

# 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](#).

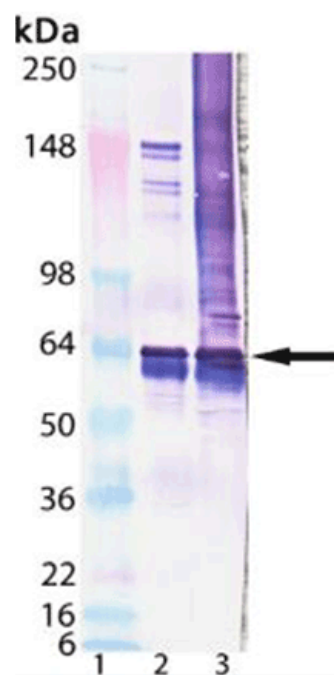
## Ordering Information

[Order Online »](#)

ENZ-ABS714-0200	200µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of ENZ-ABS714: Lane 1: MW Marker, Lane 2: GroEL (*E. coli*), (recombinant) (Prod. No. ADI-SPP-610), Lane 3: *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 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.
Host	Mouse
Immunogen	E. coli GroEL.
Isotype	IgG1κ
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from hybridoma tissue culture supernatant.
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](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)