
mpt64 Polyclonal Antibody

(Catalog # A55719)

Description

mpt64 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 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, MTCY39.39

Application

ELISA
