

HSP60 (bacterial) monoclonal antibody (A57-B9)

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

Citations: 16

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-072-R100	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	ICC, IP, WB
Application Notes	Detects a band of ~60kDa by Western blot.
Clone	A57-B9
Crossreactivity	Minimally cross-reacts with <i>Borrelia burgdorfi</i> HSP60. Does not cross-react with <i>E. coli</i> GroEL or human HSP60.
Formulation	Liquid. Ascites diluted in PBS containing 0.05% sodium azide.
Host	Mouse
Immunogen	Recombinant <i>Chlamydia trachomatis</i> HSP60 (heat shock protein 60).
Isotype	IgG1
Recommendation Dilutions/Conditions	Immunocytochemistry (1:500)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Chlamydia
Specificity	Recognizes all three genera of <i>Chlamydia</i> HSP60. Epitope mapping studies suggest binding of aa 517-522.
UniProt ID	P0C0Z7

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)