

AZIN1 Polyclonal Antibody

(Catalog # A62142)

Background

Antizyme inhibitor protein that positively regulates ornithine decarboxylase (ODC) activity and polyamine uptake by counteracting the negative effect of antizyme OAZ1, OAZ2 and OAZ3 on ODC1 activity (PubMed:17900240). Inhibits antizyme-dependent ODC degradation by binding to antizymes. Releases ODC1 from its inactive complex with antizymes, leading to formation of the catalytically active ODC1.

Description

AZIN1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

O14977

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Antizyme inhibitor 1 protein (101-400AA)

Storage

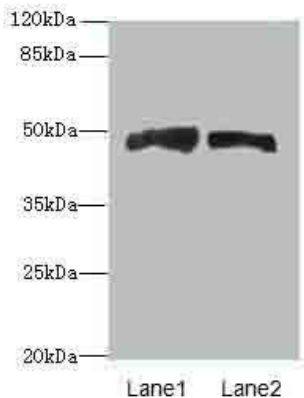
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Antizyme inhibitor 1, AZI, Ornithine decarboxylase antizyme inhibitor, AZIN1, OAZI, OAZIN

Application

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



Western blot

All lanes: AZIN1 Polyclonal Antibody at 4ug/ml

Lane 1: Mouse liver tissue

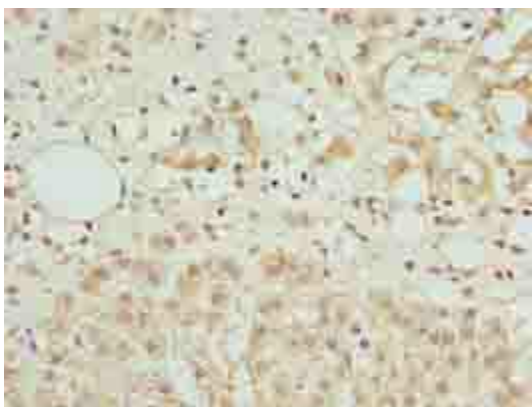
Lane 2: PC-3 whole cell lysate

Secondary

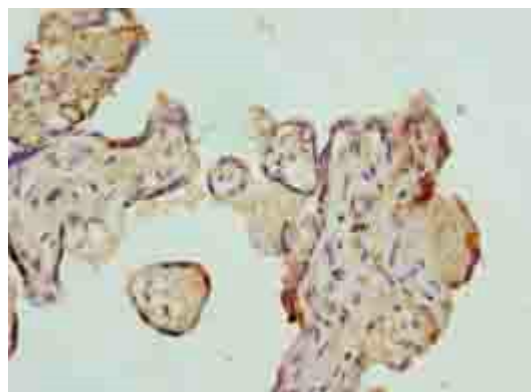
Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 50 kDa

Observed band size: 50 kDa



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



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