

CMTR2 Polyclonal Antibody

(Catalog # A58606)

Background

S-adenosyl-L-methionine-dependent methyltransferase that mediates mRNA cap2 2'-O-ribose methylation to the 5'-cap structure of mRNAs. Methylates the ribose of the second nucleotide of a m(7)GpppG-capped mRNA and small nuclear RNA (snRNA) (cap0) to produce m(7)GpppRmpNm (cap2). Recognizes a guanosine cap on RNA independently of its N(7) methylation status. Display cap2 methylation on both cap0 and cap1. Displays a preference for cap1 RNAs.

Description

CMTR2 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

Q8IYT2

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Cap-specific mRNA (nucleoside-2'-O-)-methyltransferase 2 protein (301-600AA)

Storage

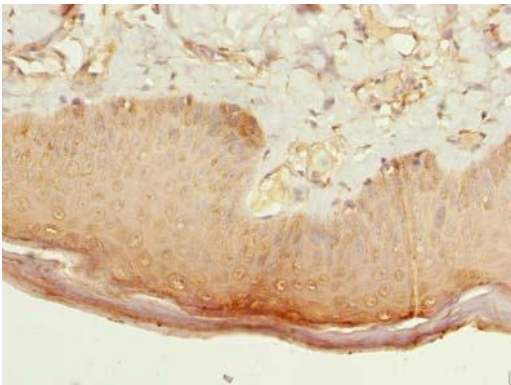
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

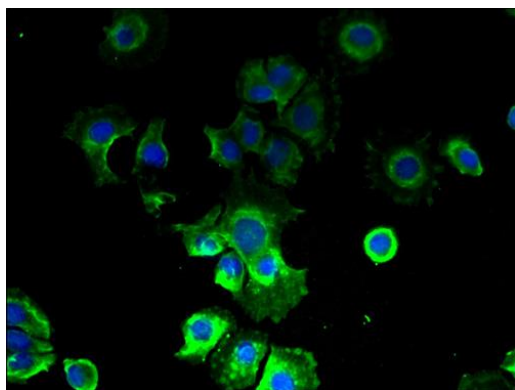
Cap-specific mRNA (nucleoside-2'-O-)-methyltransferase 2 (EC:2.1.1.296), CMTR2, AFT, FTSJD1, Cap methyltransferase 2, Cap2 2'-O-ribose methyltransferase 2, HMTr2, MTr2, FtsJ methyltransferase domain-containing protein 1, Protein adrift homolog

Application

ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human skin tissue using CMTR2 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of MCF-7 cells using CMTR2 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)