
ST8SIA4 Polyclonal Antibody

(Catalog # A66266)

Background

Catalyzes the polycondensation of alpha-2,8-linked sialic acid required for the synthesis of polysialic acid (PSA), which is present on the embryonic neural cell adhesion molecule (N-CAM), necessary for plasticity of neural cells.

Description

ST8SIA4 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

Q92187

Purification

>95%, Protein G purified

Immunogen

Recombinant Human CMP-N-acetylneuraminate-poly-alpha-2,8-sialyltransferase protein (21-168AA)

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

CMP-N-acetylneuraminate-poly-alpha-2, 8-sialyltransferase, Alpha-2, 8-sialyltransferase 8D, Polysialyltransferase-1, Sialyltransferase 8D, SIAT8-D, Sialyltransferase St8Sia IV, ST8SiaIV, ST8SIA4, PST, PST1, SIAT8D

Application

ELISA
