
ASGR1 Polyclonal Antibody, Biotin Conjugated

(Catalog # A52385)

Background

Mediates the endocytosis of plasma glycoproteins to which the terminal sialic acid residue on their complex carbohydrate moieties has been removed. The receptor recognizes terminal galactose and N-acetylgalactosamine units. After ligand binding to the receptor, the resulting complex is internalized and transported to a sorting organelle, where receptor and ligand are disassociated. The receptor then returns to the cell membrane surface.

Description

ASGR1 Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P07306

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Asialoglycoprotein receptor 1 protein (68-291AA)

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

C-type lectin domain family 4 member H1, Hepatic lectin H1, ASGR1, CLEC4H1

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
