

Sialidase-3 Polyclonal Antibody

(Catalog # A57815)

Background

Plays a role in modulating the ganglioside content of the lipid bilayer at the level of membrane-bound sialyl glycoconjugates.

Description

Sialidase-3 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

Q9UQ49

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Sialidase-3 protein (2-428AA)

Storage

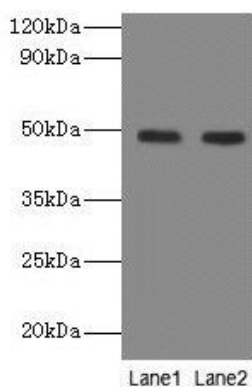
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Ganglioside sialidase, Membrane sialidase, N-acetyl-alpha-neuraminidase 3, NEU3

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: Sialidase-3 Polyclonal Antibody at 2ug/ml

Lane 1: A549 whole cell lysate

Lane 2: U251 whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 49, 52 kDa

Observed band size: 49 kDa