

α A/ α B-Crystallin polyclonal antibody

Alpha-crystallins composed of alpha A (~20 kDa) and alpha B (~20 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. The two other crystallin families, beta and gamma, are homologous to each other but not to the alpha family or the sHsp's. Alpha-crystallins are also referred to as small heat shock proteins, since they are induced by increased temperature in a variety of organisms. The alpha-crystallins have sequence homology as well as structural and functional similarities with the small Hsp's such as Hsp25/27. Most small heat shock proteins have four common structural and functional features: (i) molecular weight between 12 and 43kDa; (ii) the formation of large oligomeric complexes composed of alphaA-crystallin, alphaB-crystallin and Hsp25/27; (iii) the moderately conserved alpha-crystallin domain in the central region of the protein; and (iv) molecular chaperone activity. The alpha-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 800kDa 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.

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

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

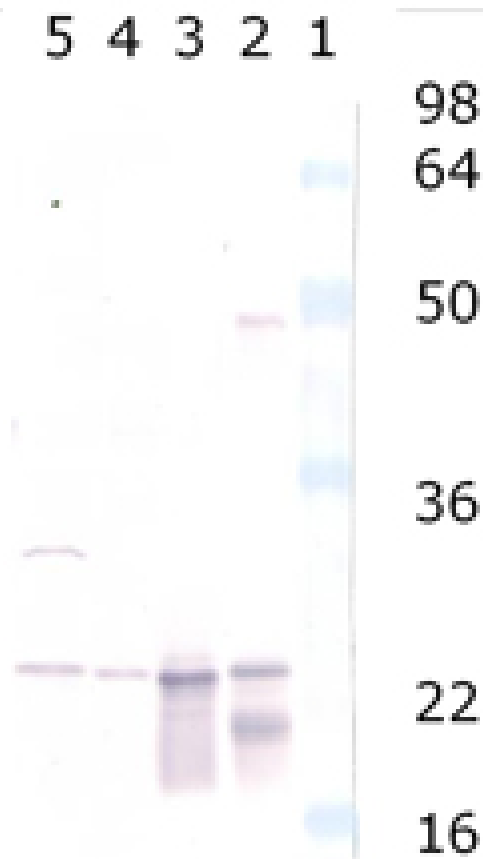
ADI-SPA-224-D

50 μ g

ADI-SPA-224-F	200µg
---------------	-------

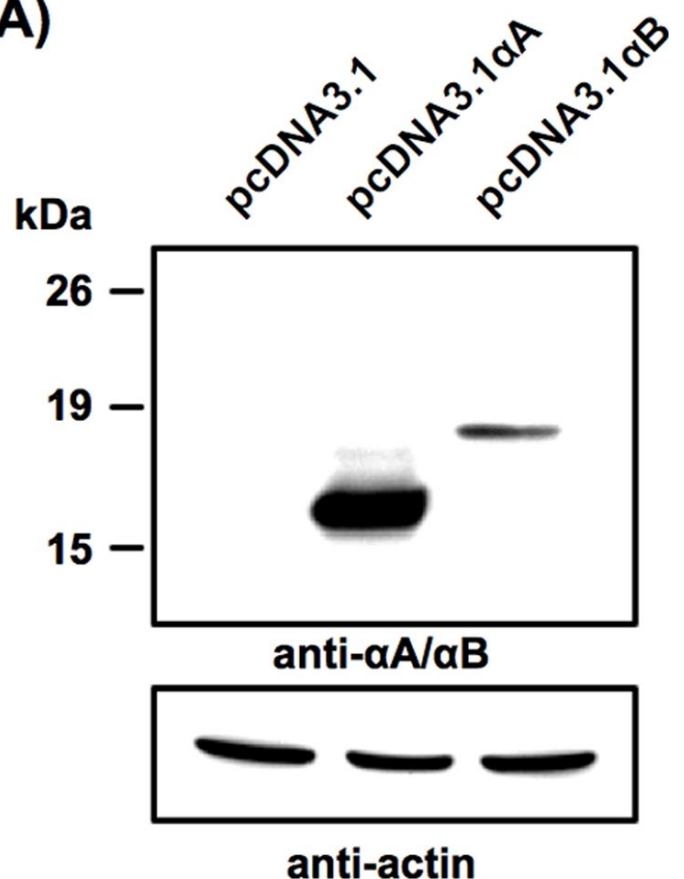
Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: Lane 1: MWM, Lane 2: Prod. No. ADI-SPP-225, Lane 3: Prod. No. ADI-SPP-227, Lane 4: HeLa HS, Lane 5: Mouse Heart.

