
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; Recommended dilution: ELISA: 1:1000
