

lepB Polyclonal Antibody

(Catalog # A64506)

Background

integral component of membrane, integral component of plasma membrane, plasma membrane, endopeptidase activity, peptidase activity, serine-type endopeptidase activity, toxic substance binding, protein processing, proteolysis, signal peptide processing

Description

lepB 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

P00803

Purification

>95%, Protein G purified

Immunogen

Recombinant Escherichia coli Signal peptidase I protein (78-324AA)

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

Signal peptidase I, SPase I, Leader peptidase I, lepB, b2568, JW2552

Application

ELISA
