

CNIH4 Polyclonal Antibody

(Catalog # A58610)

Background

Involved in G protein-coupled receptors (GPCRs) trafficking from the endoplasmic reticulum to the cell surface; it promotes the exit of GPCRs from the early secretory pathway, likely through interaction with the COPII machinery (PubMed:24405750).

Description

CNIH4 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q9P003

Purification

Affinity Purified

Immunogen

Recombinant Human Protein cornichon homolog 4 protein (78-117AA)

Storage

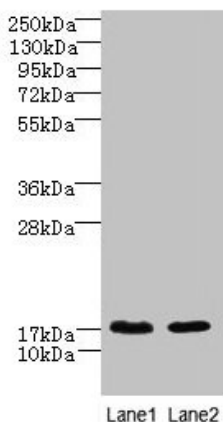
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Protein cornichon homolog 4, CNIH-4, CNIH4, HSPC163, Cornichon family AMPA receptor auxiliary protein 4

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200, IF:1:50-1:200



Western blot

All Lanes: CNIH4 Polyclonal Antibody at 0.3ug/ml

Lane 1: A431 whole cell lysate

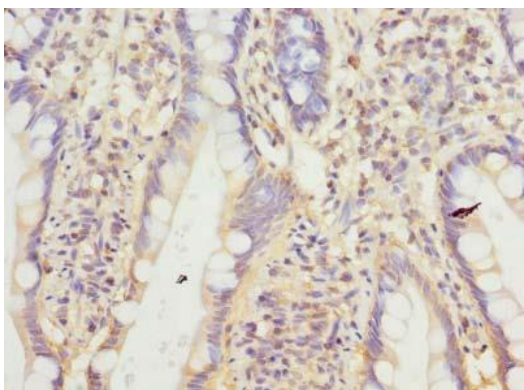
Lane 2: U251 whole cell lysate

Secondary

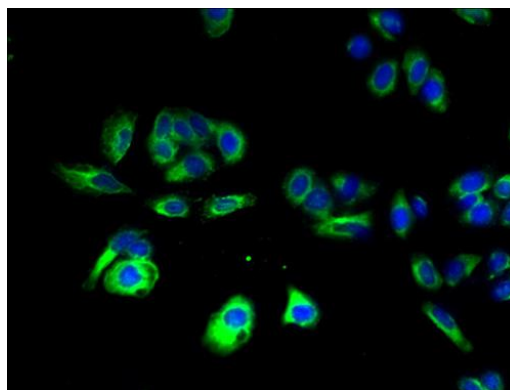
Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 16 kDa

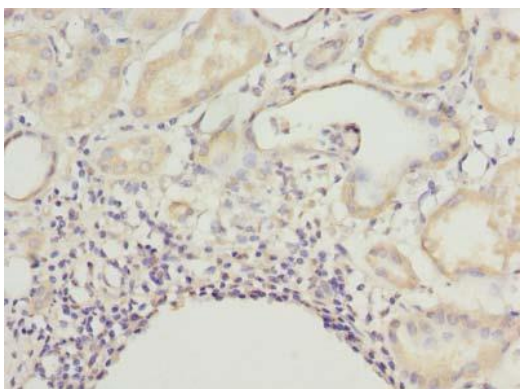
Observed band size: 16 kDa



Immunohistochemistry of paraffin-embedded human small intestine tissue using CNIH4 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using CNIH4 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human kidney tissue using CNIH4 Polyclonal Antibody at dilution of 1:100