

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

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
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



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)