

SLC52A1 Polyclonal Antibody

(Catalog # A67622)

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). In case of infection by retroviruses, acts as a cell receptor to retroviral envelopes similar to the porcine endogenous retrovirus (PERV-A).

Description

SLC52A1 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

Q9NWF4

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Solute carrier family 52, riboflavin transporter, member 1 protein (212-279AA)

Storage

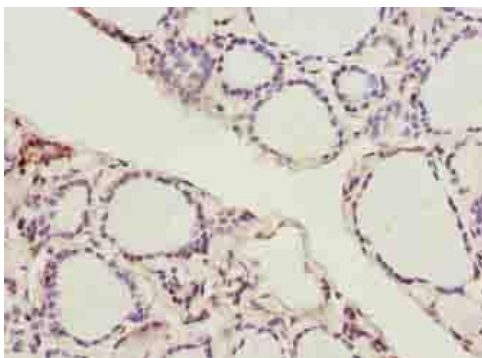
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

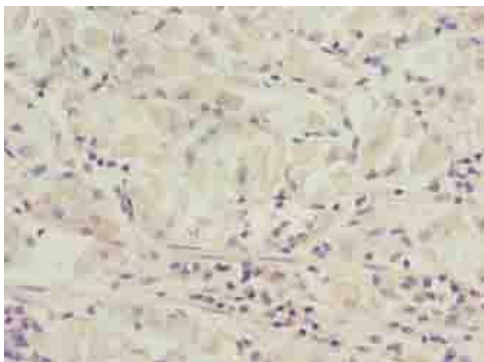
Solute carrier family 52, riboflavin transporter, member 1, Porcine endogenous retrovirus A receptor 2, PERV-A receptor 2, Protein GPR172B, Riboflavin transporter 1, hRFT1, SLC52A1, GPR172B, PAR2, RFT1

Application

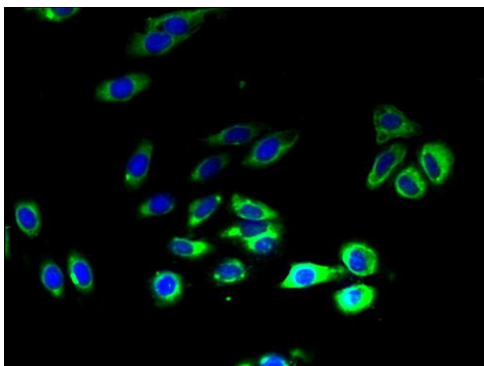
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human thyroid tissue using SLC52A1 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human gastric cancer using SLC52A1 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using SLC52A1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)
cancer using NINJ1 Antibody at dilution of 1:100