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.

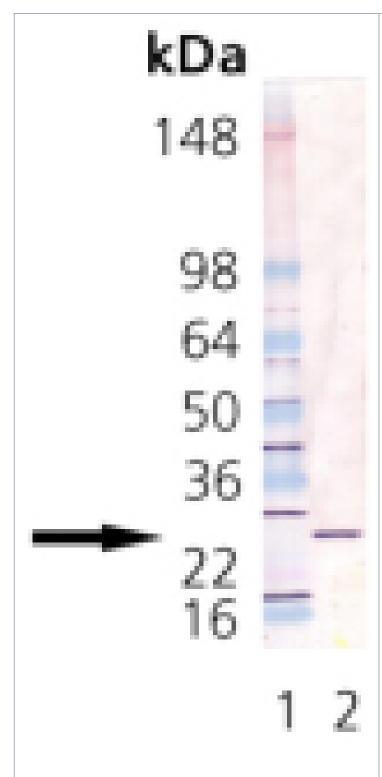
Ordering Information

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 NotesDetects 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/ConditionsWestern 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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com