

ST6GALNAC6 Polyclonal Antibody

(Catalog # A61162)

Background

Alpha-2,6-sialyltransferase involved in the synthesis of alpha-series gangliosides. Has activity toward GD1a, GT1b and GM1b. Has no activity toward glycoproteins. Responsible for the biosynthesis of DSGG (disialylgalactosylgloboside) from MSGG (monosialylgalactosylgloboside) in normal and malignant kidney. Participates in the synthesis of disialyl Lewis a (Le(a)).

Description

ST6GALNAC6 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

Q969X2

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Alpha-N-acetylgalactosaminide alpha-2,6-sialyltransferase 6 protein (65-333AA)

Storage

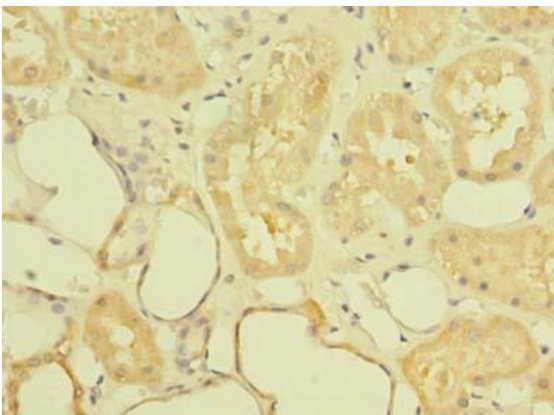
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Alpha-N-acetylgalactosaminide alpha-2, 6-sialyltransferase 6 (EC:2.4.99.-), ST6GALNAC6, SIAT7F, UNQ708/PRO1359, GalNAc alpha-2, 6-sialyltransferase VI, ST6GalNAc VI, ST6GalNAcVI, hST6GalNAc VI, Sialyltransferase 7F, SIAT7-F

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human kidney tissue using ST6GALNAC6 Polyclonal Antibody at dilution of 1:100