

## SLC35B4 Polyclonal Antibody

(Catalog # A60978)

### Background

Sugar transporter that specifically mediates the transport of UDP-xylose (UDP-Xyl) and UDP-N-acetylglucosamine (UDP-GlcNAc) from cytosol into Golgi.

### Description

SLC35B4 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

Q969S0

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human UDP-xylose and UDP-N-acetylglucosamine transporter protein (174-200AA)

### Storage

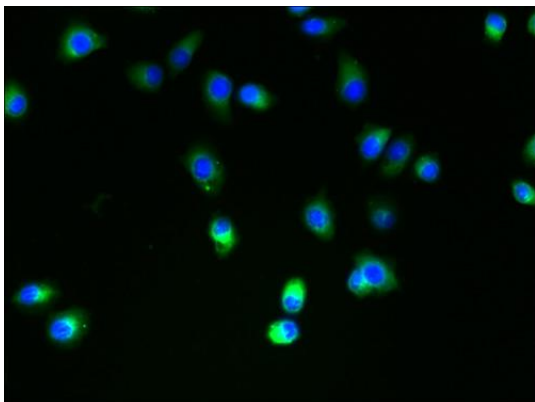
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

UDP-xylose and UDP-N-acetylglucosamine transporter, SLC35B4, YEA4, PSEC0055, Solute carrier family 35 member B4, YEA4 homolog

### Application

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



Immunofluorescent analysis of A2780 cells using SLC35B4 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)