

# HSP47 monoclonal antibody (M16.10A1)

Hsp47, also known as colligin and gp46, is an ER-resident collagen-specific stress glycoprotein involved in the synthesis and assembly of various collagens as a molecular chaperone. Hsp47 is a member of the serpin superfamily, but localizes to the ER via binding of its C-terminal RDEL sequence to KDEL receptors in the ER. Hsp47 binds procollagen cotranslationally as it enters the ER, and dissociates prior to its secretion.

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

Citations: 67

[View Online »](#)

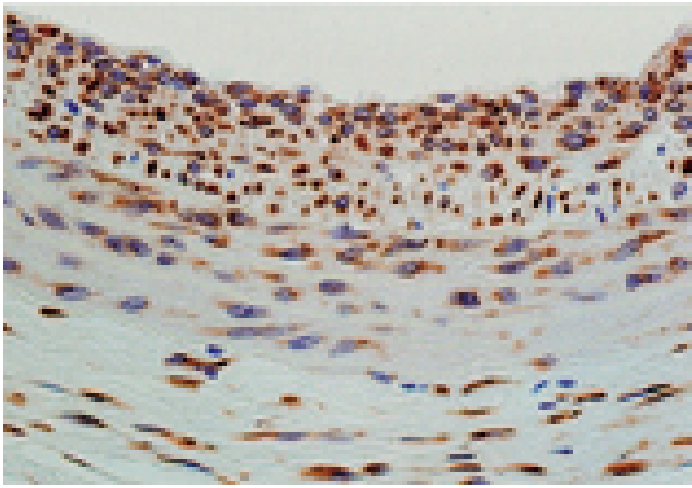
## Ordering Information

[Order Online »](#)

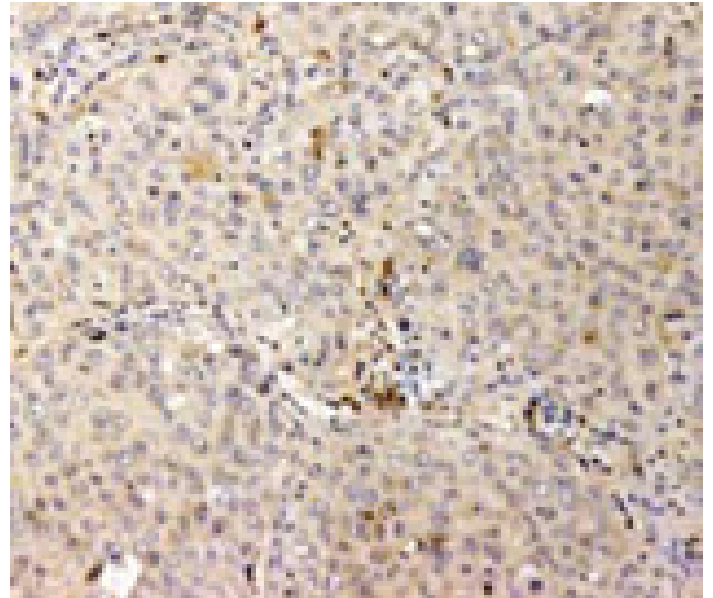
ADI-SPA-470-D	50µg
ADI-SPA-470-F	200µg

Manuals, SDS & CofA

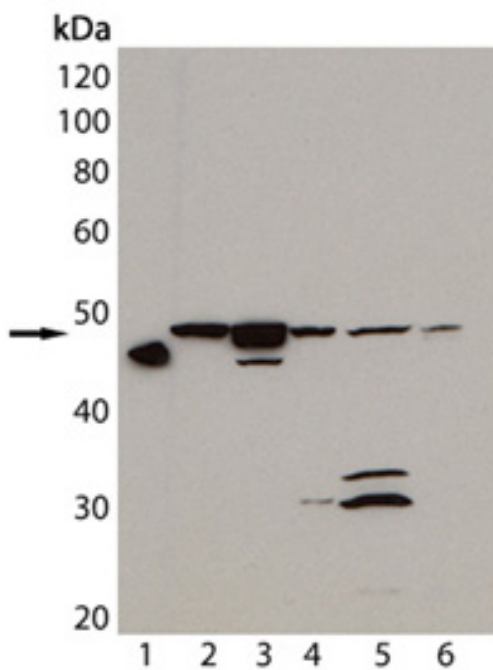
[View Online »](#)



Immunohistochemistry analysis of paraffin-embedded tissue section of rat carotid artery 14 days after balloon-injury to the artery, stained using Hsp47 (Colligin) mAb (M16.10A1).



Immunohistochemistry analysis of human breast cancer tissue immunohistochemically stained using Hsp47 mAb (M16.10A1).



Western blot analysis of HSP47, mAb (M16.10A1) (Prod. no. ADI-SPA-470): Lane 1: Rat recombinant HSP47; Lane 2: Rat-2 (heat shocked); Lane 3: L929 (heat shocked); Lane 4: 3T3 (heat shocked); Lane 5: CHO (heat shocked); Lane 6: HeLa (heat shocked).

## 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** SerpinH1, Colligin, Gp46, Heat shock protein 47, Collagen-binding protein

**Application** IHC, WB

**Application Notes** Detects a band of ~47kDa by Western blot.

**Clone** M16.10A1

**Formulation** Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

**GenBank ID** X61598

**Host** Mouse

**Immunogen** Native rat HSP47.

**Isotype** IgG2b $\lambda$

**Purity Detail** Protein G affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (1:50)Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Source** Purified from ascites.

**Species Reactivity** Bovine, Chicken, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep

Technical Info / Product Notes

Cited samples:

[For an overview on cited samples please click here.](#)

UniProt ID

P50454

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-  
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)