
ST6GALNAC3 Polyclonal Antibody

(Catalog # A66514)

Background

Involved in the biosynthesis of ganglioside GD1A from GM1B. Transfers CMP-NeuAc with an alpha-2,6-linkage to GalNAc residue on NeuAc-alpha-2,3-Gal-beta-1,3-GalNAc of glycoproteins and glycolipids. ST6GalNAcIII prefers glycolipids to glycoproteins.

Description

ST6GALNAC3 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

Q8NDV1

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Alpha-N-acetylgalactosaminide alpha-2,6-sialyltransferase 3 protein (40-305AA)

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 3, GalNAc alpha-2, 6-sialyltransferase III, ST6GalNAc III, ST6GalNAcIII, STY, Sialyltransferase 7C, SIAT7-C, ST6GALNAC3, SIAT7C, UNQ2787/PRO7177

Application

ELISA
