

---

## MT1B Polyclonal Antibody

(Catalog #A64188)

---

### 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

MT1B 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

P07438

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Metallothionein-1B 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-1B, MT-1B, Metallothionein-IB, MT-IB, MT1B, MT1Q

### Application

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

---