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 toriodal 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 »

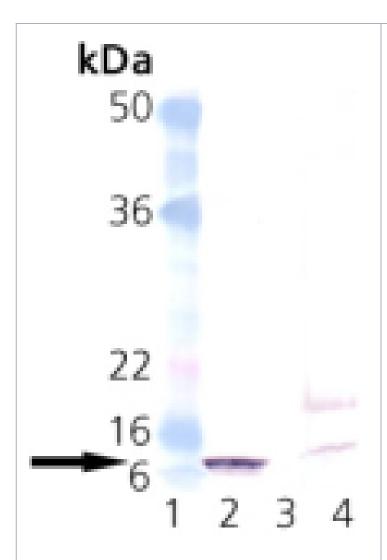
Ordering Information

Order Online »

ADI-SPA-210-D	50µI
ADI-SPA-210-F	200μΙ

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 NotesDetects 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

