

HSP110 polyclonal antibody

Hsp 110 belongs to a family of large stress proteins known as the Hsp110/SSE family, the mostly overlooked relatives of the well studied Hsp70 family. Members of this group, while significantly larger and containing sequences not present in the Hsp70 family, share 30-33% amino acid identity with members of the Hsp70 family, mostly in the conserved ATP-binding domain. Presently, all identified members of the Hsp70-like sequences belong to either the Hsp70 family or the Hsp110/SSE family. Hsp110 appears to be a normal constituent of mammalian cells with highest expression in the brain, and its induction by heat shock strongly correlates with cellular thermotolerance. Hsp110 may contribute to RNA stability, and its capacity to bind peptide has made Hsp110 a promising immunoadjuvant in anti-tumor vaccine development.

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

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-SPA-1101-E	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

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

Alternative Name	HSP105, HspH1, Heat shock protein 105, Heat shock protein 110
Application	WB
Application Notes	Detects a band of ~110kDa by Western blot.
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
GenBank ID	Z47807
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of hamster HSP110. The sequence is completely conserved in human and mouse.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	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 rabbit serum.
Species Reactivity	Bovine, Hamster, Human, Monkey, Mouse, Rat, Sheep, Yeast
UniProt ID	Q60446

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

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

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

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

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

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