

MT1G Polyclonal Antibody

(Catalog #A57673)

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

MT1G Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P13640

Purification

Protein G purified

Immunogen

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

Storage

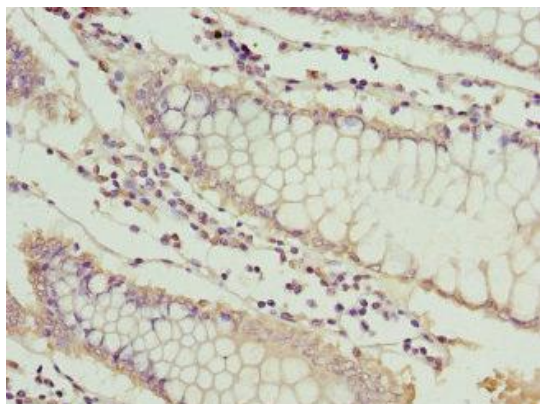
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Metallothionein-1K, Metallothionein-IG, MT1G, MT1K, MT1M

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human colon cancer using MT1G Antibody at dilution of 1: 100