

---

## ompC Polyclonal Antibody

(Catalog # A64454)

---

### Background

Forms pores that allow passive diffusion of small molecules across the outer membrane.

### Description

ompC Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Escherichia coli

### Isotype

IgG

### Uniprot ID

P06996

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Escherichia coli Outer membrane protein C protein (22-367AA)

### 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

Outer membrane protein C, Outer membrane protein 1B, Porin OmpC, ompC, meoA, par, b2215, JW2203

### Application

ELISA; Recommended dilution: ELISA 1:500-1:2000

---