

Recombinant NSUN6 protein

Catalog No: 81174

Lot No: 19218001

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.5 µg/µl

Source: Human

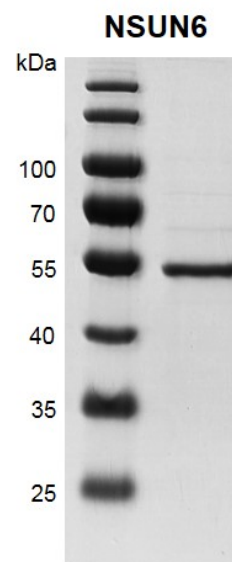
Buffer Contents: Recombinant NSUN6 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

Background: NSUN6 (NOP2/Sun Domain Family, Member 6), also called as NOPD1, is a member of NSUN family. It may have S-adenosyl-L-methionine-dependent methyltransferase activity. Recently, researches found that human NSUN6 is associated with tRNAs and acts as a tRNA methyltransferase. It may play a major role in bone metastasis through the methylation of Hippo/MST1 and consequent activation of YAP, implying that NSUN6 could be a valuable target for bone metastasis and therapy-resistant oncology research

Protein Details: Recombinant NSUN6 protein was expressed in a baculovirus expression system as the full length protein (accession number NP_872349.1) with an N-terminal FLAG tag. The molecular weight of the protein is 53 kDa.

Application Notes: Recombinant NSUN6 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 NSUN6 protein
10% SDS-PAGE Coomassie staining
MW: 53 kDa
Purity: >90%