
NIT1 Polyclonal Antibody

(Catalog #A65766)

Background

Plays a role in cell growth and apoptosis: loss of expression promotes cell growth and resistance to DNA damage stress. Has tumor suppressor properties that enhances the apoptotic responsiveness in cancer cells; this effect is additive to the tumor suppressor activity of FHIT. It is also a negative regulator of primary T-cells. Has apparently no omega-amidase activity such as NIT2.

Description

NIT1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q86X76

Purification

Protein G purified

Immunogen

Recombinant Human Deaminated glutathione amidase protein (1-243AA)

Storage

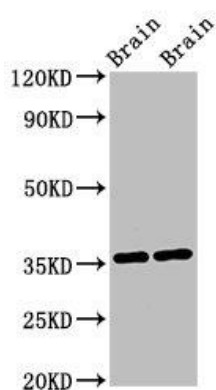
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Nitrilase homolog 1, NIT1

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:1000-1:2000, IF:1:200-1:500



Western Blot

Positive WB detected in: Mouse brain tissue, Rat brain tissue

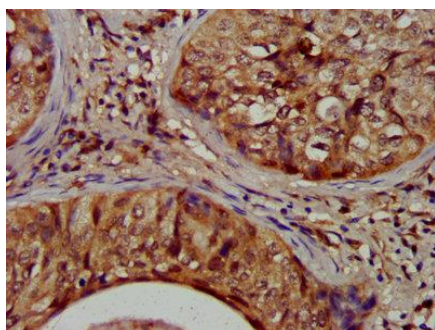
All lanes: NIT1 antibody at 3ug/ml

Secondary

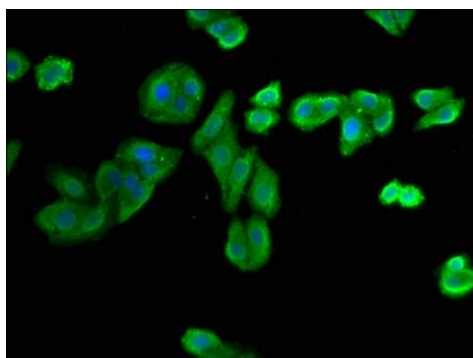
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 36, 32, 41, 35, 27 kDa

Observed band size: 36 kDa



IHC image of NIT1 Antibody diluted at 1:1200 and staining in paraffin-embedded human cervical cancer performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Immunofluorescence staining of HepG2 cells with NIT1 Antibody at 1:400, 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).