

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 *Borrelia burgdorfi* HSP60. Does not cross-react with *E. coli* GroEL or 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 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 *Chlamydia* HSP60. Epitope mapping studies suggest binding of aa 517-522.



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-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@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

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee.