

CHIP polyclonal antibody

CHIP has been shown to interact both with the constitutive form of hsc70 and the stress inducible form, hsp70. This novel 35 kDa cytoplasmic protein has been shown to be highly expressed in striated muscle in vivo. Additional studies have shown that this protein is expressed over a broad range of cultured tissues. Through immunoprecipitation experiments, CHIP has been shown to directly bind to the carboxyl terminus of hsc70 and hsp70 where it decreases ATPase activity and reduces overall chaperone efficiency.

CHIP has also been identified as an important protein in the ubiquitin-proteasome system. CHIP contains a U-box domain and acts as an E3 ubiquitin-ligase in conjunction with hsc70 and hsp90.

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

Ordering Information

[Order Online »](#)

ENZ-ABS273-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Carboxyl terminus of Hsc70-interacting protein

Application WB

Application Notes Detects a band of ~35kDa by Western blot.

Formulation Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 218-232 of human CHIP.

Isotype IgG

Purity Detail Antigen affinity chromatography

Species Reactivity Human, Mouse

UniProt ID Q9UNE7

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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@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)