

GroEL (*E. coli*) monoclonal antibody (9A1/2)

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: 12

[View Online »](#)

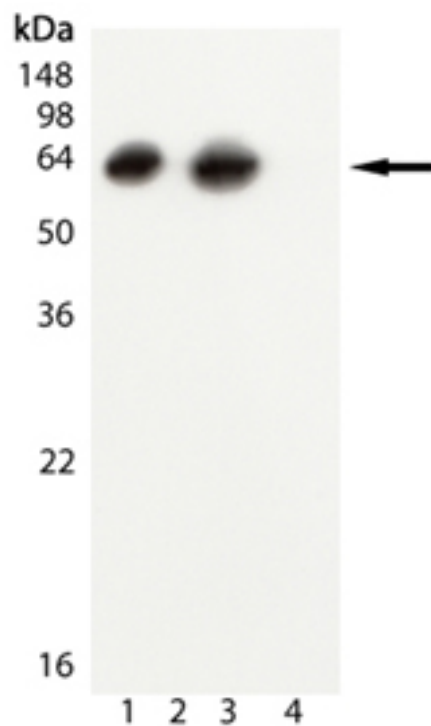
Ordering Information

[Order Online »](#)

ADI-SPS-870-D	50µg
ADI-SPS-870-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of GroEL, mAb (9A1/2) (Prod. No. ADI-SPS-870): Lane 1: GroEL (E. coli) (rec) (Prod. No. ADI-SPP-610) Protein, Lane 2: HSP60 (Human) (rec) (negative control) (Prod. No. ADI-SPP-540), Lane 3: E. coli Lysate, Lane 4: HeLa Lysate, heat shocked (negative control) (Prod. No. ADI-LYC-HL101)

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	IP, WB
--------------------	--------

Application Notes	Detects a band of ~60kDa by Western blot.
--------------------------	---

Clone 9A1/2

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

GenBank ID X07899

Host Mouse

Immunogen E. coli GroEL.

Isotype IgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions	Western Blot (1:10,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
--	---

Source	Purified from mouse ascites.
---------------	------------------------------

Species Reactivity E. coli

UniProt ID P0A6F5 (strain K12)

Worry-free Guarantee

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com