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

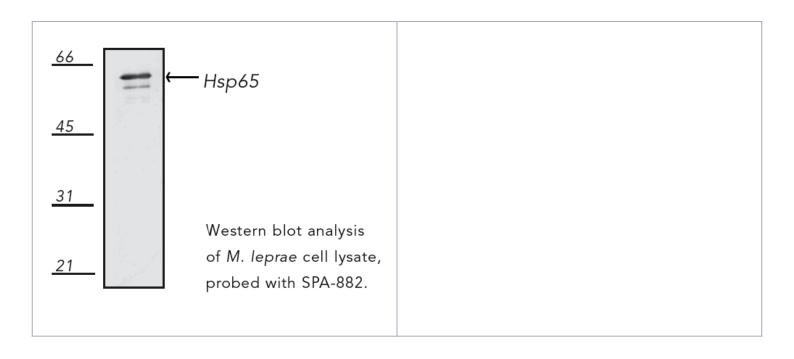
Order Online »

ADI-SPA-882-E

100µg

Manuals, SDS & CofA

View Online »



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

lsotype lgG1

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com