
IpxC Polyclonal Antibody

(Catalog #A53321)

Background

The key enzyme in the biosynthesis of lipid A, a phosphorylated glycolipid that anchors the lipopolysaccharide to the outer membrane of the cell. Degraded by FtsH; when the activity of FtsH is reduced, too much lipid A and not enough phospholipids are made (both pathways use the same precursor), which is lethal.

Description

IpxC 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

P0A725

Purification

>95%, Protein G purified

Immunogen

Recombinant Escherichia coli UDP-3-O-[3-hydroxymyristoyl] N-acetylglucosamine deacetylase protein (1-305AA)

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

Protein EnvA, UDP-3-O-acyl-GlcNAc deacetylase, IpxC, asmB, envA, b0096, JW0094

Application

ELISA; Recommended dilution: ELISA 1:1000
