

trxA Polyclonal Antibody

(Catalog # A57734)

Background

Participates in various redox reactions through the reversible oxidation of its active center dithiol to a disulfide and catalyzes dithiol-disulfide exchange reactions.

Description

trxA 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

P0AA25

Purification

>95%, Protein G purified

Immunogen

Recombinant Escherichia coli Thioredoxin-1 protein (2-109AA)

Storage

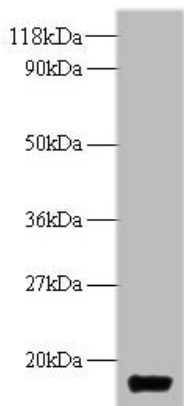
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

trxA, fipA, tsnC

Application

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



Western blot

All lanes: trxA Polyclonal Antibody at 2ug/ml+DH5a whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 12kDa

Observed band size: 12kDa