
NREP Polyclonal Antibody

(Catalog # A68871)

Background

May have roles in neural function. Ectopic expression augments motility of gliomas. Promotes also axonal regeneration. May also have functions in cellular differentiation. Induces differentiation of fibroblast into myofibroblast and myofibroblast amoeboid migration. Increases retinoic-acid regulation of lipid-droplet biogenesis. Down-regulates the expression of TGFB1 and TGFB2 but not of TGFB3. May play a role in the regulation of alveolar generation.

Description

NREP 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

Q16612

Purification

Protein G purified

Immunogen

Recombinant Human Neuronal regeneration-related protein (1-68AA)

Storage

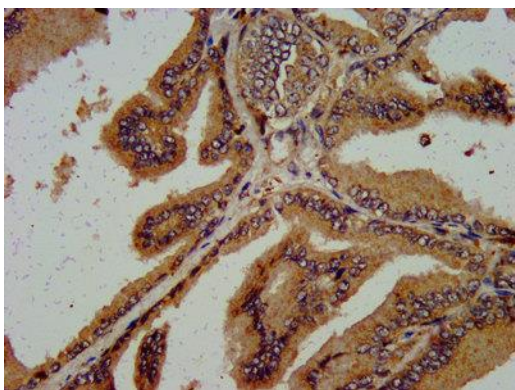
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

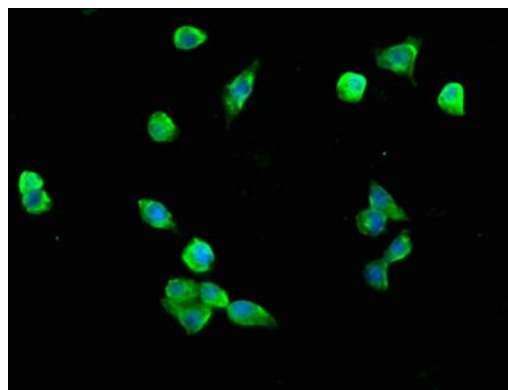
Neuronal regeneration-related protein, Neuronal protein 3.1, Protein p311, NREP, C5orf13, P311

Application

ELISA, IHC, IF; Recommended dilution: IHC:1:500-1:1000, IF:1:200-1:500



IHC image of NREP Polyclonal Antibody diluted at 1:600 and staining in paraffin-embedded human prostate tissue performed on a Leica Bond™ system.



Immunofluorescence staining of SH-SY5Y cells with NREP 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).