GroEL (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: 11

View Online »

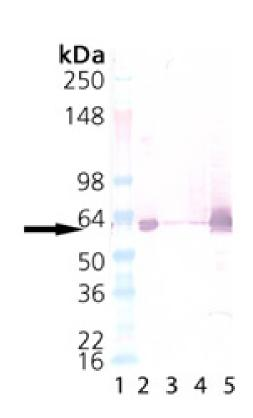
Ordering Information

Order Online »

ADI-SPS-875-D	50µg
ADI-SPS-875-F	200μg

Manuals, SDS & CofA

View Online »



Western Blot Analysis of GroEL: Lane 1: MW Marker, Lane 2: GroEL (E. coli), (recombinant) (Prod. No. ADI-SPP-610), Lane 3: HSP60 (human), (recombinant) (Prod. No. ADI-NSP-540), Lane 4: HeLa (Heat Shocked), Cell Lysate (Prod. No. ADI-LYC-HL101), Lane 5: E. coli Cell Lysate all probed with GroEL (*E. coli*), pAb (Prod. No. ADI-SPS-875).

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 60

Application WB

Application NotesDetects a band of ~60kDa by Western blot.

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

azide.

Host Rabbit

Immunogen E. coli GroEL.

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 P0A6F5 (strain K12)

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 infouk@enzolifesciences.com