

# HSP65

## (mycobacterial)

## monoclonal antibody

## (4H11)

Hsp65 is a member of the Hsp60 family of heat shock proteins isolated from *Mycobacterium bovis* BCG. Hsp65 from *M. bovis* is identical to that of *M. tuberculosis*, and similar to that of *M. leprae*, two highly pathogenic strains of mycobacterium. Hsp65 is the immunodominant antigen during mycobacterial infection and vaccination, and has been linked to the development of autoimmune adjuvant arthritis in rats.

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

Citations: 2

[View Online »](#)

### Ordering Information


[Order Online »](#)

ADI-SPA-882-E	100µg
---------------	-------

### Manuals, SDS & CofA

[View Online »](#)

66  
45  
31  
21



← *Hsp65*

Western blot analysis  
of *M. leprae* cell lysate,  
probed with SPA-882.

## 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 65

**Application** in vitro Assay, WB

**Clone** 4H11

**Formulation** Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

**Host** Mouse

**Immunogen** Recombinant M. bovis BCG Hsp65

**Isotype** IgG1

**Purity Detail** Protein G affinity purified.

**Recommendation Dilutions/Conditions** Western Blot (1µg/ml, colorimetric) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from mouse ascites.

**Species Reactivity** Bovine, Mycobacterium bovis, Mycobacterium leprae

**UniProt ID** P0A521

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



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)