
NETO1 Polyclonal Antibody

(Catalog #A70083)

Background

Involved in the development and/or maintenance of neuronal circuitry. Accessory subunit of the neuronal N-methyl-D-aspartate receptor (NMDAR) critical for maintaining the abundance of GRIN2A-containing NMDARs in the postsynaptic density. Regulates long-term NMDA receptor-dependent synaptic plasticity and cognition, at least in the context of spatial learning and memory.

Description

NETO1 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

Q8TDF5

Purification

Protein G purified

Immunogen

Recombinant Human Neuropilin and tolloid-like protein 1 protein (367-507AA)

Storage

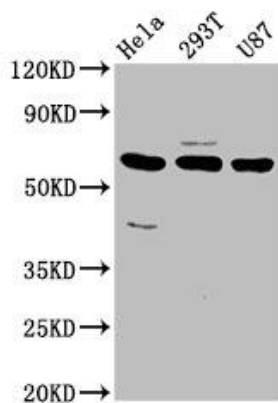
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Neuropilin and tolloid-like protein 1, Brain-specific transmembrane protein containing 2 CUB and 1 LDL-receptor class A domains protein 1, NETO1, BTCL1

Application

ELISA, WB, IF; Recommended dilution: WB 1:500 - 1:5000, IF 1:50 - 1:200



Western Blot

Positive WB detected in: HeLa whole cell lysate, 293T whole cell lysate, U87 whole cell lysate

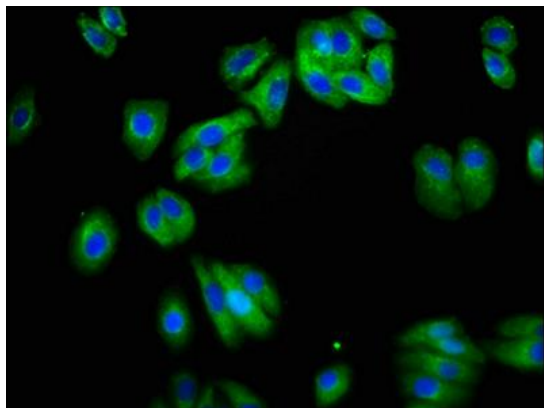
All lanes: NETO1 antibody at 3.2ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 61, 18 KDa

Observed band size: 61 KDa



Immunofluorescence staining of HepG2 cells with NETO1 Polyclonal Antibody at 1:200, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).