

surA Polyclonal Antibody

(Catalog # A57221)

Background

Chaperone involved in the correct folding and assembly of outer membrane proteins, such as OmpA, OmpF and LamB. Recognizes specific patterns of aromatic residues and the orientation of their side chains, which are found more frequently in integral outer membrane proteins. May act in both early periplasmic and late outer membrane-associated steps of protein maturation. Essential for the survival of E. coli in stationary phase. Required for pilus biogenesis.

Description

surA 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

P0ABZ6

Purification

>95%, Protein G purified

Immunogen

Recombinant Escherichia coli Chaperone SurA protein (21-428AA)

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

Peptidyl-prolyl cis-trans isomerase SurA, PPIase SurA, Rotamase SurA, Survival protein A, b0053, JW0052, surA

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Lane 1: 293T whole cell lysate
Lane 2: DH5α whole cell lysate
Secondary
Goat polyclonal to Rabbit IgG at 1/10000 dilution
Predicted band size: 47 kDa
Observed band size: 47 kDa

Western blot

All Lanes: surA Polyclonal Antibody at 12ug/ml