## HSP40/Hdj1 (human), (recombinant)

Recombinant human HSP40/Hdj1 that acts as a co-chaperone to Hsp70, facilitating protein folding and preventing aggregation.

Hsp40/Hdj1 is a cytosolic co-chaperone belonging to the class DnaJ, named after its homolog in *E. coli*. Hsp40 functions in protein folding by binding nascent peptides and unfolded substrates and facilitating substrate interaction with Hsp70. Binding of Hsp40 to Hsp70 increases Hsp70 ATP hydrolysis and substrate binding, which is reversed by the action of nucleotide exchange factors (GrpE in *E. coli*). Repeated cycles of Hsp40/Hsp70 peptide binding and ATP hydrolysis prevent premature folding and aggregation, promoting protein maturation throughout the cell.

Citations: 9

**View Online »** 

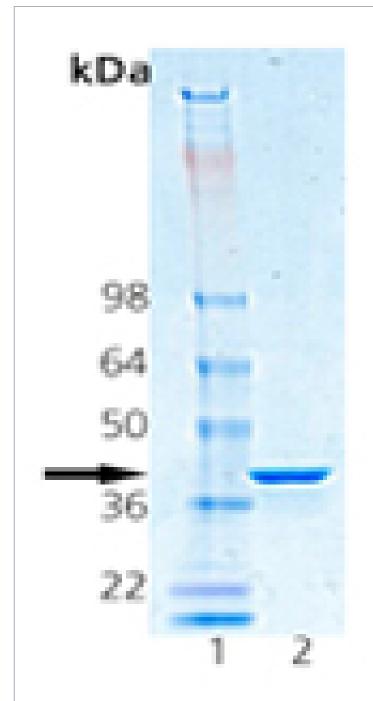
## **Ordering Information**

**Order Online** »

ADI-SPP-400-D	50µg
ADI-SPP-400-F	200µg

Manuals, SDS & CofA

**View Online** »



1 2 kDa 

SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 2.0µg of purified Human HSP40 Protein (Prod. No. ADI-SPP-400).

SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 2.0µg of purified Human HSP40 Protein (Prod. No .ADI-SPP-400).

## **Handling & Storage**

**Long Term Storage** -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

**Product Details** 

Alternative Name Heat shock protein 40, DnaJ protein homolog 1

Application Notes Western blot control.

**Formulation** Liquid. In Dulbecco's PBS.

MW ~50kDa

Purity ≥80% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Source Produced in E. coli. Contains N-terminal His tag.

UniProt ID P25685

Last modified: October 9, 2025

