

---

## **GALNT16 Polyclonal Antibody**

(Catalog # A66946)

---

### **Background**

Catalyzes the initial reaction in O-linked oligosaccharide biosynthesis, the transfer of an N-acetyl-D-galactosamine residue to a serine or threonine residue on the protein receptor.

### **Description**

GALNT16 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### **Formulation**

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

### **Specificity**

Human, Mouse

### **Isotype**

IgG

### **Uniprot ID**

Q8N428

### **Purification**

>95%, Protein G purified

### **Immunogen**

Recombinant Human Polypeptide N-acetylgalactosaminyltransferase 16 protein (301-558AA)

### **Storage**

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

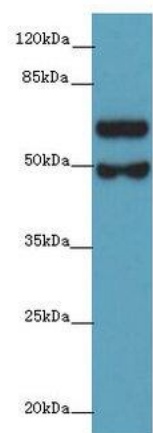
### **Alternative Names**

Polypeptide N-acetylgalactosaminyltransferase 16, Polypeptide GalNAc transferase 16, GalNAc-T16, Polypeptide GalNAc transferase-like protein 1, GalNAc-T-like protein 1, pp-GaNTase-like protein 1, Polypeptide N-acetylgalactosaminyltransferase-like protein 1, Protein-UDP acetylgalactosaminyltransferase-like protein 1, polypeptide N-acetylgalactosaminyltransferase-like protein 1, GALNT16, GALNTL1, KIAA1130

### **Application**

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

---



Western blot

All lanes: GALNT16 Polyclonal Antibody at 12 ug/ml+Mouse kidney tissue

Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 63 kDa

Observed band size: 63,49 kDa