

NonO antibody (pAb)

Catalog Nos: 61247, 61248

RRID: AB_2793567

Isotype: IgG

Application(s): IP, WB

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Rabbit

Concentration: 1 µg/µl

Molecular Weight: 60 kDa

Background: NonO (non-POU domain containing, octamer-binding, p54(nrb), NRB54, NMT55) is a DNA- and RNA binding protein, involved in several nuclear processes. Binds the conventional octamer sequence in double-stranded DNA. Also binds single-stranded DNA and RNA at a site independent of the duplex site. This gene also encodes an RNA-binding protein which plays various roles in the nucleus, including transcriptional regulation and RNA splicing. A rearrangement between this gene and the transcription factor E3 gene has been observed in papillary renal cell carcinoma.

Immunogen: This NonO antibody was raised against a peptide corresponding to the C-terminus of human NonO.

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:

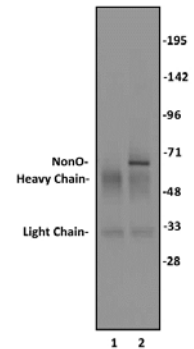
IP: 5 µg per IP

WB: 0.5 - 2 µg/ml dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

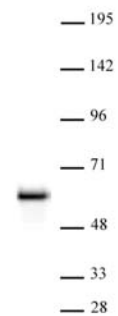
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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



NonO antibody (pAb) tested by Immunoprecipitation.

5 µg of NonO antibody was used to immunoprecipitate NonO from 400 µg of HeLa whole cell extract (lane 2). 5 µg of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the NonO antibody at a dilution of 1 µg/ml.



NonO antibody (pAb) tested by Western blot.

Whole cell extract of HeLa cells (20 µg) probed with NonO antibody (1 µg/ml dilution).