

Crystallin, α B (phospho-Ser45) Polyclonal Antibody

Product Specifications

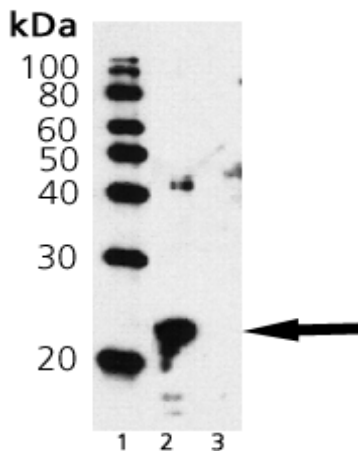
Catalog Number:	SPA-226										
Host:	Rabbit										
Species Reactivity:	Human, Mouse, Rat, Bovine, Canine, Chicken, Guinea Pig, Hamster, Monkey, Pig, Rabbit, Scallop, Xenopus Other species not tested.										
Application:	WB (ECL, 1:500), IP Other applications not tested. <i>The optimal dilution for a specific application must be determined by the investigator</i>										
Predicted MW:	~22 kDa										
Concentration:	See product label										
Purification:	Affinity Purified										
Format:	0.1M sodium phosphate buffer, pH 7.0, 0.1% azide, 1.5% BSA										
Storage:	-20°C <i>Shipping conditions may differ from the recommended storage temperature</i>										
Immunogen:	Synthetic phospho-peptide derived from the sequence of human α B Crystallin										
Related Products:	<table border="0"> <tr> <td>LYC-HL101</td> <td>HeLa Cell Lysate, Heat Shocked</td> </tr> <tr> <td>LYC-PC100</td> <td>PC-12 Cell Lysate</td> </tr> <tr> <td>LYC-3T100</td> <td>3T3 Cell Lysate</td> </tr> <tr> <td>SPA-227</td> <td>Crystallin, αB (phospho-Ser59) Polyclonal</td> </tr> <tr> <td>960-074</td> <td>Crystallin, αB ImmunoSet</td> </tr> </table>	LYC-HL101	HeLa Cell Lysate, Heat Shocked	LYC-PC100	PC-12 Cell Lysate	LYC-3T100	3T3 Cell Lysate	SPA-227	Crystallin, α B (phospho-Ser59) Polyclonal	960-074	Crystallin, α B ImmunoSet
LYC-HL101	HeLa Cell Lysate, Heat Shocked										
LYC-PC100	PC-12 Cell Lysate										
LYC-3T100	3T3 Cell Lysate										
SPA-227	Crystallin, α B (phospho-Ser59) Polyclonal										
960-074	Crystallin, α B ImmunoSet										

Background:

Alpha-crystallins composed of ~20 kDa α A and α B 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 tissues¹. Crystallin families β and γ share homology with each other but not the α family or the small heat shock protein (sHsp) family, while Alpha-crystallins, induced by heat in a variety of organisms, belong to the sHsp family². The α -crystallins possess structural and functional similarities and share sequence homology with Hsp25/27³. Most small heat shock proteins exhibit four common structural and functional features: molecular weight between 12 and 43kDa; the formation of large oligomeric complexes composed of α A-crystallin, α B-crystallin and Hsp25/27; a moderately conserved α -crystallin domain in the central region of the protein; and molecular chaperone activity^{2,4}. The α -crystallin domain bounded by variable N-terminal and C-terminal extensions contains approximately 90 residues and participates in oligomer assembly. Oligomers, potentially 800kDa or more, exhibit dynamic subunit exchanges and organizational plasticity, which may promote functional diversity. Phosphorylation of serine residues occurs in response to stress during development, typically decreasing oligomer size⁴. Chaperone activity requires oligomerization (which, in turn, modulates the chaperone activity) and is confined to binding unfolded intermediates to prevent irreversible aggregation^{2,4}, even though productive release and refolding of denatured proteins requires close cooperation with other chaperones. Other proposed functions include a role in membrane stabilization² and modulation of intermediate filament organization during physiological stress and neurodegenerative disease⁵. Alpha B Crystallin is phosphorylated on serine 45 by Erk1/2 and serine 59 by MAPKAPK-2^{6,7}. When isolated, mono-phosphorylated α B-crystallin contains only α B-crystallin phosphorylated at serine 19⁸.

References:

1. Augusteyn, R.C., *et al.* (1998) *Prog in Polymer Sci.* **23**, 375-413.
2. Narberhaus, F. (2002) *Microbiol Mol Biol Rev.* **66**, 64-93.
3. Merck, K.B., *et al.* (1993) *J Biol Chem.* **268**, 1046-1052.
4. MacRae, T.H. (2000) *Cell Mol Life Sci.* **57**, 899-913.
5. Head, M.W., *et al.* (2000) *Neuropathol Appl Neurobiol.* **26**, 304-312.
6. Armstrong, S.C., *et al.* (2000) *J Mol Cell Cardiol.* **32**, 1301-1314.
7. Ito H, *et al.* (1999) *FEBS Lett.* **446**, 269-272.
8. Kamei, A., *et al.* (2001) *Biol Pharm Bull.* **24**, 96-99.



Western Blot Analysis of Crystallin α B (phospho-Ser45):

Lane 1: MWM, Lane 2: α B1 Crystallin Protein (phospho), Lane 3: α B2 Crystallin Protein (non-phospho)

Generally reagents are good for one year from the date of receipt, except for conjugates which are good for six months and reagents with an expiration date indicated on the label or other supporting document.

FOR RESEARCH USE ONLY; NOT FOR THERAPEUTIC OR DIAGNOSTIC USE

5777 Hines Drive • Ann Arbor, MI • 48108 | Tel: 800-833-8651 or 800-668-6113 | Fax: 734-668-2793
www.assaydesigns.com | orders@assaydesigns.com | technical@assaydesigns.com