
SLC52A2 Polyclonal Antibody

(Catalog # A61018)

Background

Riboflavin transporter. Riboflavin transport is Na(+)-independent but moderately pH-sensitive. Activity is strongly inhibited by riboflavin analogs, such as lumiflavin. Weakly inhibited by flavin adenine dinucleotide (FAD) and flavin mononucleotide (FMN). In case of infection by retroviruses, acts as a cell receptor to retroviral envelopes similar to the porcine endogenous retrovirus (PERV-A).

Description

SLC52A2 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

Q9HAB3

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Solute carrier family 52, riboflavin transporter, member 2 protein (217-276AA)

Storage

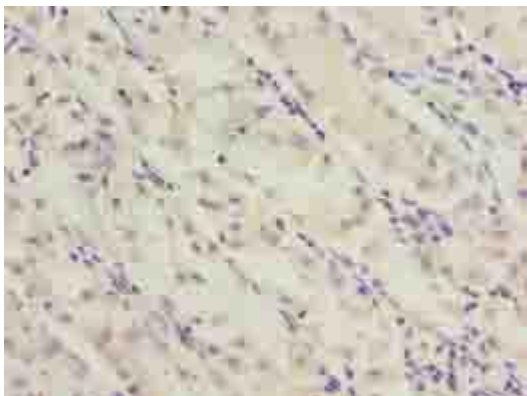
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

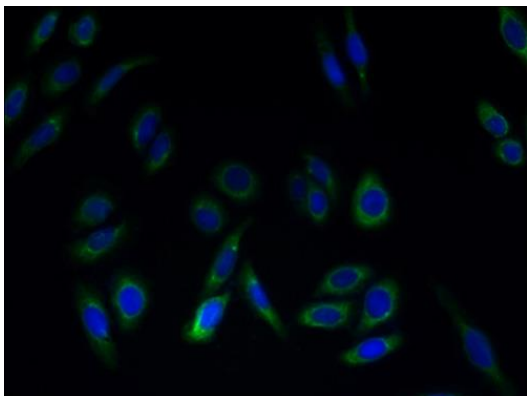
Solute carrier family 52, riboflavin transporter, member 2, SLC52A2, GPR172A, PAR1, RFT3, Porcine endogenous retrovirus A receptor 1, PERV-A receptor 1, Protein GPR172A, Riboflavin transporter 3, hRFT3

Application

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



Immunohistochemistry of paraffin-embedded human gastric cancer using SLC52A2 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using SLC52A2 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)