

deoD Polyclonal Antibody

(Catalog # A53072)

Background

Cleavage of guanosine or inosine to respective bases and sugar-1-phosphate molecules.

Description

deoD 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

P0ABP8

Purification

>95%, Protein G purified

Immunogen

Recombinant Escherichia coli Purine nucleoside phosphorylase DeoD-type protein (2-239AA)

Storage

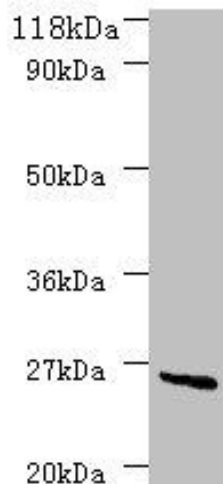
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

deoD, pup b4384, JW4347

Application

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



Western blot
All lanes: deoD Polyclonal Antibody at 2ug/ml+DH5a whole cell lysate
Secondary
Goat polyclonal to Rabbit IgG at 1/10000 dilution
Predicted band size: 26kDa
Observed band size: 26kDa