

GroES (*E. coli*) polyclonal antibody

GroEL and GroES are *E. coli* chaperonins, which are homologs of mitochondrial chaperonins Hsp60 and Hsp10. GroEL is a double toroidal assembly of 14 identical subunits which form two heptameric rings stacked back-to-back, with a cavity at each end. GroEL and its co-chaperonin GroES facilitate protein folding with an ATP-dependent mechanism.

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

Citations: 4

[View Online »](#)

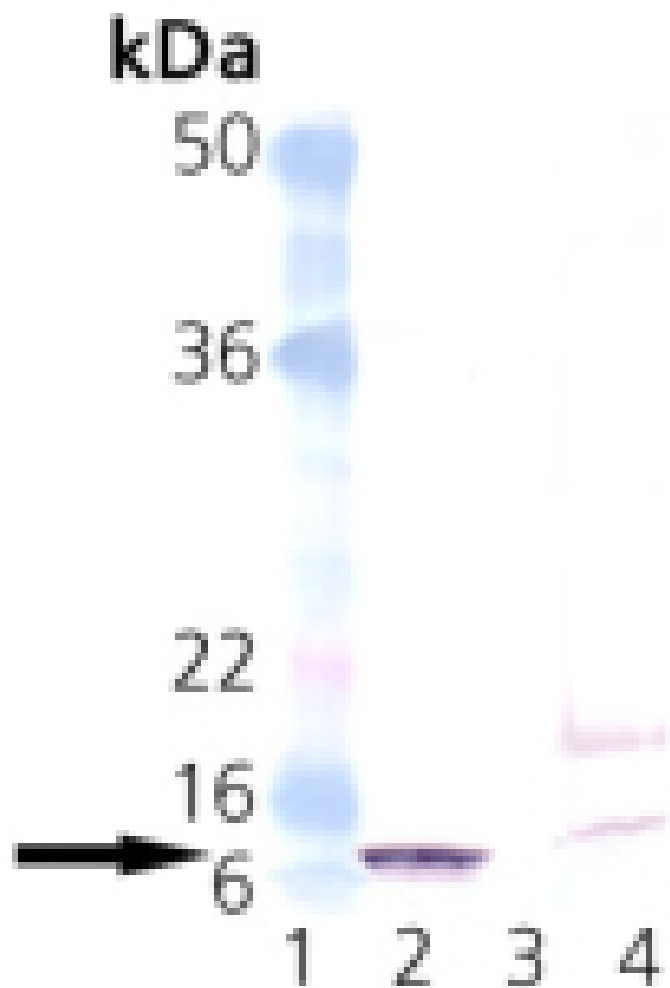
Ordering Information

[Order Online »](#)

ADI-SPA-210-D	50µl
ADI-SPA-210-F	200µl

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of GroES: Lane 1: MW marker, Lane 2: GroES Recombinant *E. coli* Protein, Lane 3: (Empty), Lane 4: *E. coli* Lysate.

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 Chaperonin 10

Application WB

Application Notes Detects a band of ~10kDa by Western blot.

Formulation Liquid. Whole rabbit serum.

GenBank ID X07850

Host Rabbit

Immunogen E. coli GroES.

Recommendation Western Blot (1:5,000, colorimetric) Suggested dilutions/conditions may not be available
Dilutions/Conditions for all applications. Optimal conditions must be determined individually for each application.

Source Whole rabbit serum.

Species Reactivity E. coli

UniProt ID P0A6F9 (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)