
MT1H Polyclonal Antibody

(Catalog # A59938)

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

MT1H 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

P80294

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Metallothionein-1H protein (1-59AA)

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-1H, MT-1H, MT1H, Metallothionein-0, MT-0, Metallothionein-IH, MT-IH

Application

ELISA
