

SAT1 Polyclonal Antibody

(Catalog # A53921)

Description

SAT1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P21673

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Diamine acetyltransferase 1 protein (5-171AA)

Storage

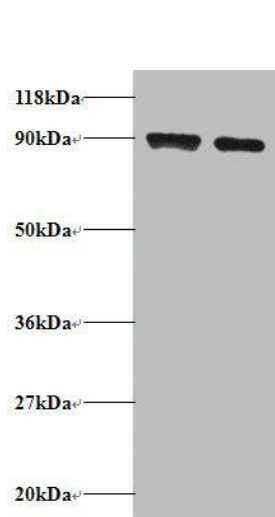
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Polyamine N-acetyltransferase 1, Putrescine acetyltransferase, Spermidine/spermine N(1)-acetyltransferase 1, SSAT, SSAT-1

Application

ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200



Western blot

All lanes: SAT1 Polyclonal Antibody at 2/ml

Lane 1: 293T whole cell lysate

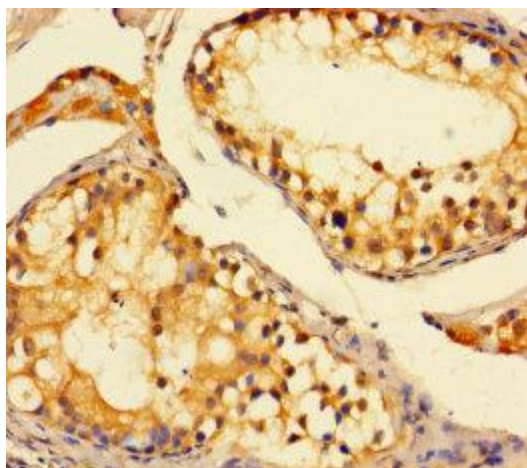
Lane 2: EC109 whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/15000 dilution

Predicted band size: 19 kDa

Observed band size: 90 kDa



Immunohistochemistry of paraffin-embedded human testis tissue at dilution of 1:100