
ompC Polyclonal Antibody, FITC Conjugated

(Catalog #A64456)

Background

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

Description

ompC Polyclonal Antibody, FITC Conjugated. FITC. 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 1:100-1:500
