## HSP65 (low endotoxin) (M. bovis), (recombinant)

Recombinant M. bovis HSP65 (low endotoxin) that serves as an

Hep65 an 65 kDa protein isolated from Myropaeterium boyisi BCG, is a member of the hsp60 family of heat shock proteins. Sequence comparison of Hsp65 from different mycobacterium strains showed that the protein sequence of M. bovis BCG is identical to that of M. tuberculosis, and very similar to that of M. leprae, the pathogens that cause tuberculosis and tuberculoid leprosy, respectively. M. bovis BCG Hsp65 was identified as the immunodominant antigen during mycobacterial diseases and vaccination.

Citations: 1

**View Online »** 

## **Ordering Information**

**Order Online** »

ADI-ESP-581-D	50µg
ADI-ESP-581-F	200µg

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

**Long Term Storage** -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

**Product Details** 

Alternative Name Heat shock protein 65

**Application Notes** ATPase activity assay (positive). Western blot control.

**Endotoxin Content** <50EU/mg purified protein (LAL test)

**Formulation** Liquid. In 1x Dulbecco's PBS.

MW ~65kDa

Purity ≥95% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

**Source** Produced in *E. coli*.

UniProt ID P0A521

Last modified: October 9, 2025

