
MTR Polyclonal Antibody

(Catalog #A59958)

Background

Catalyzes the transfer of a methyl group from methyl-cobalamin to homocysteine, yielding enzyme-bound cob(I)alamin and methionine. Subsequently, remethylates the cofactor using methyltetrahydrofolate.

Description

MTR 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

Q99707

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Methionine synthase protein (923-1265AA)

Storage

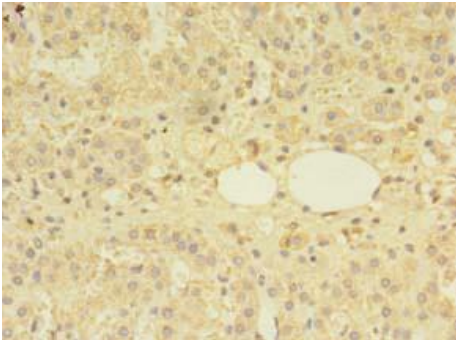
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

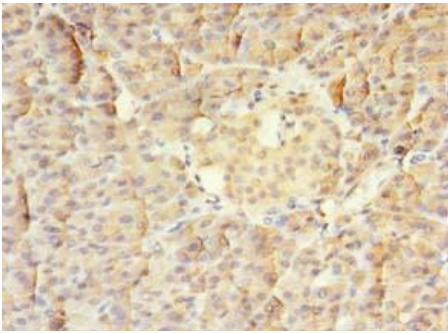
Methionine synthase (EC:2.1.1.13), MTR, 5-methyltetrahydrofolate--homocysteine methyltransferase, Vitamin-B12 dependent methionine synthase, MS

Application

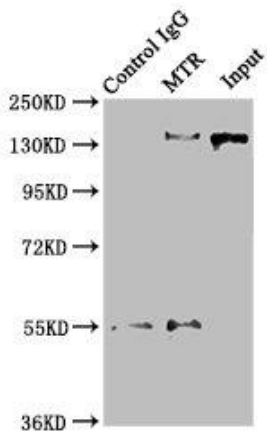
ELISA, IHC, IP; Recommended dilution: IHC 1:20 - 1:200, IP 1:200 - 1:2000



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using MTR Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human pancreatic tissue using MTR Antibody at dilution of 1:100



Immunoprecipitating MTR in HeLa whole cell lysate
Lane 1: Rabbit control IgG instead of MTR Antibody in HeLa whole cell lysate. For western blotting, a HRP-conjugated Protein G antibody was used as the secondary antibody (1/2000)
Lane 2: MTR Antibody (8ug) + HeLa whole cell lysate (500ug)
Lane 3: HeLa whole cell lysate (10ug)