
NMNAT3 Polyclonal Antibody

(Catalog # A57005)

Background

Catalyzes the formation of NAD⁺ from nicotinamide mononucleotide (NMN) and ATP. Can also use the deamidated form; nicotinic acid mononucleotide (NaMN) as substrate with the same efficiency. Can use triazofurin monophosphate (TrMP) as substrate. Can also use GTP and ITP as nucleotide donors. Also catalyzes the reverse reaction, i. e. the pyrophosphorolytic cleavage of NAD⁺. For the pyrophosphorolytic activity, can use NAD (+), NADH, NAAD, nicotinic acid adenine dinucleotide phosphate (NHD), nicotinamide guanine dinucleotide (NGD) as substrates. Fails to cleave phosphorylated dinucleotides NADP⁺, NADPH and NAADP⁺. Protects against axonal degeneration following injury.

Description

NMNAT3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q96T66

Purification

Protein G purified

Immunogen

Recombinant Human Nicotinamide/nicotinic acid mononucleotide adenylyltransferase 3 protein (1-215AA)

Storage

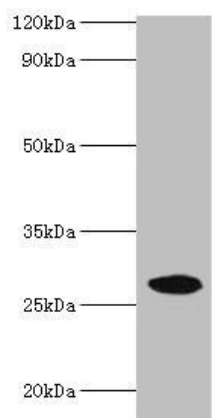
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

NMN adenylyltransferase 3, Nicotinate-nucleotide adenylyltransferase 3, NaMN adenylyltransferase 3, Pyridine nucleotide adenylyltransferase 3, PNAT-3FKSG76, NMNAT3

Application

ELISA, WB, IHC; Recommended dilution: WB:1:200-1:1000, IHC:1:20-1:200



Western blot

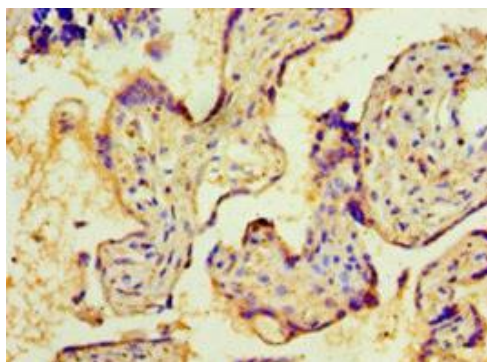
All lanes: NMNAT3 antibody at 2ug/ml + Mouse brain tissue

Secondary

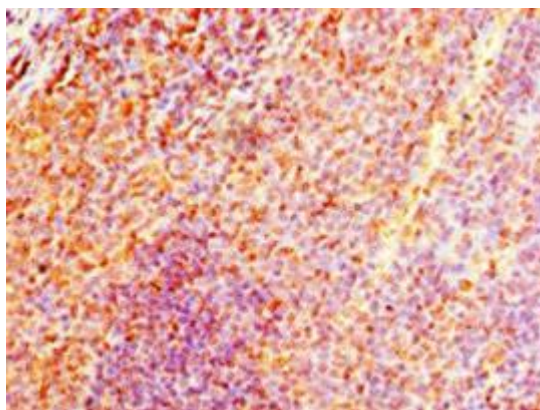
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 29, 25, 19 kDa

Observed band size: 29 kDa



Immunohistochemistry of paraffin-embedded human placenta tissue using NMNAT3 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human tonsil tissue using NMNAT3 Antibody at dilution of 1:100