

GLS Polyclonal Antibody

(Catalog # A62646)

Background

Catalyzes the first reaction in the primary pathway for the renal catabolism of glutamine. Plays a role in maintaining acid-base homeostasis. Regulates the levels of the neurotransmitter glutamate in the brain. Isoform 2 lacks catalytic activity.

Description

GLS 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

O94925

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Glutaminase kidney isoform, mitochondrial protein (616-669AA)

Storage

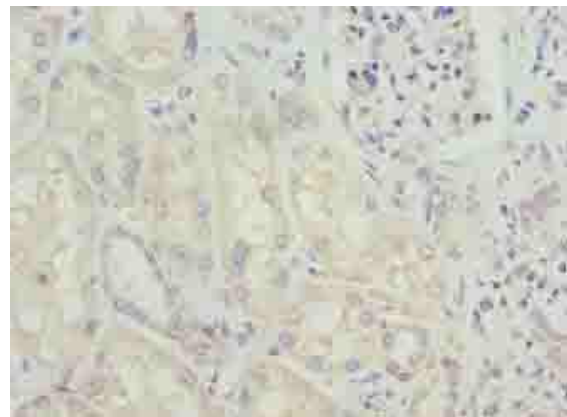
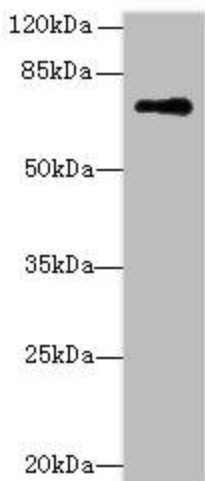
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Glutaminase kidney isoform, mitochondrial, GLS, K-glutaminase, L-glutamine amidohydrolase, GLS, GLS1, KIAA0838

Application

ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human kidney tissue at dilution of 1:100

Western blot all lanes: GLS Polyclonal Antibody at 5ug/ml+Mouse brain tissue
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 74, 18, 66 kDa
Observed band size: 73 kDa