

NSUN2

CATALOG NO.: MT-21-592

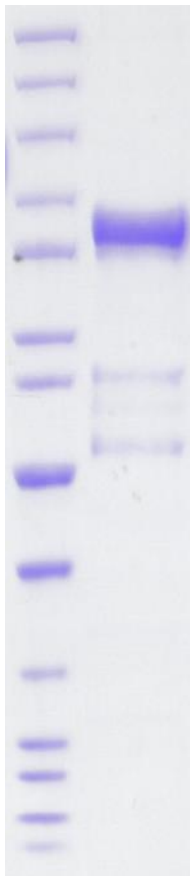
LOT:

DESCRIPTION: Human recombinant NSUN2 (residues 2-767; Genbank Accession # NP_001193455.2; MW = 87.89 kDa) expressed as a C-terminal StrepII-tag protein in Sf9 insect cells.

PURITY: > 80% by SDS-PAGE.

SUPPLIED AS: _____ $\mu\text{g}/\mu\text{L}$ in 50 mM Tris, pH 8.0, 500 mM NaCl, 10% glycerol, 0.25 mM TCEP, as determined by AU_{280} .

STORAGE: -70°C . Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice/ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots ($<5 \mu\text{l}$) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 μg of RBC NSUN2. MW markers (left) are, from top, 200, 150, 120, 100, 85, 70, 60, **50**, 40, 30, 25, **20**, 15, & 10 kDa.

Enzymatic activity has not been tested.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com