
mpt64 Polyclonal Antibody

(Catalog # A57629)

Background

The present invention relates to fusion proteins containing at least two Mycobacterium species antigens. In particular, it relates to nucleic acids encoding fusion proteins that include two or more individual M. tuberculosis antigens, which increase serological sensitivity of sera from individuals infected with tuberculosis, and methods for their use in the diagnosis, treatment, and prevention of tuberculosis infection.

Description

mpt64 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300. 50% Glycerol, 0.01M PBS, pH 7.4.

Specificity

Mycobacterium tuberculosis

Isotype

IgG

Uniprot ID

P9WIN9

Purification

Protein G purified

Immunogen

Recombinant Mycobacterium tuberculosis Immunogenic protein MPT64 protein (24-228AA)

Storage

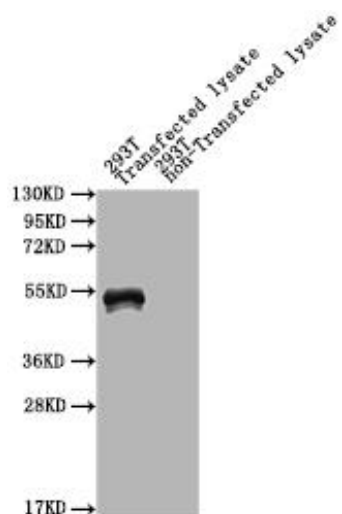
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Antigen MPT64, mpt64, Rv1980c, MT2032, MTCY39.39

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western Blot

Positive WB detected in: mpt64 293T Transfected lysate, 293T non-Transfected lysate

All lanes: mpt64 antibody at 1:2000

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 51 kDa

Observed band size: 51 kDa