

# $\alpha$ A-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: 14

[View Online »](#)

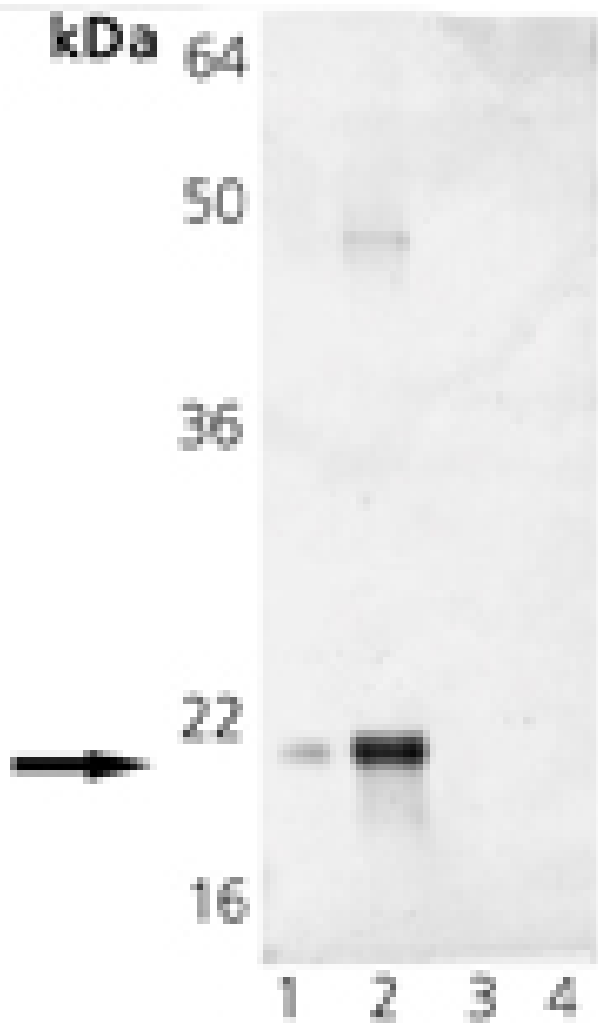
### Ordering Information

[Order Online »](#)

ADI-SPA-221-D	50µg
ADI-SPA-221-F	200µg

### Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Crystallin, alpha-A pAb: Lane 1: Crystallin, alpha Native Bovine Protein, Lane 2: Crystallin, alpha A Native Bovine Protein, Lane 3: Crystallin, alpha B Native Bovine Protein (Negative Control), Lane 4: Crystallin, beta Native Bovine Protein (Negative Control).

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name CRYAA, HSPB4

Application IF, IP, WB

Application Notes Detects a band of ~20kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

GenBank ID U05569

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the C-terminus of human  $\alpha$ A-Crystallin. The sequence is completely conserved in frog and chicken.

Purity Detail Protein A affinity purified.

**Recommendation** Western Blot (1:1,000, colorimetric) Suggested dilutions/conditions may not be available  
**Dilutions/Conditions** for all applications. Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Human, Mouse

UniProt ID P02489

**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-  
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)