

α B-Crystallin monoclonal antibody (1B6.1-3G4)

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: 1

[View Online »](#)

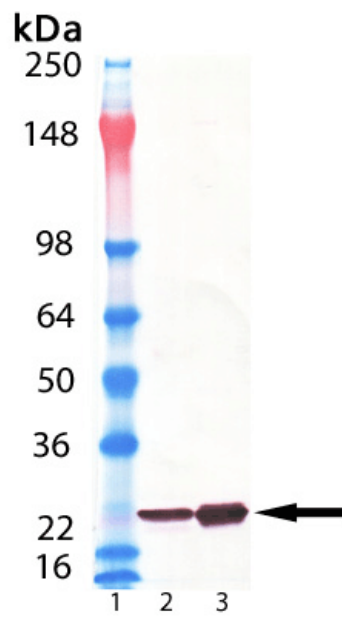
Ordering Information

[Order Online »](#)

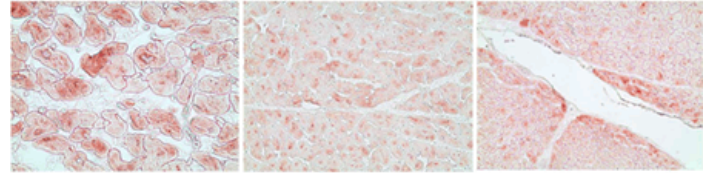
| | |
|-----------------|-------|
| ENZ-ABS706-0200 | 200µg |
|-----------------|-------|

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of α B-Crystallin mAb (1B6.1-3G4) (Prod. No. ENZ-ABS706): Lane 1: MW Marker, Lane 2: α -Crystallin (bovine), (native) (Prod. No. ADI-SPP-225), Lane 3: α B-Crystallin (bovine), (native) (Prod. No. ADI-SPP-227).



Immunohistochemistry analysis of human heart tissue with alpha B-Crystallin, mAb (1B6.1-3G4), (Prod. No. ENZ-ABS706), at 20 μ g/ml.

| | |
|--------------------------|---------------------------|
| Handling | Avoid freeze/thaw cycles. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

| | |
|-------------------------------------|--|
| Alternative Name | CRYAB, HSPB5 |
| Application | ICC, IF, IHC (PS), WB |
| Application Notes | Detects a band of ~20kDa by Western blot. |
| Clone | 1B6.1-3G4 |
| Formulation | Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide. |
| Host | Mouse |
| Immunogen | Bovine αB-crystallin. |
| Isotype | IgG1 |
| Purity Detail | Protein G affinity purified. |
| Recommendation Dilutions/Conditions | Immunohistochemistry (paraffin sections, 10µg/ml)Western Blot (1:1,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| Source | Purified from hybridoma tissue culture supernatant. |
| Species Reactivity | Bovine, Chicken, Human, Mouse, Rat |
| UniProt ID | P02510 |

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-
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)