

---

## ompA Polyclonal Antibody

(Catalog # A56362)

---

### Background

In elementary bodies (EBs, the infectious stage, which is able to survive outside the host cell) provides the structural integrity of the outer envelope through disulfide cross-links with the small cysteine-rich protein and the large cysteine-rich periplasmic protein. It has been described in publications as the Sarkosyl-insoluble COMC (Chlamydia outer membrane complex), and serves as the functional equivalent of peptidoglycan. By similarity. Permits diffusion of specific solutes through the outer membrane.

### Description

ompA Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Chlamydophila pneumoniae

### Isotype

IgG

### Uniprot ID

P27455

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Chlamydophila pneumoniae Major outer membrane porin protein (24-389AA)

### 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

ompA, omp1, CPn\_0695, CP\_0051, CpB0722

### Application

ELISA, Recommended dilution: ELISA 1:1000

---