

Phosphothreonine 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 »](#)

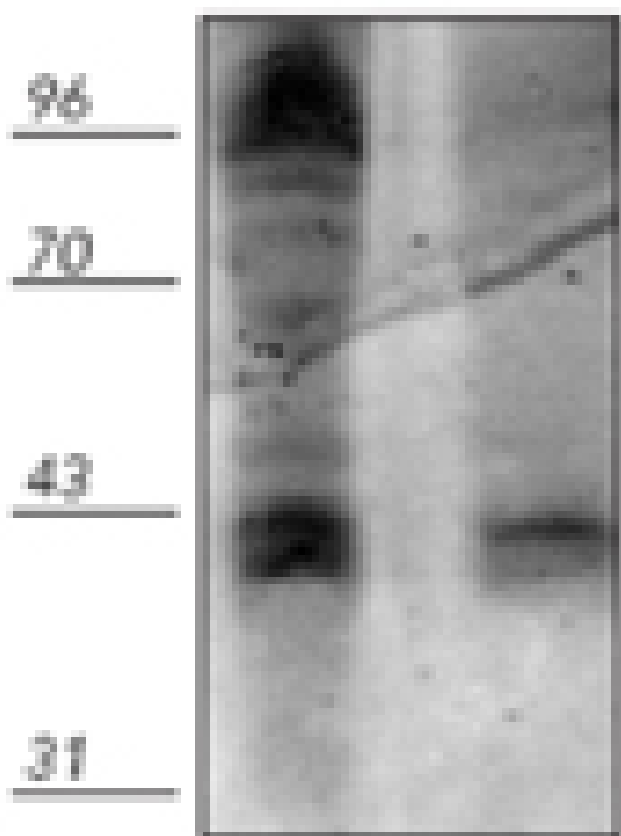
Ordering Information

[Order Online »](#)

ADI-KAP-ST211-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of mouse spleen extract probed with Prod. No. ADI-KAP-ST211. Left Lane: Vanadium treated; Right lane: Untreated.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, IP, WB

Formulation Liquid. In PBS, pH 7.0 containing 50% glycerol.

Host Rabbit

Immunogen Phosphothreonine conjugated to KLH.

Purity Detail Immunoaffinity purified.

Recommendation Dilutions/Conditions ELISA (0.5µg/ml)Immunoprecipitation (10µg/mg protein)Western Blot (2µg/ml, colorimetric)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)