## 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 collagenspecific stress glycoprotein involved in the synthesis and assembly of varoius 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.

Citations: 1

View Online »

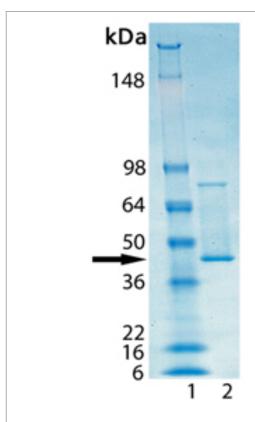
## **Ordering Information**

**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  $\mu g$  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

