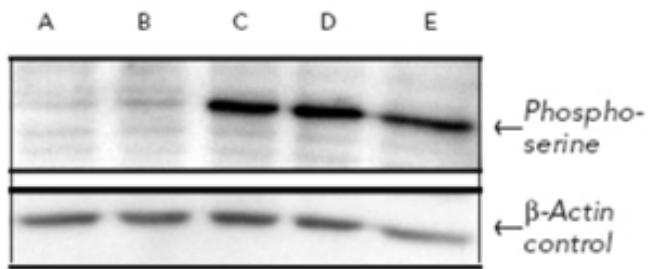
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of human MMRU melanoma cells probed with Phosphoserine, pAb. Top (from the left): MMRU cells treated with 0 (A), 50 (B), 200 (C), 400 (D) long wavelength UV, or 0.1 $\mu$ M of okadaic acid (E). Bottom: Same samples probed with  $\beta$ -actin antibody (control).

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application ELISA, IHC, IP, WB

Formulation Liquid. In PBS with 50% glycerol.

Host Rabbit

Immunogen Phosphoserine conjugated to KLH.

Purity Detail Peptide affinity purified.

**Recommendation Dilutions/Conditions** ELISA (1µg/ml)Immunohistochemistry (5µg/ml)Immunoprecipitation (10µg/mg protein)Western Blot (1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Species independent

**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)