

NF-1A antibody (pAb)

Catalog Nos: 39397, 39398

RRID: AB_2314931

Isotype: Serum

Application(s): ICC, IF, IHC, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 60 kDa

Background: NF-1A (NF-1, Nuclear factor-1) is a ubiquitous 47 kDa dimeric DNA-binding protein whose recognition sequence, TGG(C/A)N(5)GCCAA, is found in the genomes of a number of DNA viruses. The same sequence occurs frequently in the human genome. The NF-1 protein stimulates initiation of adenovirus DNA replication *in vitro* and is capable of stimulating the transcription of genes in cooperation with other factors, such as the estrogen receptor. The Nuclear factor-1 (NF-1) family of transcription factors are encoded by the genes NF-1A, NF-1B2, NF-1C and NF-1X and feature a highly conserved N-terminal DNA binding region and a variable C-terminal transcriptional activation or repression domain. NF-1 binding controls the expression of many target genes, including insulin, TGF-β, nucleotide, steroid hormone and vitamin synthetic genes, TNFα and FSH. NF-1 transcripts can be differentially spliced and the resultant protein variants homo- and heterodimerize, allowing for multiple potential transcriptional activation and repression activities.

Immunogen: This NF-1A antibody was raised against a synthetic peptide corresponding to amino acid residue(s) 478-492 of human NF-1A.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Validated Applications:

ICC/IF: 1:250 - 1:500 dilution

WB: 1:1,000 - 1:5,000 dilution

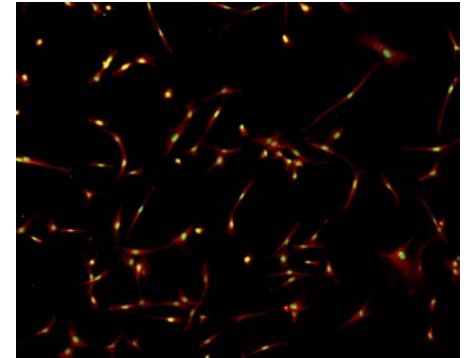
Published Applications:

IF

IHC

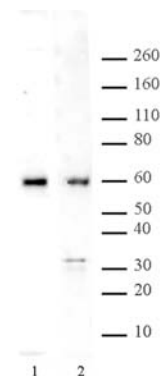
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



NF-1A antibody (pAb) tested by immunofluorescence.

Detection of NF-1A antibody by immunofluorescence. Human Astrocytes-cerebellar (HA-c) cells were stained with NF-1A antibody at a 1:500 dilution. Green: NF-1A antibody (pAb)
Red: Vimentin



NF-1A pAb tested by Western blot.

Lane 1: Nuclear extract (15 µg) of MCF-7 cells was probed with NF-1A pAb (1:1,000).
Lane 2: Nuclear extract (20 µg) of HeLa cells was probed with NF-1A pAb (1:2,500).