

HSP47 (human), (recombinant)

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 »](#)

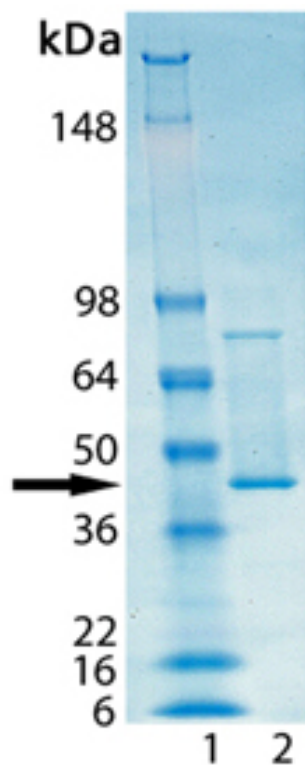
Ordering Information

[Order Online »](#)

| | |
|---------------|-------|
| ADI-SPP-535-D | 50µg |
| ADI-SPP-535-F | 200µg |

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis. Lane 1: MW marker, Lane 2: 1.0 μ 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



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-
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)