

NINJ1 Polyclonal Antibody

(Catalog # A67086)

Background

Homophilic cell adhesion molecule that promotes axonal growth. May play a role in nerve regeneration and in the formation and function of other tissues. Cell adhesion requires divalent cations.

Description

NINJ1 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

Q92982

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Ninjurin-1 protein (1-80AA)

Storage

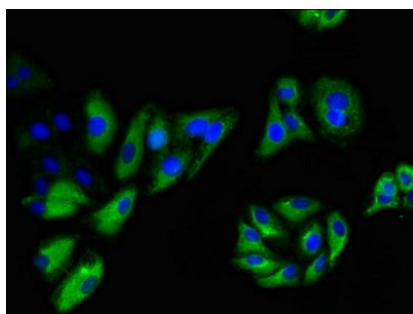
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

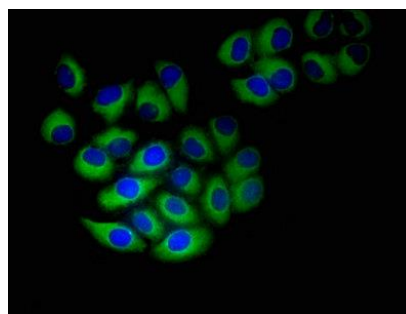
Ninjurin-1, Nerve injury-induced protein 1, NINJ1

Application

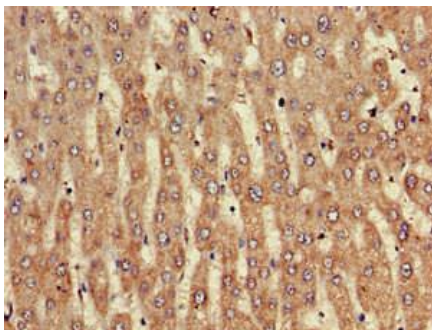
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



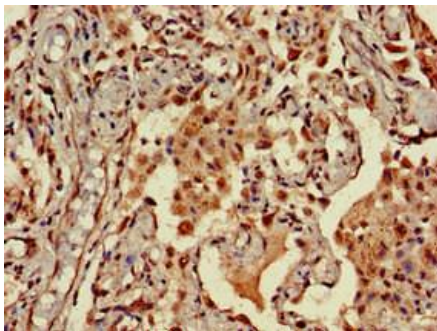
Immunofluorescent analysis of HepG2 cells using NINJ1 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunofluorescent analysis of A549 cells using NINJ1 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human liver tissue using NINJ1 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human lung cancer using NINJ1 Antibody at dilution of 1:100