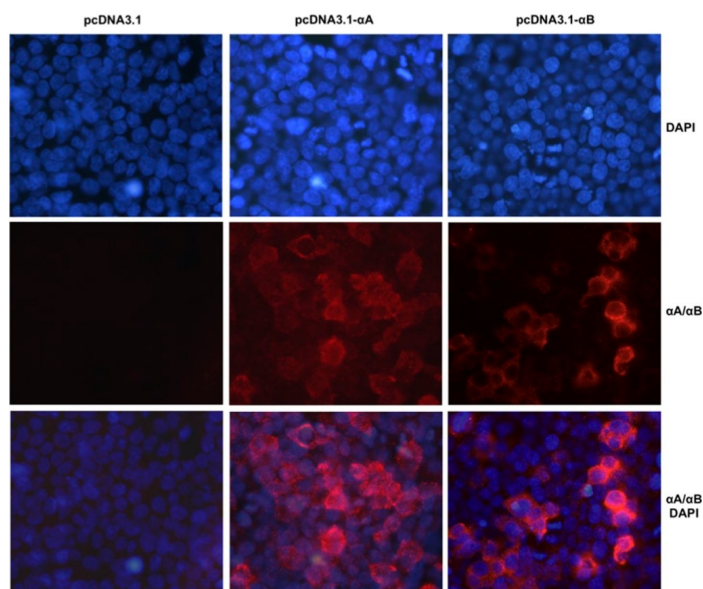
(A)



Expression of α A- and α B-crystallins in transiently transfected 293T cells. (A) Western blot analysis of α A- and α B-crystallin levels 24 h post-transfection. Fifty micrograms of total proteins from cell extracts were subjected to 12% SDS-PAGE and immunoassayed with anti- α A/ α B-crystallin to detect the overexpressed α -crystallins and with anti- β -actin as a control of equal protein loading. (B) Immunofluorescence analysis with anti- α A/ α B-crystallin showing cytoplasmic expression of α A- (pcDNA3.1- α A) and α B- (pcDNA3.1- α B) crystallins 24 h post-transfection, while no detection was observed in cells transfected with the empty plasmid (pcDNA3.1).

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Analysis of the cytoprotective role of α -crystallins in cell survival and implication of the α A-crystallin C-terminal extension domain in preventing Bax-induced apoptosis. *PLoS One* (2013)

(B)



Expression of α A- and α B-crystallins in transiently transfected 293T cells. (A) Western blot analysis of α A- and α B-crystallin levels 24 h post-transfection. Fifty micrograms of total proteins from cell extracts were subjected to 12% SDS-PAGE and immunoassayed with anti- α A/ α B-crystallin to detect the overexpressed α -crystallins and with anti- β -actin as a control of equal protein loading. (B) Immunofluorescence analysis with anti- α A/ α B-crystallin showing cytoplasmic expression of α A- (pcDNA3.1- α A) and α B- (pcDNA3.1- α B) crystallins 24 h post-transfection, while no detection was observed in cells transfected with the empty plasmid (pcDNA3.1).

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Analysis of the cytoprotective role of α -crystallins in cell survival and implication of the α A-crystallin C-terminal extension domain in preventing Bax-induced apoptosis. *PLoS One* (2013)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CRYAA/CRYAB, HSPB4/HSPB5

Application ICC, WB

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

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

GenBank ID U05569 (alphaA-crystallin), M28638 (alphaB-crystallin)

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the C-terminus of human α A- and α B-Crystallin.

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 (alphaA-crystallin), P02511 (alphaB-crystallin)

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)