

NAALADL2 Polyclonal Antibody

(Catalog # A56611)

Background

May be catalytically inactive.

Description

NAALADL2 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

Q58DX5

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Inactive N-acetylated-alpha-linked acidic dipeptidase-like protein 2 protein (143-795AA)

Storage

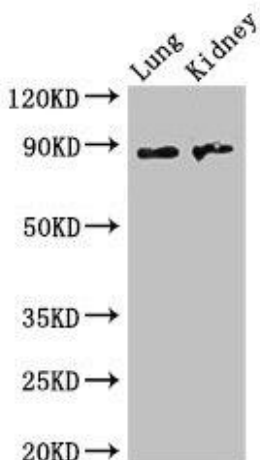
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

NAALADL2, NAALADase L2

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:2000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

Positive WB detected in: Mouse lung tissue, Mouse kidney tissue

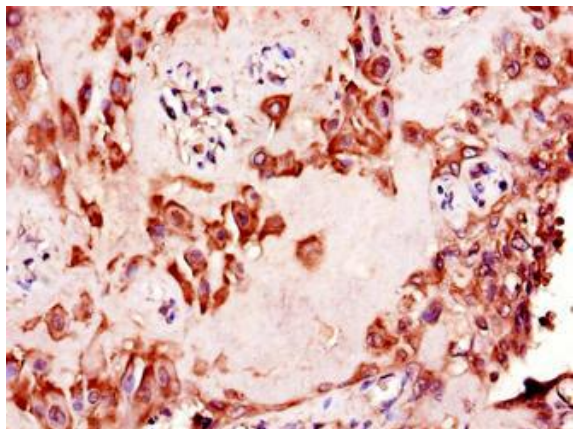
All lanes: NAALADL2 Polyclonal Antibody at 3.5µg/ml

Secondary

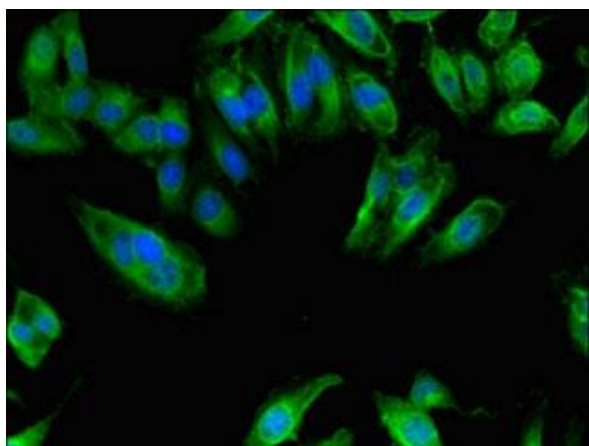
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 89, 37 kDa

Observed band size: 89 kDa



Immunohistochemistry of paraffin-embedded human placenta tissue using NAALADL2 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HeLa cells using NAALADL2 Polyclonal Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)