
scaIM Polyclonal Antibody

(Catalog # A56851)

Background

This methylase recognizes the double-stranded sequence AGTACT, causes specific methylation on C-5 on both strands, and protects the DNA from cleavage by the ScaI endonuclease.

Description

scaIM Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Streptomyces caespitosus

Isotype

IgG

Uniprot ID

O52692

Purification

>95%, Protein G purified

Immunogen

Recombinant Streptomyces caespitosus Modification methylase ScaI protein (1-304AA)

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

M.ScaI, N-4 cytosine-specific methyltransferase ScaI, scaIM

Application

ELISA
