

[pSer¹⁹]αB-Crystallin polyclonal antibody

Alpha-crystallins, which are part of the small Heat shock family members, are major water-soluble proteins present in the lens of the mammalian eye. Phosphorylation of serine residues which occurs during development and in response to stress, is intimately linked with its function. Chaperone activity requires, and is modulated by, oligomerization and is limited to binding unfolded intermediates to prevent irreversible aggregation.

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

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

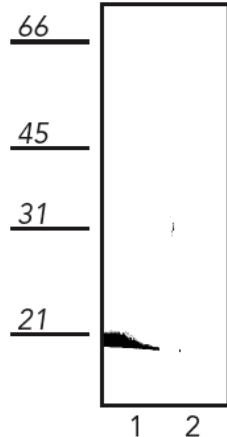
ADI-SPA-225-C	25µg
ADI-SPA-225-E	100µg

Manuals, SDS & CofA

[View Online »](#)

phosphorylated
 α B-Crystallin

non-phosphorylated
 α B-Crystallin



Western blot analysis
of phosphorylated
 α B-Crystallin and
non-phosphorylated
 α B-Crystallin proteins,
probed with SPA-225.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CRYAB, HSPB5
Application	IP, WB
Formulation	Liquid. In sodium phosphate buffer, pH 7.0, containing 3% BSA.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human α B-crystallin. The sequence is completely conserved in mouse, rat and bovine phosphorylated at Ser19.
Recommendation Dilutions/Conditions	Immunoprecipitation (12.5 μ g/ml)Western Blot (1 μ g/ml, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Bovine, Human, Mouse
UniProt ID	P02511
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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@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)