

Recombinant NSUN2 protein

Catalog No: 81171, 81871

Lot No: 15818001

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.3 µg/µl

Source: Human

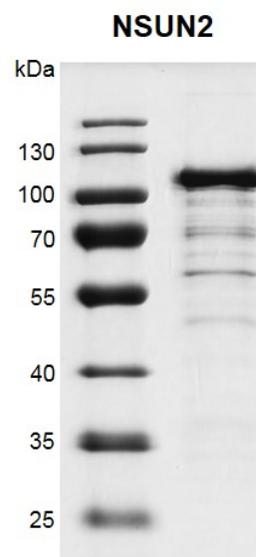
Buffer Contents: Recombinant NSUN2 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

Background: NSUN2 (NOP2/Sun Domain Family, Member 2), also called as TRM4 (TRNA Methyltransferase 4 Homolog), is a member of NSUN family. It's a RNA methyltransferase that methylates tRNAs, and possibly RNA polymerase III transcripts. NSUN2 also methylates cytosine to 5-methylcytosine (m5C) at positions 34 and 48 of introncontaining tRNA(Leu)(CAA) precursors, and at positions 48, 49 and 50 of tRNA(Gly)(GCC) precursors. It may act downstream of Myc to regulate epidermal cell growth and proliferation. NSUN2 is required for proper spindle assembly and chromosome segregation, independently of its methyltransferase activity.

Protein Details: Recombinant NSUN2 protein was expressed in a baculovirus expression system as the full length protein (accession number NP_060225.4) with an N-terminal FLAG tag. The molecular weight of the protein is 88.1 kDa.

Application Notes: Recombinant NSUN2 protein is suitable for use in enzyme kinetics, inhibitor screening, and selectivity profiling.

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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



Recombinant NSUN2 protein
10% SDS-PAGE Coomassie staining
MW: 88.1 kDa
Purity: >80%