

## A4GNT Polyclonal Antibody

(Catalog # A72325)

---

### Background

This gene encodes a protein from the glycosyltransferase 32 family. The enzyme catalyzes the transfer of N-acetylglucosamine (GlcNAc) to core 2 branched O-glycans. It forms a unique glycan, GlcNAc $\alpha$ 1 $\rightarrow$ 4Gal $\beta$ 1 $\rightarrow$ 6R and is largely associated with the Golgi apparatus membrane.

### Description

A4GNT Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q9UNA3

### Purification

Affinity Purification

### Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-340 of human A4GNT (NP\_057245.1).

### Storage

Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

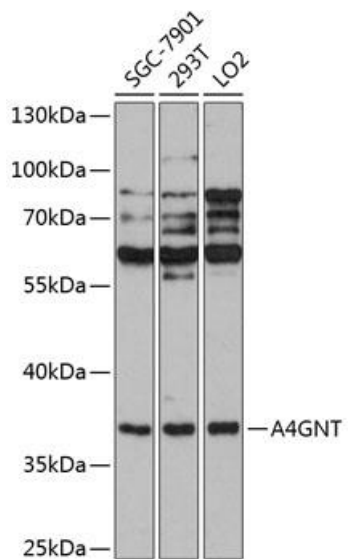
### Alternative Names

A4GNT; alpha4GnT; alpha-1,4-N-acetylglucosaminyltransferase

### Application

WB; Recommended dilution: WB 1:500 - 1:2000

---



Western blot analysis of extracts of various cell lines, using A4GNT Polyclonal Antibody at 1:1000 dilution.  
Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.  
Lysates/proteins: 25ug per lane.  
Blocking buffer: 3% nonfat dry milk in TBST.  
Exposure time: 90s.