

## **ply Polyclonal Antibody**

(Catalog # A50732)

---

### **Background**

Sulfhydryl-activated toxin that causes cytolysis by forming pores in cholesterol containing host membranes. After binding to target membranes, the protein undergoes a major conformation change, leading to its insertion in the host membrane and formation of an oligomeric pore complex. Cholesterol may be required for binding to host membranes, membrane insertion and pore formation. Can be reversibly inactivated by oxidation.

### **Description**

ply Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### **Formulation**

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

### **Specificity**

Streptococcus pneumoniae serotype 4

### **Isotype**

IgG

### **Uniprot ID**

P0C2J9

### **Purification**

>95%, Protein G purified

### **Immunogen**

Recombinant Streptococcus pneumoniae serotype 4 Pneumolysin protein (2-471AA)

### **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**

Thiol-activated cytolysin, SP\_1923, ply

### **Application**

ELISA

---