DnaJ (E. coli), (recombinant) (Histag)

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

Escharichia poti that shock protein Protein Protein the molecular chaperone class of proteins. Located in an operon with dnaK, theE. coli dnaJ gene shares sequence identity with eukaryotic cytosolic and **translania tinulan drafoldings Alspin folded** in protein folding, membranestranslocation 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 asE. 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 Ssal 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 »

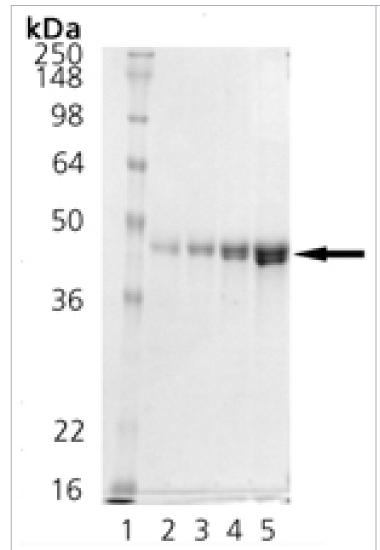
Ordering Information

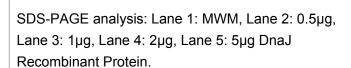
Order Online »

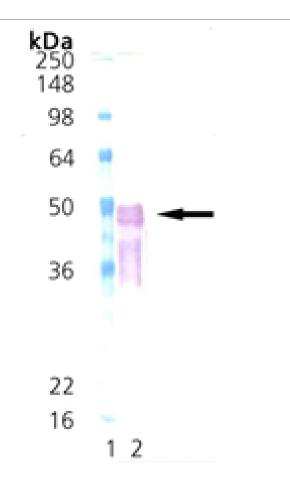
ADI-SPP-640-J	1mg
ADI-SPP-640-D	50µg
ADI-SPP-640-F	200µg

Manuals, SDS & CofA

View Online »







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.

Produced in E. coli. E. coli DnaJ is fused at the N-terminus Source

to a His-tag.

P08622 **UniProt ID**

Last modified: October 9, 2025



info-

eu@enzolifesciences.com