

N-Myc antibody (pAb)

Catalog Nos: 61185, 61186

RRID: AB_2793543

Isotype: IgG

Application(s): ChIP-Seq, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 60 kDa

Background: N-Myc (MYCN) is a member of the Myc family of proto-oncogenes. Myc family members play crucial roles in regulating cell proliferation, size and differentiation. N-Myc is required for proper development of the nervous system. Myc family proteins can induce malignant transformation inappropriately modulating gene transcription, leading to unchecked cell proliferation. Misexpression of N-Myc is often associated with pediatric neural cancers such as neuroblastoma, medulloblastoma and retinoblastoma.

Immunogen: This N-Myc antibody was raised against a peptide derived from the C-terminal region of human N-Myc.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

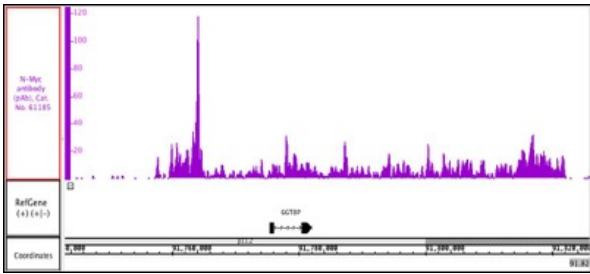
ChIP-Seq: 10 - 20 µl per ChIP

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

For optimal results in Western Blot, primary antibody incubations should be performed at 4°C. 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.



N-Myc antibody (pAb) tested by ChIP-Seq

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT® High Sensitivity Kit (Cat. No. 53040) with 200 µg of chromatin from human cell line chromatin and 20 µg N-Myc antibody. ChIP DNA was sequenced on the Illumina HiSeq and 9.2 million sequence tags were mapped to identify N-Myc binding sites on chromosome 3. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, [here](#).

N-Myc antibody (pAb) tested by Western blot.

Nuclear extract (20 µg) of Kelly cells probed with N-Myc antibody at a dilution of 1:1,000

