

# GrpE (*E. coli*) polyclonal antibody

GrpE (Glucose Regulated Protein *E. coli*) is encoded by a monocistronic *E. coli* gene, which is under the control of both s32 and s70 transcription factors. The GrpE protein is a so-called co-chaperone because it is known to assist the DnaK (Hsp70) protein to effectively carry out DnaK-dependent chaperone activity (i.e. protein folding, protein transport, disaggregation of heat inactivated proteins, activation of mutant protein).

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

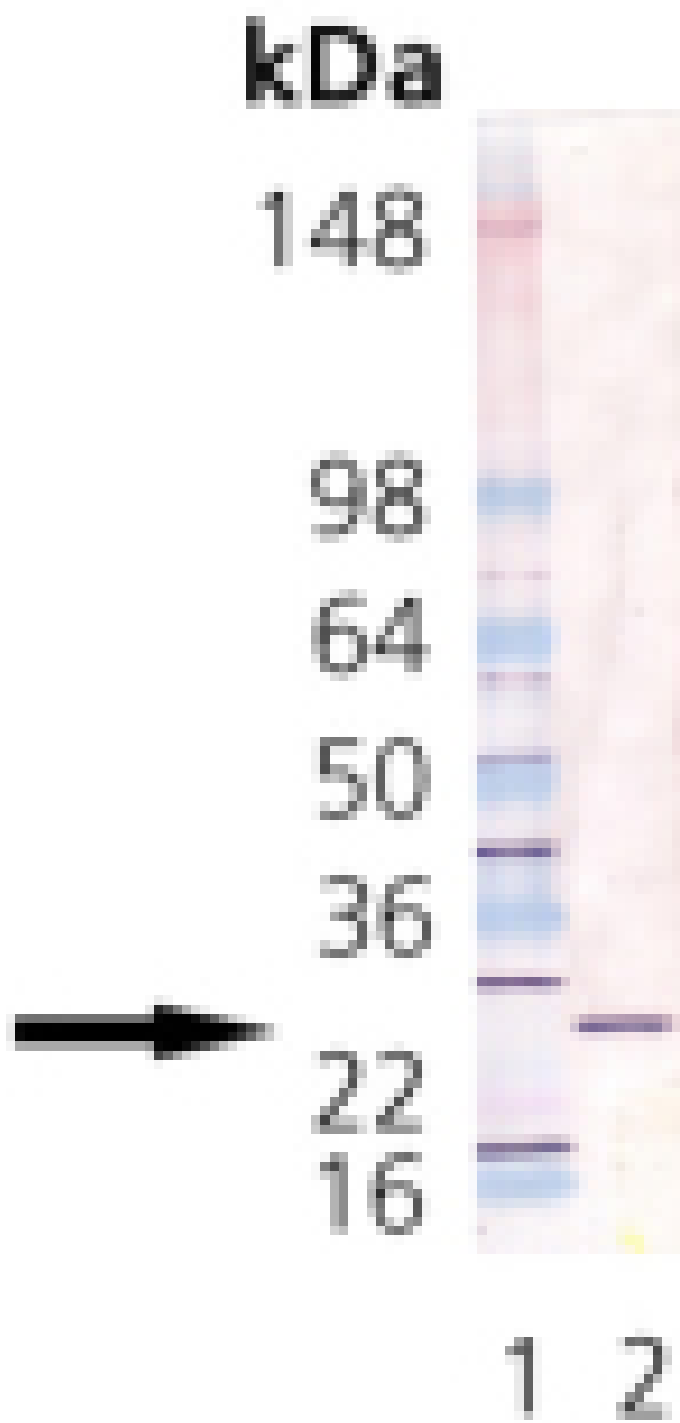
## Ordering Information

[Order Online »](#)

ADI-SPA-240-J	1mg
ADI-SPA-240-0200	200µg

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of GrpE: Lane 1: MW marker,  
Lane 2: GrpE 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** HSP70 cofactor, HSP24

**Application** WB

**Application Notes** Detects a band of ~24kDa by Western blot.

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

**GenBank ID** X07863

**Host** Rabbit

**Immunogen** Purified E. coli GrpE.

**Purity Detail** Protein A affinity purified.

**Recommendation Dilutions/Conditions** Western Blot (1:1,000, colorimetric) Suggested 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** P09372 (strain K12)

**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)