

GALNT15 Polyclonal Antibody

(Catalog #A66358)

Background

Catalyzes the initial reaction in O-linked oligosaccharide biosynthesis, the transfer of an N-acetyl-D-galactosamine residue to a serine or threonine residue on the protein receptor. Although it displays a much weaker activity toward all substrates tested compared to GALNT2, it is able to transfer up to seven GalNAc residues to the Muc5AC peptide, suggesting that it can fill vicinal Thr/Ser residues in cooperation with other GALNT proteins. Prefers Muc1a as substrate.

Description

GALNT15 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

Q8N3T1

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Polypeptide N-acetylgalactosaminyltransferase 15 protein (35-335AA)

Storage

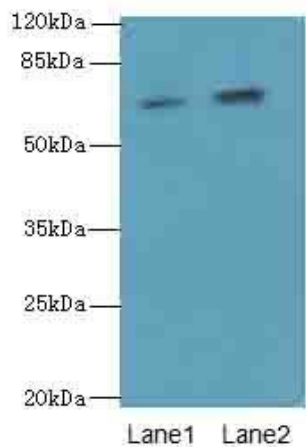
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Polypeptide N-acetylgalactosaminyltransferase 15, Polypeptide GalNAc transferase-like protein 2, GalNAc-T-like protein 2, pp-GaNTase-like protein 2, Polypeptide N-acetylgalactosaminyltransferase-like protein 2, Protein-UDP acetylgalactosaminyltransferase-like protein 2, polypeptide N-acetylgalactosaminyltransferase-like protein 2, GALNT15, GALNTL2, UNQ770/PRO1564

Application

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



Western blot

All Lanes: GALNT15 Polyclonal Antibody at 5ug/ml

Lane 1: Mouse kidney tissue

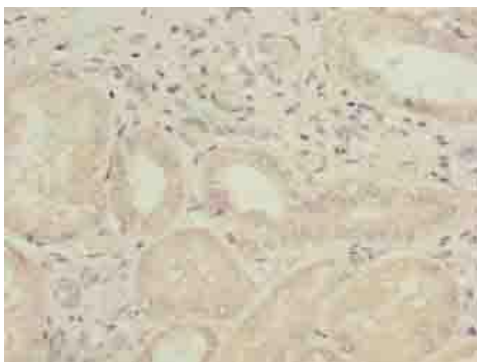
Lane 2: Mouse liver tissue

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 73 kDa

Observed band size: 73 kDa



Immunohistochemistry of paraffin-embedded human kidney tissue using GALNT15 Polyclonal Antibody at dilution of 1:100