

# 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](#).

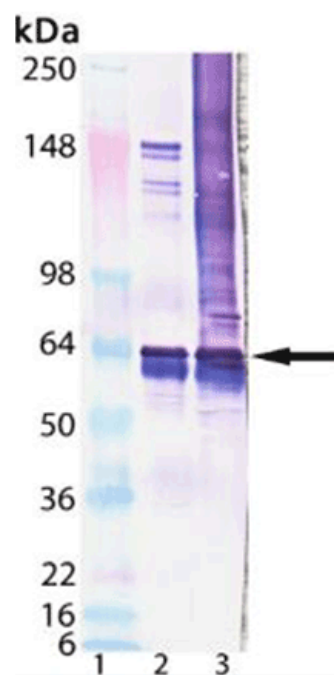
## Ordering Information

[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** Western Blot (1:1000, colorimetric) Suggested dilutions/conditions may not be available  
**Dilutions/Conditions** 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)



ENZO LIFE SCIENCES,  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 1 42 42 42 42  
[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)

**This antibody is covered by our Worry Free Guarantee**