

# [pSer19] $\alpha$ B-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 »](#)

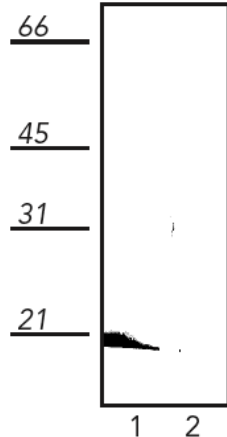
ADI-SPA-225-C	25 $\mu$ g
ADI-SPA-225-E	100 $\mu$ g

## Manuals, SDS & CofA

[View Online »](#)

phosphorylated  
 $\alpha$ B-Crystallin

non-phosphorylated  
 $\alpha$ B-Crystallin



Western blot analysis  
of phosphorylated  
 $\alpha$ B-Crystallin and  
non-phosphorylated  
 $\alpha$ 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  $\alpha$ B-crystallin. The sequence is completely conserved in mouse, rat and bovine phosphorylated at Ser19.

Recommendation Dilutions/Conditions      Immunoprecipitation (12.5 $\mu$ g/ml)Western Blot (1 $\mu$ 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)