
bar Polyclonal Antibody, Biotin Conjugated

(Catalog # A54540)

Background

This enzyme is an effector of phosphinothricin tripeptide (PTT or bialaphos) resistance. Inactivates PTT by transfer of an acetyl group.

Description

bar Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

Formulation

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

Specificity

Streptomyces hygrosopicus

Isotype

IgG

Uniprot ID

P16426

Purification

>95%, Protein G purified

Immunogen

Recombinant Streptomyces hygrosopicus Phosphinothricin N-acetyltransferase protein (1-183AA)

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

Phosphinothricin-resistance protein bar

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
