

[pSer⁴⁵]αB-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: 5

[View Online »](#)

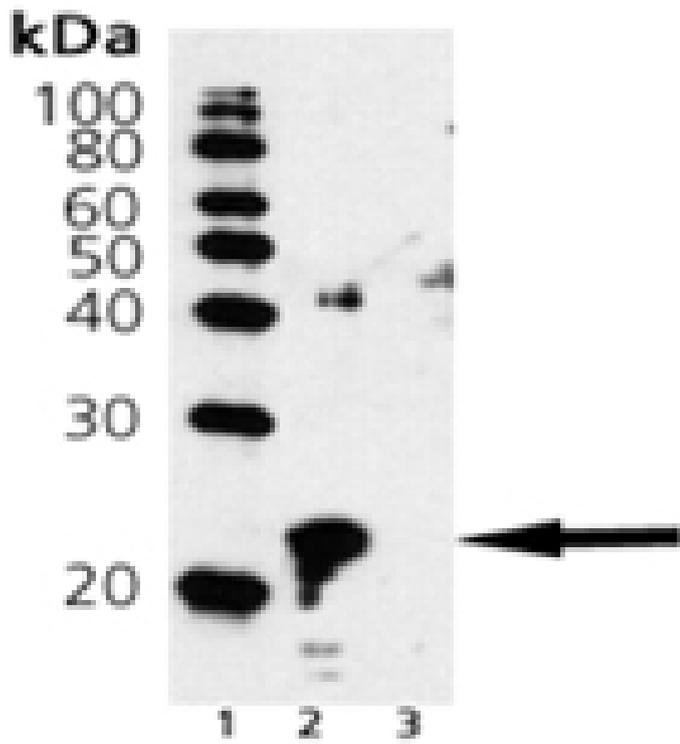
Ordering Information

[Order Online »](#)

ADI-SPA-226-D	50µg
ADI-SPA-226-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Crystallin alpha-B (pSer45):
Lane 1: MW marker, Lane 2: alpha-B1 Crystallin Protein (phospho), Lane 3: alpha-B2 Crystallin Protein (non-phospho)

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

Application Notes

Detects a band of ~22kDa by Western blot.

Formulation

Liquid. In 0.1M sodium phosphate buffer, pH 7.0, containing 1.5% BSA and 0.1% sodium azide.

Host

Rabbit

Immunogen

Synthetic peptide corresponding to a portion of human α B-crystallin phosphorylated at Ser45.

Purity Detail

Affinity purified.

Recommendation Dilutions/Conditions

Western Blot (1:500, ECL) 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, Chicken, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Scallop, Xenopus

UniProt ID

P02511

Worry-free Guarantee

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



ENZO LIFE SCIENCES,
INC.
Phone: +1 949 948 9120
info@enzolifesciences.com
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com
uk@enzolifesciences.com