

GALT Polyclonal Antibody

(Catalog #A54772)

Background

Galactose-1-phosphate uridylyl transferase (GALT) catalyzes the second step of the Leloir pathway of galactose metabolism, namely the conversion of UDP-glucose + galactose-1-phosphate to glucose-1-phosphate + UDP-galactose. The absence of this enzyme results in classic galactosemia in humans and can be fatal in the newborn period if lactose is not removed from the diet. The pathophysiology of galactosemia has not been clearly defined.

Description

GALT Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P07902

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Galactose-1-phosphate uridylyltransferase protein (1-379AA)

Storage

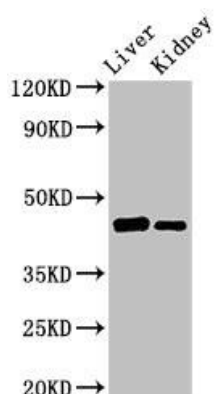
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

UDP-glucose--hexose-1-phosphate uridylyltransferase GALT

Application

ELISA, WB, IHC, IF, IP; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200, IP:1:200-1:2000



Western Blot

Positive WB detected in: Rat liver tissue, Mouse kidney tissue

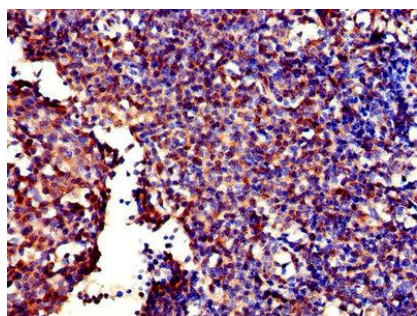
All lanes: GALT Polyclonal Antibody at 3µg/ml

Secondary

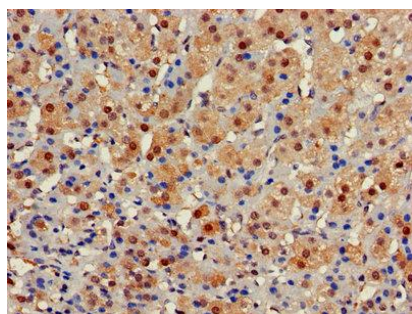
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 44, 32 kDa

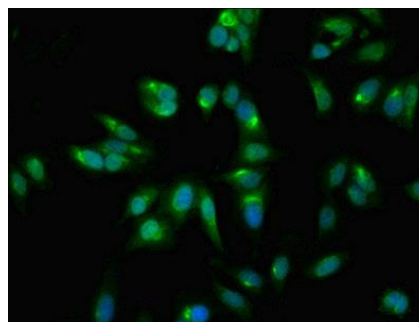
Observed band size: 44 kDa



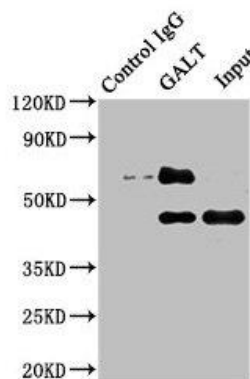
Immunohistochemistry of paraffin-embedded human tonsil tissue using GALT Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using GALT Polyclonal Antibody at dilution of 1:100



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



Immunoprecipitating GALT in HepG2 whole cell lysate

Lane 1: Rabbit control IgG instead of GALT Polyclonal Antibody in HepG2 whole cell lysate. For western blotting, a HRP-conjugated Protein G antibody was used as the secondary antibody (1/2000)

Lane 2: GALT Polyclonal Antibody (8µg) + HepG2 whole cell lysate (500µg)

Lane 3: HepG2 whole cell lysate (10µg)