

HSP40/Hdj1 (human), (recombinant)

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 »](#)

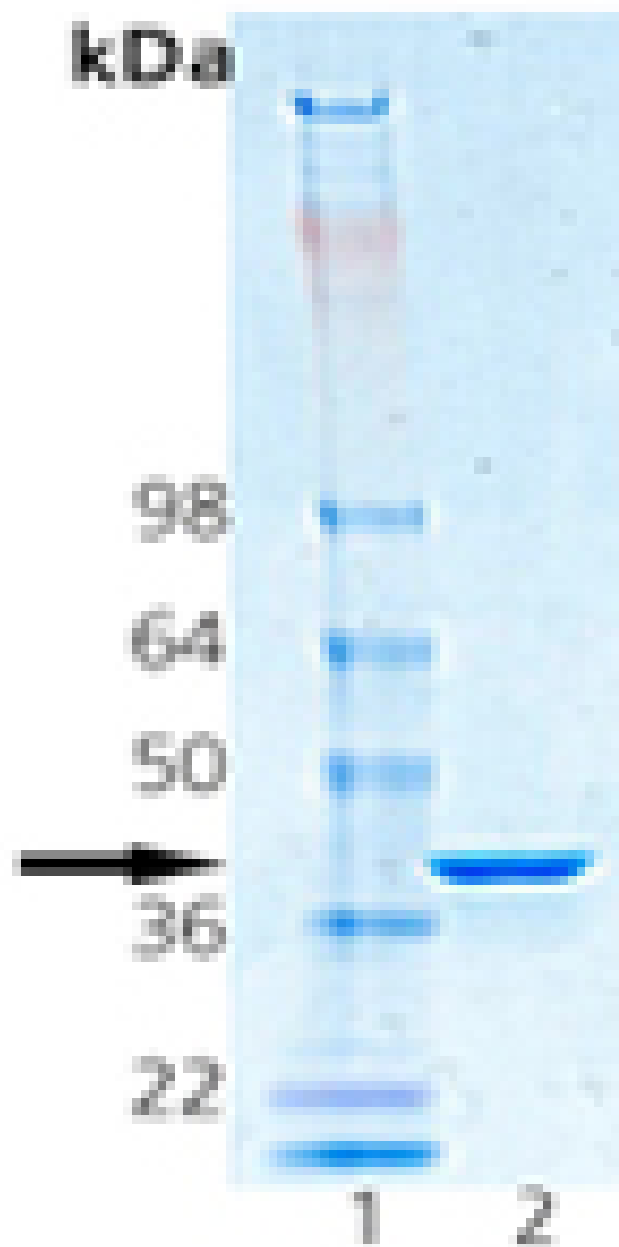
Ordering Information

[Order Online »](#)

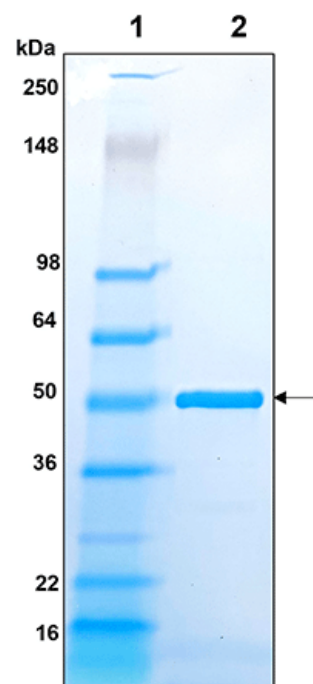
ADI-SPP-400-D	50µg
ADI-SPP-400-F	200µg

Manuals, SDS & CofA

[View Online »](#)



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: 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)