

alaC Polyclonal Antibody

(Catalog # A55431)

Background

Involved in the biosynthesis of alanine.

Description

alaC 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

P77434

Purification

>95%, Protein G purified

Immunogen

Recombinant Escherichia coli Glutamate-pyruvate aminotransferase AlaC protein (1-412AA)

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

alaC, yfdZ, b2379, JW2376

Application

ELISA
