

## PQLC2 Polyclonal Antibody

(Catalog #A60398)

### Background

Amino acid transporter that specifically mediates the pH-dependent export of the cationic amino acids arginine, histidine and lysine from lysosomes.

### Description

PQLC2 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

Q6ZP29

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Lysosomal amino acid transporter 1 homolog protein (91-150AA)

### Storage

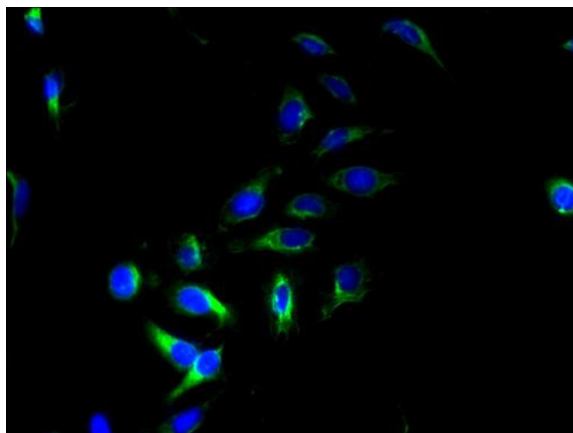
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Lysosomal amino acid transporter 1 homolog, PQLC2, PQ-loop repeat-containing protein 2

### Application

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



Immunofluorescent analysis of HeLa cells using PQLC2 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)