

SLC7A7 Polyclonal Antibody

(Catalog # A68118)

Background

Involved in the sodium-independent uptake of dibasic amino acids and sodium-dependent uptake of some neutral amino acids. Requires coexpression with SLC3A2/4F2hc to mediate the uptake of arginine, leucine and glutamine. Plays a role in nitric oxide synthesis in human umbilical vein endothelial cells (HUVECs) via transport of L-arginine. Involved in the transport of L-arginine in monocytes.

Description

SLC7A7 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9UM01

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Y+L amino acid transporter 1 protein (325-382AA)

Storage

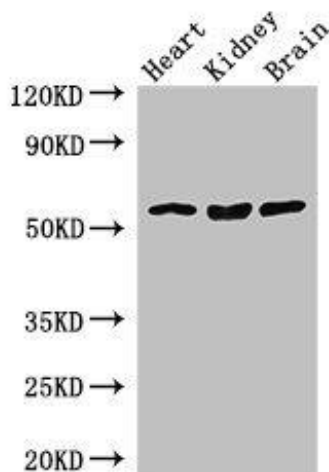
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Y+L amino acid transporter 1, Monocyte amino acid permease 2, MOP-2, Solute carrier family 7 member 7, y(+)L-type amino acid transporter 1, Y+LAT1, y+LAT-1, SLC7A7

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

Positive WB detected in: Mouse heart tissue, Mouse kidney tissue, Mouse brain tissue

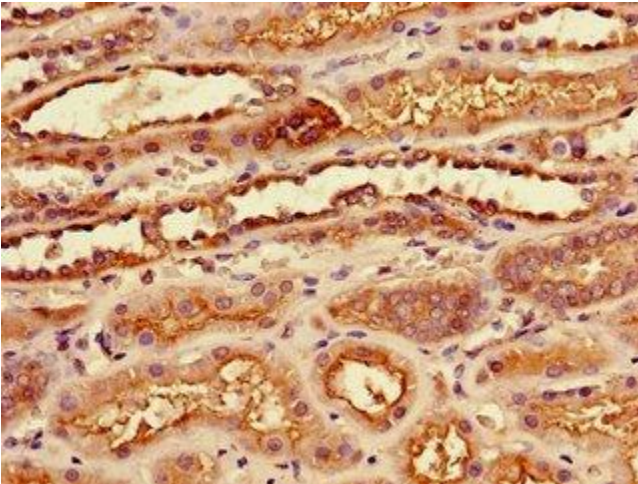
All lanes: SLC7A7 Polyclonal Antibody at 3.5ug/ml

Secondary

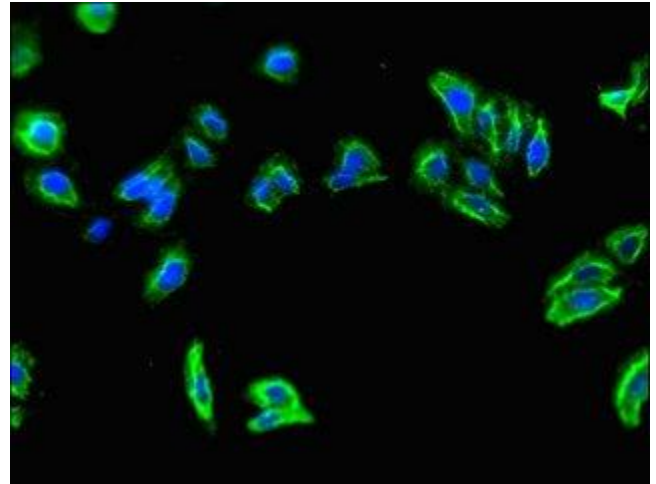
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 56 kDa

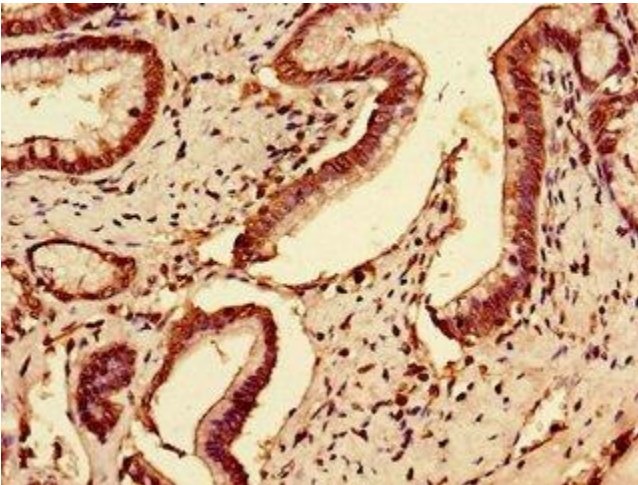
Observed band size: 56 kDa



Immunohistochemistry of paraffin-embedded human kidney tissue using SLC7A7 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using SLC7A7 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)



Immunohistochemistry of paraffin-embedded human pancreatic tissue using SLC7A7 Polyclonal Antibody at dilution of 1:100