

# HSP47 (human), (recombinant)

**Recombinant human HSP47 that acts as an ER-resident chaperone, assisting in collagen synthesis and assembly.**

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.

Citations: 1

[View Online »](#)

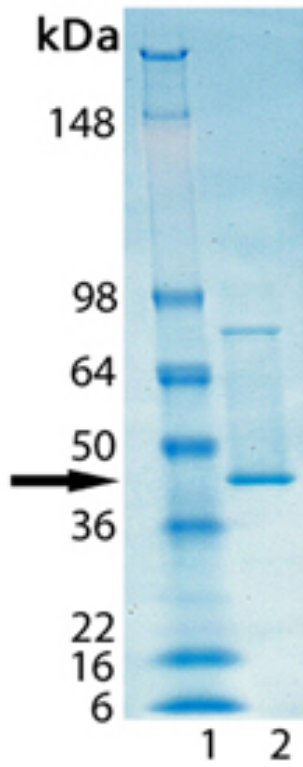
**Ordering Information**

[Order Online »](#)

ADI-SPP-535-D	50µg
---------------	------

**Manuals, SDS & CofA**

[View Online »](#)



SDS-PAGE Analysis. Lane 1: MW marker, Lane 2: 1.0  $\mu\text{g}$  of purified Hsp47 (Prod. No. ADI-SPP-535)

## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** SerpinH1, Colligin, Gp46, Heat shock protein 47, Collagen binding protein

**Application Notes** Western blot control.

**Endotoxin Content** < 50 EU/mg determined by Limulus Amebocyte Lysate (LAL) gel clot assay.

**Formulation** Liquid. In 30mM TRIS, pH 8.0, containing 100mM sodium chloride, 1mM DTT, and 0.4M arginine.

**MW** ~47kDa

**Purity** ≥90% (SDS-PAGE)

**Purity Detail** Purified by multi-step chromatography.

**Source** Produced in *E. coli*.

**UniProt ID** P50454

**Last modified: October 9, 2025**



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)