

HSP60 (bacterial) monoclonal antibody (A57-E4)

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-0710	100µl
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

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 Heat shock protein 60 (bacterial)

Application WB

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

Clone A57-E4

Crossreactivity Does not cross-react with human HSP60.

Formulation Liquid. Ascites diluted in PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Recombinant *Chlamydia trachomatis* HSP60 (heat shock protein 60).

Isotype IgG1

Recommendation Dilutions/Conditions Western Blot (1:5,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity *Chlamydia*, *Coxiella burnetti*, *E. coli*, *Neisseria gonorrhoeae*, *Salmonella typhimurium*

Specificity Recognizes bacterial HSP60 from *Chlamydia*, *E. coli* (GroEL), *Salmonella typhimurium*, *Neisseria gonorrhoeae*, and *Coxiella burnetti*.

UniProt ID P0C0Z7

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)