

NRMT1 (N-terminal Xaa-Pro-Lys N-methyltransferase 1) (GST)

CATALOG NO.: HMT-11-319

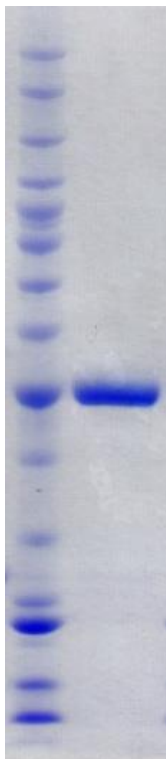
LOT NO.:

DESCRIPTION: Human recombinant NRMT1 (residues 2-223; Genbank Accession # NM_014064.3; MW = 52.3 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

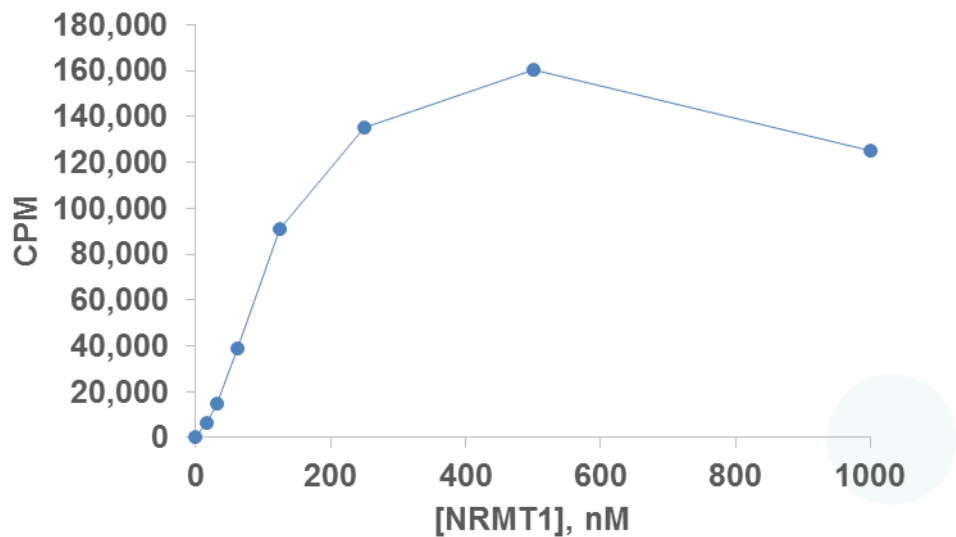
PURITY: >95% by SDS-PAGE

SUPPLIED AS: $_ \mu\text{g}/\mu\text{L}$ in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol

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 μl) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (12% acrylamide) of 4 μg of RBC NRMT1 (GST)
 MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, **20**, 15, 10 kDa.



Methyltransferase Activity of NRMT1 with RCC1 (HMT-11-321).
 Methylation determined as TCA-precipitable counts in a scintillation/filter plate assay. Reactions were 25 μL , 60 min., 30°C, with 1 μM [^3H]-SAM and 0.6 μM RCC1.

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