

SLC25A32 Polyclonal Antibody

(Catalog # A67486)

Background

Transports folate across the inner membranes of mitochondria.

Description

SLC25A32 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

Q9H2D1

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Mitochondrial folate transporter/carrier protein (44-88AA)

Storage

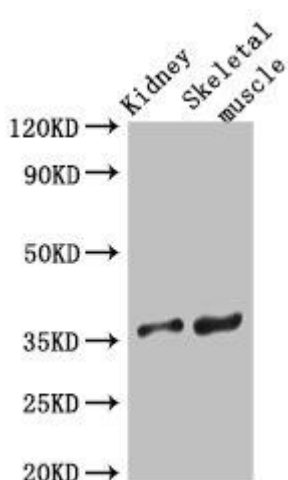
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Mitochondrial folate transporter/carrier, Solute carrier family 25 member 32, SLC25A32, MFT, MFTC

Application

ELISA, WB; Recommended dilution: WB:1:500-1:5000



Western Blot

Positive WB detected in: Mouse kidney tissue, Mouse skeletal muscle tissue

All lanes: SLC25A32 Polyclonal Antibody at 3ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 36 kDa

Observed band size: 36 kDa