

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: 5

[View Online »](#)

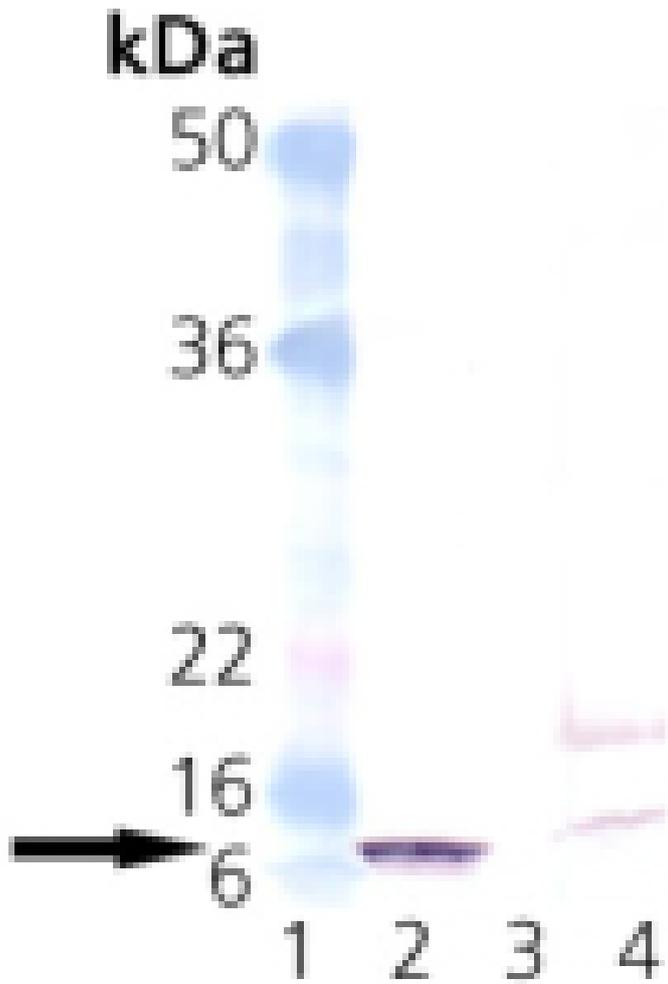
Ordering Information

[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 Dilutions/Conditions	Western Blot (1:5,000, colorimetric)Suggested dilutions/conditions may not be available 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](#)