

# [pSer<sup>59</sup>]αB-Crystallin polyclonal antibody

α-Crystallins composed of αA (~19 kDa) and αB (~19.2 kDa) subunits, are major water-soluble proteins accounting for almost 50% of total protein in the mammalian transparent eye lens and they are also found in a variety of other tissues. Alpha-crystallins are also referred to as small heat shock proteins, since they are induced by increased temperature in a variety of organisms. The α-crystallins have sequence homology as well as structural and functional similarities with the small Hsp's such as Hsp25/27. Most α-crystallins have four common structural and functional features: (i) molecular weight between 12 and 43kDa; (ii) the formation of large oligomeric complexes composed of αA-crystallin, αB-crystallin and Hsp25/27; (iii) the moderately conserved α-crystallin domain in the central region of the protein; and (iv) molecular chaperone activity. The α-crystallin domain comprises approximately 90 residues, is bounded by variable N-terminal and C-terminal extensions and is involved in oligomer assembly. Oligomers can reach 800 kDa or more and are dynamic, exhibiting subunit exchanges and organizational plasticity, possibly leading to functional diversity. Phosphorylation of serine residues occurs during development and in response to stress, and usually decreases oligomer size. Chaperone activity requires, and is modulated by, oligomerization and is limited to binding unfolded intermediates to prevent irreversible aggregation. Although productive release and refolding of denatured proteins requires close cooperation with other chaperones. Other proposed functions include a role in membrane stabilization and modulation of intermediate filament organization during physiological stress and neurodegenerative disease. Alpha B Crystallin is phosphorylated on serine 45 by ERK1/2 and serine 59 by MAPKAPK-2. When mono-phosphorylated αB-crystallin is isolated it contains only αB-crystallin phosphorylated at serine 19.

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

Citations: 16

[View Online »](#)

Ordering Information

[Order Online »](#)

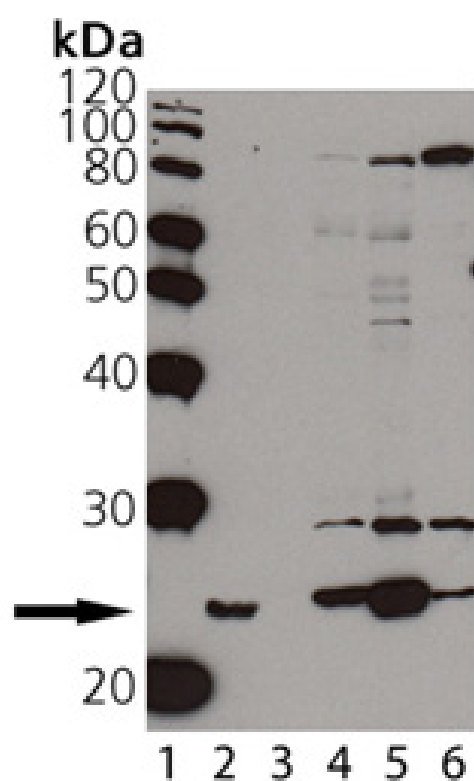
ADI-SPA-227-D

50µg

ADI-SPA-227-F	200µg
---------------	-------

**Manuals, SDS & CofA**

**[View Online »](#)**



Western blot analysis of [pSer59]  $\alpha$ B-Crystallin: Lane 1: MWM, Lane 2:  $\alpha$ -Crystallin (bovine), (native) (Prod No. ADI-SPP-225), Lane 3:  $\beta$ -Crystallin (bovine), (native) (Prod No. ADI-SPP-235), Lane 4: Mouse Heart Tissue Lysate, Lane 5: Rat Heart Tissue Lysate, Lane 6: ESK-4 Cell Lysate

## 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 0.1% sodium azide, 3% BSA, and 50% glycerol.

Host      Rabbit

Immunogen      Synthetic peptide corresponding to aa 54-64 of human  $\alpha$ B-crystallin phosphorylated at Ser59.

Recommendation      Western Blot (ECL, 1:1,000) 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, Mouse, Porcine, Rat, Sheep

UniProt ID      P02511

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



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)