

---

## PIGL Polyclonal Antibody

(Catalog #A68218)

---

### Background

Involved in the second step of GPI biosynthesis. De-N-acetylation of N-acetylglucosaminyl-phosphatidylinositol.

### Description

PIGL 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

Q9Y2B2

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human N-acetylglucosaminyl-phosphatidylinositol de-N-acetylase protein (23-252AA)

### Storage

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

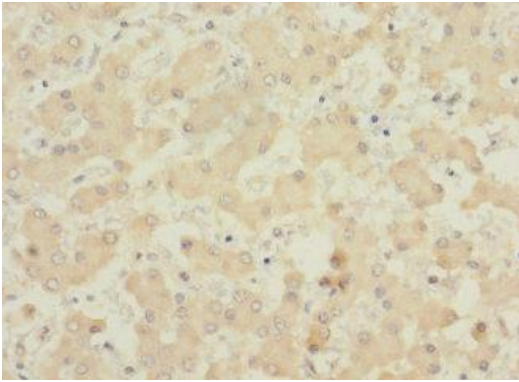
### Alternative Names

N-acetylglucosaminyl-phosphatidylinositol de-N-acetylase, Phosphatidylinositol-glycan biosynthesis class L protein, PIG-L, PIGL

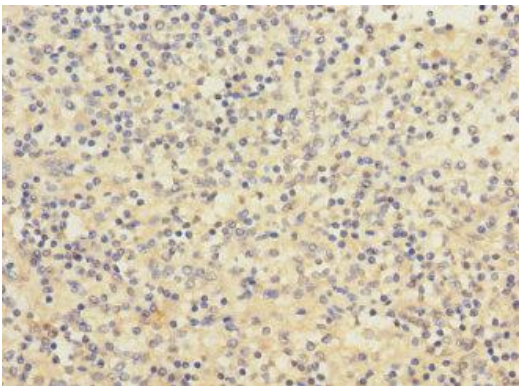
### Application

ELISA, IHC; Recommended dilution: IHC 1:20 - 1:200

---



Immunohistochemistry of paraffin-embedded human liver tissue using PIGL Polyclonal Antibody at dilution 1:100



Immunohistochemistry of paraffin-embedded human spleen tissue using PIGL Polyclonal Antibody at dilution 1:100