

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 co-translationally as it enters the ER, and dissociates prior to its secretion.

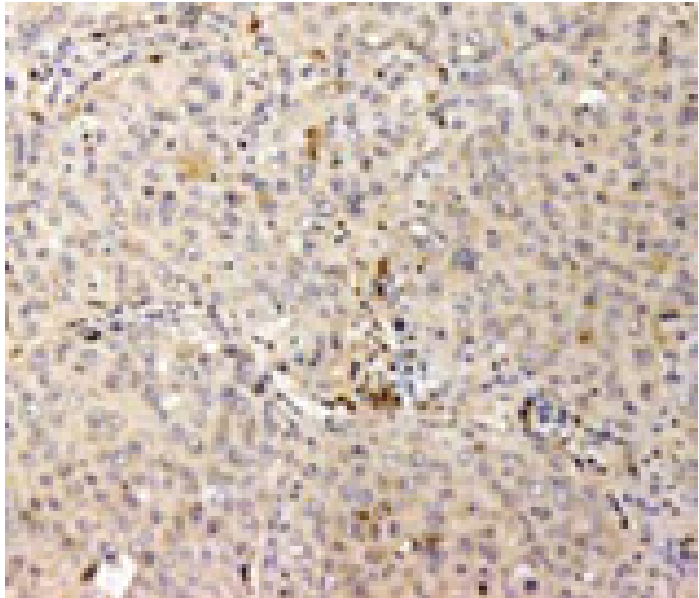
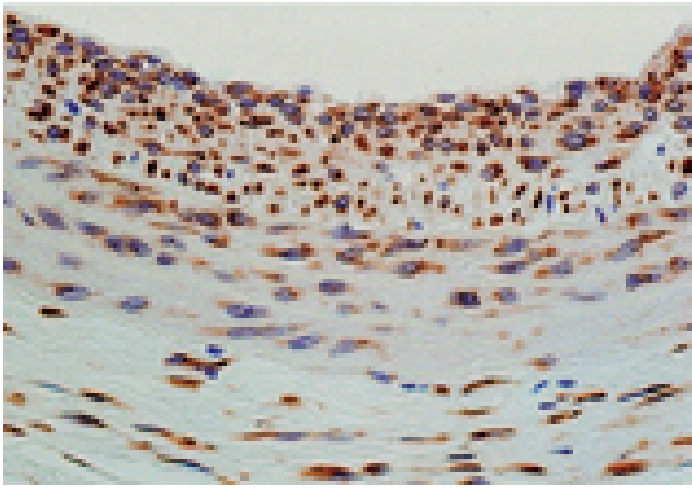
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 67 [View Online »](#)

Ordering Information [Order Online »](#)

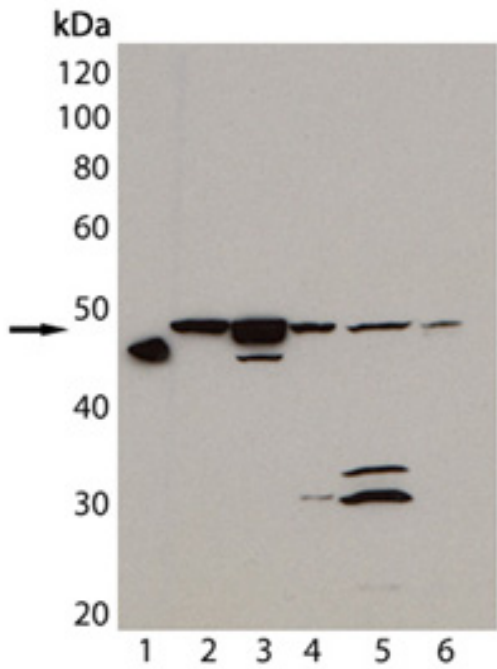
ADI-SPA-470-J	1mg
ADI-SPA-470-D	50µg
ADI-SPA-470-F	200µg

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	IgG2bλ
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)