

IMMUNOSET[®] αB-Crystallin ELISA development set

The IMMUNOSET[®] αB-Crystallin ELISA development set contains sufficient reagents

for the development of 5 x 96 well EIA plates which provides analysis of ~200 samples in duplicate.

The IMMUNOSET[®] αB-Crystallin EIA development set contains sufficient reagents for the development of 5 x 96 well EIA plates which provides analysis of ~200 samples in duplicate. Save time, money, and sample by developing your own EIA kits. Simply coat your plates and go! This kit provides superior accuracy with fully quantitative results compared to Western blot analysis and has low reactivity with related molecules.

α-crystallins composed of ~20kDa αA and αB subunits function as major water-soluble proteins accounting for almost 50% of total protein in the mammalian transparent eye lens, also existing in a variety of other tissues. The α-crystallin proteins are members of the small heat shock protein (sHsp) family, as their expression can be induced by heat and other stress insults in a variety of organisms. The α-crystallins possess structural and functional similarities and share sequence homology with Hsp25/273. The conserved α-crystallin domain participates in oligomer assembly, which is critical to their function in the prevention of irreversible protein aggregation.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

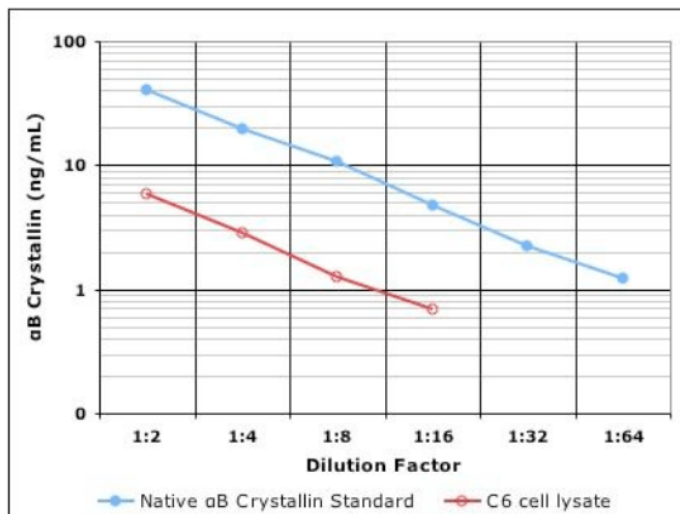
ADI-960-074

5x96 wells

Manuals, SDS & CofA

[View Online »](#)

- Higher throughput - assay ~200 samples in duplicate in a fraction of the time
- Accurate - fully quantitative results surpass semi-quantitative Western blot analysis
- Specific - low reactivity with related molecules



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CRYAB, HSPB5

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of α B-Crystallin in cell lysates of human, mouse, rat, and cow origin.

Contents Capture antibody, Standard, Detection antibody, SA-HRP

Crossreactivity Reacts with α B-Crystallin from human, mouse, rat, and bovine. Does not react with HSP10, HSP25, HSP27, HSP40, γ -crystallin, or β L-crystallin and reacts <0.4% with α A-crystallin.

Sensitivity 0.59 ng/ml (range 1.25 – 40 ng/ml)

UniProt ID P02511 (human), P02510 (bovine)

Wavelength 450 nm



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)