

# DnaJ (E. coli)

## polyclonal antibody

DnaJ is a basic 41 kDa *E. coli* Heat shock protein which belongs to the molecular chaperone class of proteins. Bacterial DnaJ protein acts synergistically with the bacterial chaperone DnaK (Hsp70 homologue) and the other co-chaperone GrpE in suppressing eukaryotic and prokaryotic polypeptide aggregation, facilitating protein translocation through intracellular compartments or protein secretion, reactivating some partially aggregated enzymes, and activating pre-priming complex during initiation of DNA replication.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 7

[View Online »](#)

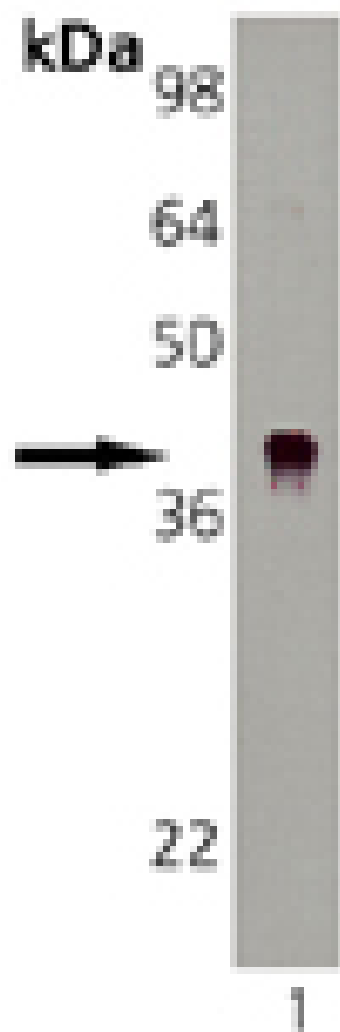
### Ordering Information

[Order Online »](#)

ADI-SPA-410-D	50µg
ADI-SPA-410-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: Lane 1: DnaJ Recombinant *E. coli* Protein.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** HSP40, Heat shock protein 40, Heat shock protein J

**Application** IP, WB

**Application Notes** Detects a band of ~41kDa by Western blot.

**Formulation** Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

**Host** Rabbit

**Immunogen** E. coli DnaJ.

**Purity Detail** Protein A affinity purified.

**Recommendation** Immunoprecipitation (1:200)Western Blot (1:1,000, ECL)Suggested dilutions/conditions  
**Dilutions/Conditions** may not be available for all applications.Optimal conditions must be determined individually for each application.

**Source** Purified from rabbit serum.

**Species Reactivity** E. coli

**UniProt ID** P08622

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).



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)