
MT1F Polyclonal Antibody

(Catalog #A64184)

Background

Metallothioneins have a high content of cysteine residues that bind various heavy metals; these proteins are transcriptionally regulated by both heavy metals and glucocorticoids.

Description

MT1F Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P04733

Purification

Protein G purified

Immunogen

Recombinant Human Metallothionein-1F protein (1-61AA)

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

Metallothionein-1F, MT-1F, Metallothionein-IF, MT-IF, MT1F, PRO0376

Application

ELISA; Recommended dilution: ELISA 1:1000
