

# UNC45a monoclonal antibody (AbS1)

GC UNC45 (general cell UNC45/unc45a) is an Hsp90 co-chaperone that has been shown to play a critical role in the chaperoning of myosin proteins as well as the progesterone receptor (PR). There are two known vertebrate isoforms of UNC45 that are functional orthologs of the *C. elegans* unc-45, an important protein in the development of organized body wall muscles. The vertebrate general cell form (GC UNC45/unc45a) is ubiquitously expressed, while the striated muscle-specific isoform (SM UNC45/unc45b) is found in cardiac and skeletal muscle. In vitro, GC UNC45 preferentially binds the Hsp90 beta isoform, and is specifically required for proper cellular distribution of Hsp90 $\beta$  but not Hsp90 $\alpha$ .

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

Citations: 7

[View Online »](#)

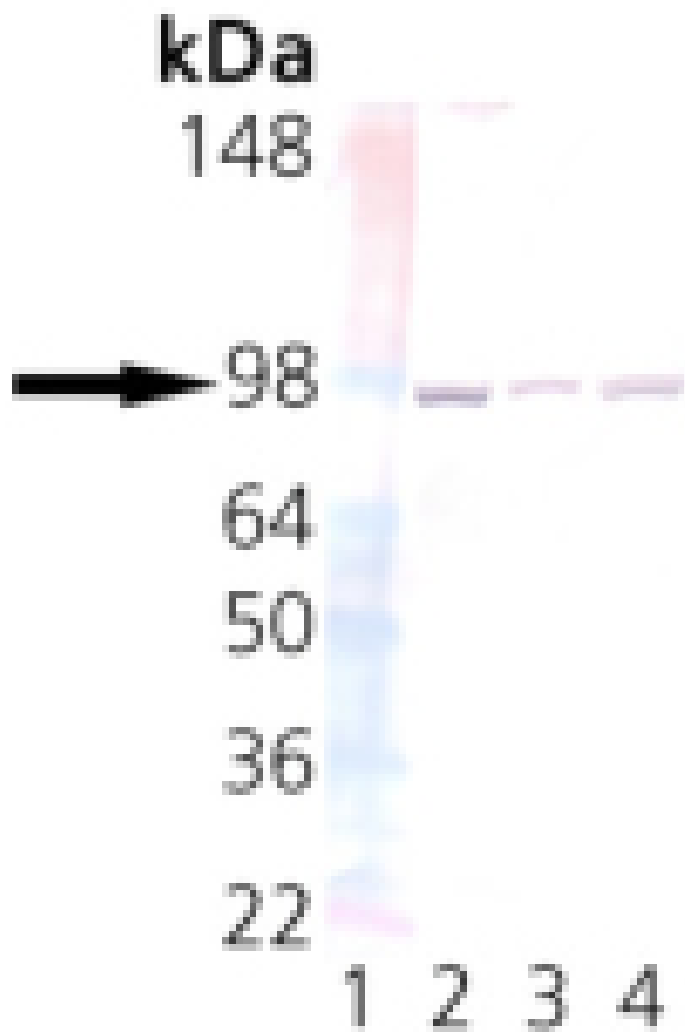
## Ordering Information

[Order Online »](#)

ADI-SRA-1800-D	50 $\mu$ g
ADI-SRA-1800-F	200 $\mu$ g

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of UNC45: Lane 1: MW marker, Lane 2: HeLa, Lane 3: 3T3, Lane 4: PC-12.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** WB

**Application Notes** Detects a band of ~98kDa by Western blot.

**Clone** AbS1

**Formulation** Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

**Host** Mouse

**Immunogen** Recombinant human UNC45.

**Isotype** IgG2b

**Purity Detail** Protein G 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 mouse ascites.

**Species Reactivity** Human, Mouse, Rat

**UniProt ID** Q9H3U1

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



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)