
SLC38A2 Polyclonal Antibody

(Catalog #A60994)

Background

Functions as a sodium-dependent amino acid transporter. Mediates the saturable, pH-sensitive and electrogenic cotransport of neutral amino acids and sodium ions with a stoichiometry of 1:1. May function in the transport of amino acids at the blood-brain barrier and in the supply of maternal nutrients to the fetus through the placenta.

Description

SLC38A2 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

Q96QD8

Purification

Protein G purified

Immunogen

Recombinant Human Sodium-coupled neutral amino acid transporter 2 protein (1-76AA)

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

Sodium-coupled neutral amino acid transporter 2, SLC38A2, ATA2, KIAA1382, SAT2, SNAT2, Amino acid transporter A2, Protein 40-9-1, Solute carrier family 38 member 2, System A amino acid transporter 2, System A transporter 1, System N amino acid transporter 2

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500 - 1:2000, IHC:1:20 - 1:200, IF:1:50 - 1:200



Western blot

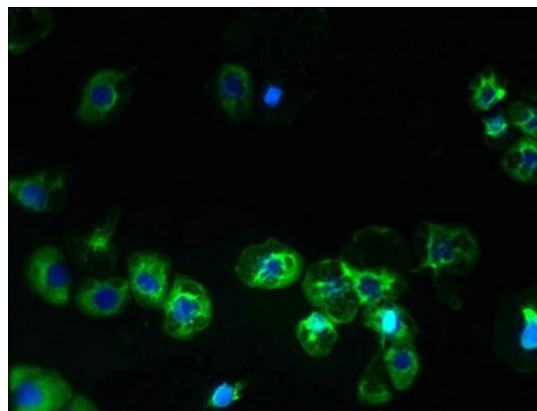
All lanes: SLC38A2 antibody at 1.5ug/ml + A431 whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

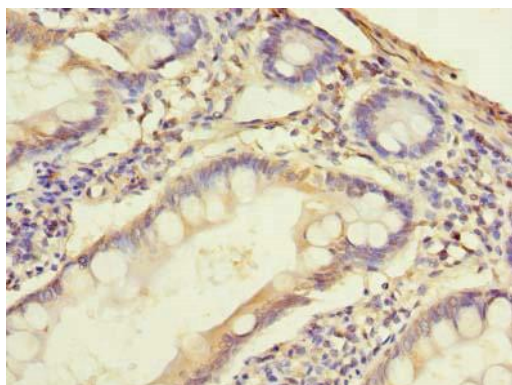
Predicted band size: 57, 46 kDa

Observed band size: 57 kDa

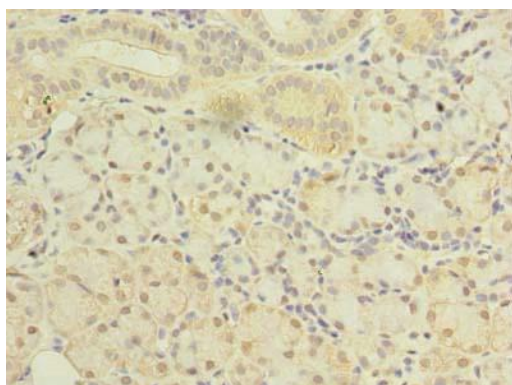


Immunofluorescent analysis of A431 cells using

SLC38A2 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human small intestine tissue using SLC38A2 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human pancreatic tissue using SLC38A2 Antibody at dilution of 1:100