
ompA Polyclonal Antibody

(Catalog # A55519)

Background

Required for the action of colicins K and L and for the stabilization of mating aggregates in conjugation. Serves as a receptor for a number of T-even like phages. Also acts as a porin with low permeability that allows slow penetration of small solutes.

Description

ompA Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Salmonella typhi

Isotype

IgG

Uniprot ID

Q8Z7S0

Purification

>95%, Protein G purified

Immunogen

Recombinant Salmonella typhi Outer membrane protein A protein (27-349AA)

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, STY1091, t1850

Application

ELISA
