

## GPC Polyclonal Antibody

(Catalog # A64474)

---

### Background

Stable signal peptide (SSP) is cleaved but is apparently retained as the third component of the GP complex. The SSP is required for efficient glycoprotein expression, post-translational cleavage of GP1 and GP2, glycoprotein transport to the cell plasma membrane, formation of infectious virus particles, and acid pH-dependent glycoprotein-mediated cell fusion.

### Description

GPC Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Lymphocytic choriomeningitis virus

### Isotype

IgG

### Uniprot ID

P09991

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Lymphocytic choriomeningitis virus Pre-glycoprotein polyprotein GP complex protein (266-498AA)

### 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

Pre-glycoprotein polyprotein GP complex  
Cleaved into the following 3 chains: Stable signal peptide, SSP  
Glycoprotein G1, GP1  
Glycoprotein G2, GP2, GPC, GP-C

### Application

ELISA

---