

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

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

**Immunodominant antigen in mycobacterial infection and vaccination.**  
Hsp65, a 65 kDa protein isolated from Mycobacterium bovis 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
---------------	------

## 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 <i>E. coli</i> .
UniProt ID	P0A521

Last modified: October 9, 2025



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)