

NF-1A antibody (pAb)

Catalog Nos: 39397, 39398

RRID: AB_2314931

Isotype: Serum

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

Reactivity: Human

Volumes: 100 μ l, 10 μ l

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, TNF α 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:

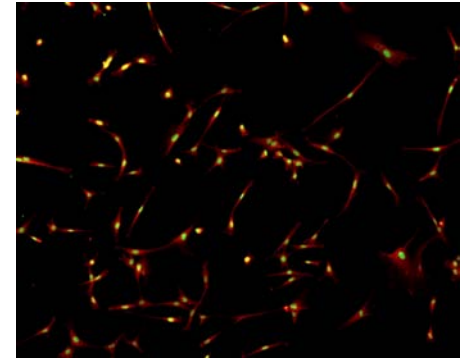
Applications Validated by Active Motif:

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

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

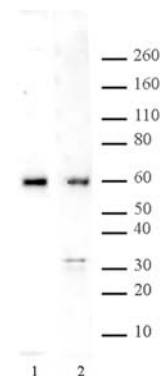
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 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 μ g) of MCF-7 cells was probed with NF-1A pAb (1:1,000).
Lane 2: Nuclear extract (20 μ g) of HeLa cells was probed with NF-1A pAb (1:2,500).