
tetR Polyclonal Antibody, FITC Conjugated

(Catalog # A64540)

Background

TetR is the repressor of the tetracycline resistance element; its N-terminal region forms a helix-turn-helix structure and binds DNA. Binding of tetracycline to TetR reduces the repressor affinity for the tetracycline resistance gene (tetA) promoter operator sites.

Description

tetR Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Escherichia coli

Isotype

IgG

Uniprot ID

P04483

Purification

>95%, Protein G purified

Immunogen

Recombinant Escherichia coli Tetracycline repressor protein class B from transposon Tn10 protein (1-207AA)

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

Tetracycline repressor protein class B from transposon Tn10, tetR

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
