

HSP65

(mycobacterial)

monoclonal antibody

(3F7)

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: 1

[View Online »](#)

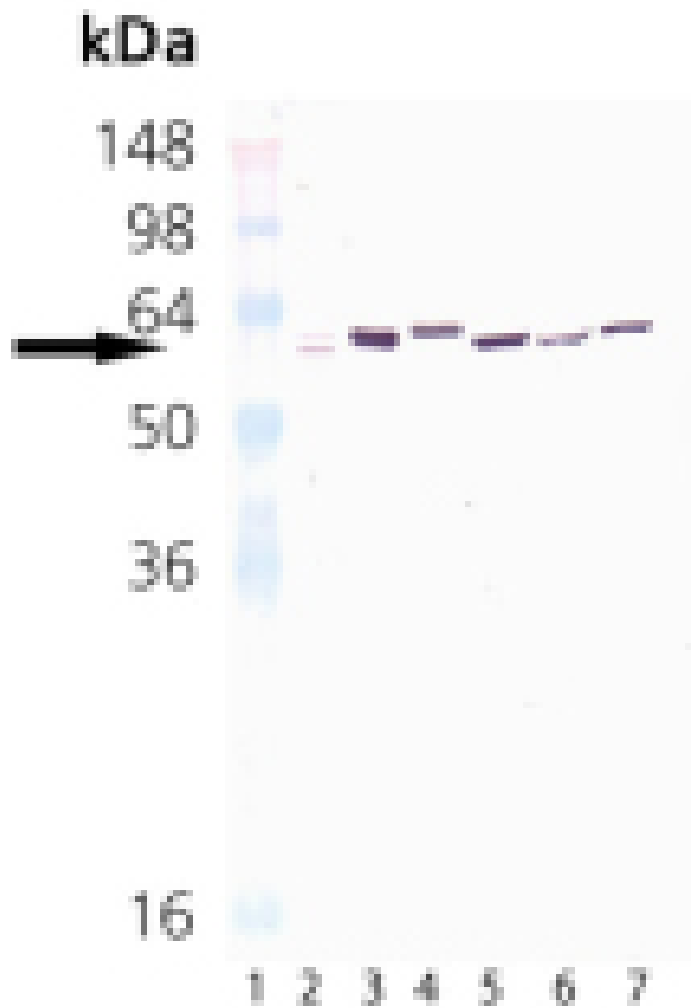
Ordering Information

[Order Online »](#)

ADI-SPA-881-D	50µl
ADI-SPA-881-F	200µl

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HSP65: Lane 1: MW marker, Lane 2: GroEL Recombinant E. coli Protein Lane 3: HSP60 Recombinant Human Protein, Lane 4: HSP65 Recombinant Mycobacterium bovis Protein, Lane 5: HeLa (HS) Cell Lysate, Lane 6: PC-12 Cell Lysate, Lane 7: 3T3 Cell Lysate.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Heat shock protein 65
Application	WB
Application Notes	Detects a band of ~65kDa by Western blot.
Clone	3F7
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Mouse
Immunogen	Recombinant M. bovis HSP65.
Isotype	IgM
Purity Detail	Ammonium sulfate precipitation purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, 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, Cryptosporidium, Drosophila, E. coli, Guinea pig, Hamster, Human, Mouse, Mycobacterium bovis, Mycobacterium leprae, Rat
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)