

# DnaJ (E. coli), (recombinant) (His- tag)

## Recombinant DnaJ (E. coli, His-tag) that acts as a molecular chaperone, cooperating with

**Escherichia coli** heat shock protein DnaJ belongs to the molecular chaperone class of proteins. Located in an operon with *dnaK*, the *E. coli* *dnaJ* gene shares sequence identity with eukaryotic cytosolic and endoplasmic reticulum DnaJ homologs (Hsp40) involved in protein folding, membrane translocation of newly synthesized proteins, and initiation of translation. The N-terminal, highly-conserved "J" domain of the DnaJ protein supports interaction with the DnaK protein and represents the signature sequence of DnaJ family members. Bacterial DnaJ protein acts synergistically with bacterial chaperones DnaK (Hsp70 homolog) and GrpE in various functions, including suppressing eukaryotic and prokaryotic polypeptide aggregation to promote protein folding, facilitating protein translocation through intracellular compartments or protein secretion, and repairing and reactivating partially aggregated enzymes such as *E. coli* RNAP and luciferase. In these reactions, DnaJ protein facilitates the binding of DnaK to its substrate by accelerating DnaK-catalyzed ATP hydrolysis to produce a DnaK-ADP form with a higher affinity for certain protein substrates. Some eukaryotic homologues of DnaJ also display an ability to stimulate the ATPase activity of Hsp70 proteins, including DnaJ homologue *S. cerevisiae* YDJ1 which interacts with DnaK homologue Ssa1 to stimulate its ATPase activity. Data suggests that DnaJ protein acting alone possesses chaperone activity, demonstrating an ability to bind tightly to denatured proteins such as rhodanese or luciferase to prevent their aggregation or misfolding.

Citations: 5

[View Online »](#)

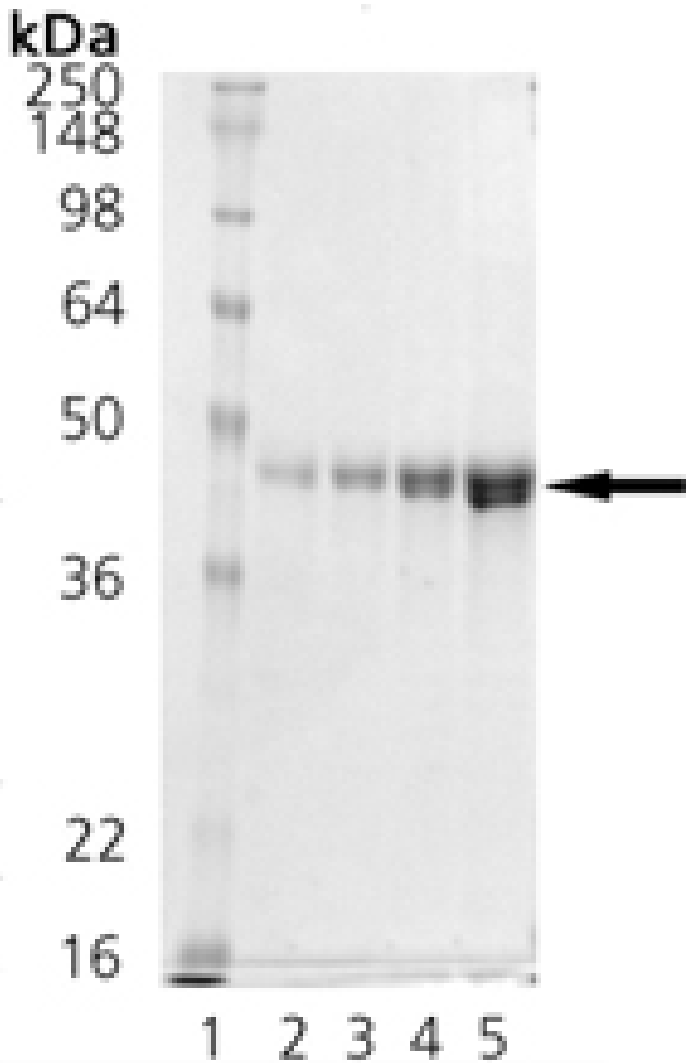
## Ordering Information

[Order Online »](#)

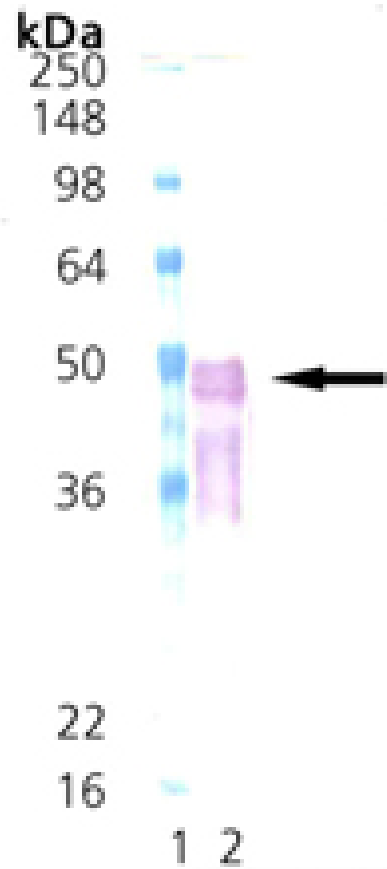
ADI-SPP-640-D	50µg
ADI-SPP-640-F	200µg

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MWM, Lane 2: 0.5μg, Lane 3: 1μg, Lane 4: 2μg, Lane 5: 5μg DnaJ Recombinant Protein.



Western Blot Analysis: Lane 1: MWM, Lane 2: 100 ng of DnaJ probed with DnaJ pAb.

## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** DNAJB2, DnaJ HSJ1, HSPF3, HSP40

**Application Notes** Western blot control.

**Formulation** Liquid. In 25mM HEPES, pH 7.5, containing 200mM potassium chloride, 10% glycerol, and 5mM DTT.

**MW** ~43kDa

**Purity** ≥90% (SDS-PAGE; Western blot)

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

**Source** Produced in *E. coli*. *E. coli* DnaJ is fused at the N-terminus to a His-tag.

**UniProt ID** P08622

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)