

## NF-1B2 antibody (pAb)

**Catalog No:** 39091

**RRID:** AB\_2314934

**Isotype:** IgG

**Application(s):** ChIP, EMSA, IF, WB

**Reactivity:** Human, Rat

**Volume:** 100 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 50 kDa

**Background:** NF-1B2 (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- $\beta$ , nucleotide, steroid hormone and vitamin synthetic genes, TNF $\alpha$  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-1B2 antibody was raised against a synthetic peptide corresponding to amino acid residues 402-415 of human NF-1B2.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

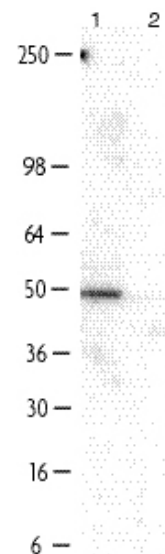
Applications Validated by Active Motif:

WB: 1:500 - 1:2,000 dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

**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-1B2 antibody (pAb) tested by Western blot.

NF-1B2 detection by Western blot analysis. NF-1B2 detected in MG-63 nuclear extract (20µg) using NF-1B2 antibody at a 1:1,000 dilution (lane 1). The signal is eliminated by adding the immunizing peptide to the reaction mixture (lane 2).