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 »

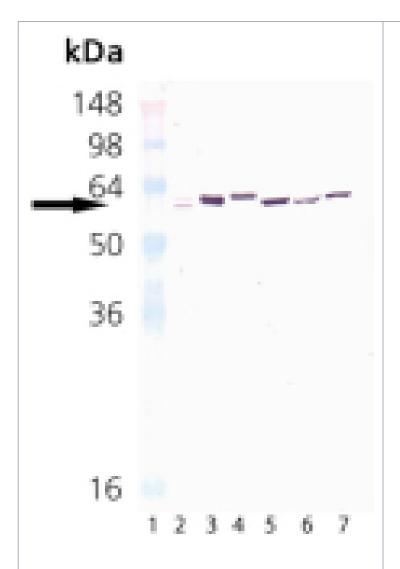
Ordering Information

Order Online »

ADI-SPA-881-D	50µI
ADI-SPA-881-F	200μΙ

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 NotesDetects 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.

lgM lgM

Purity Detail Ammonium sulfate precipitation purified.

Recommendation Dilutions/ConditionsWestern 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

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

