

## Recombinant NONO protein

**Catalog No:** 31539, 31939

**Lot No:** 19916001

**Expressed In:** Baculovirus

**Quantity:** 20 µg

**Concentration:** 0.4 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant full length NONO protein is supplied at a concentration of 0.4 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

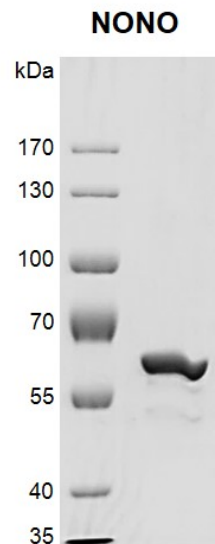
**Background:** NONO (non-POU domain containing, octamer-binding, p54(nrb), NRB54, NMT55) is a DNA- and RNA binding protein and involved in several nuclear processes. NONO binds the conventional octamer sequence in double-stranded DNA. It also binds single-stranded DNA and RNA at a site independent of 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.

**Protein Details:** Recombinant human NONO was expressed in a baculovirus expression system as the full length protein (accession number NP\_031389.3) with an N-terminal FLAG tag. The molecular weight of the protein is 57.2 kDa. The NONO protein is >90% pure by SDS-PAGE.

**Application Notes:** Recombinant NONO is suitable for use in RNA binding, assays and protein-protein interaction.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. 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.



**Recombinant NONO protein gel.**  
NONO protein was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.