

NNMT (His) (Nicotinamide N-Methyltransferase)

CATALOG NO.: MT-11-358

DESCRIPTION: Human recombinant NNMT (residues 2-264; Genbank Accession # NM_006169.2; MW = 32.3 kDa) expressed with an N-terminal His tag in *E. coli*.

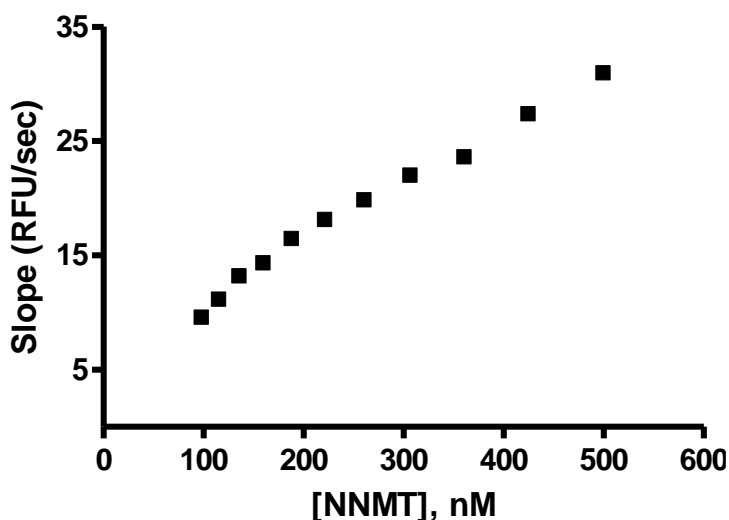
PURITY: > 95% by SDS-PAGE.

SUPPLIED AS: ___µg/µL in 50mM Tris-HCl, pH 7.5, 200mM NaCl, 5% glycerol, as determined by OD₂₈₀.

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 (4-12% acrylamide) of 2 and 4 µg of RBC NNMT. MW markers (right) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



NNMT methyltransferase activity using SAH hydrolase (SAHH)-coupled fluorescent assay. NNMT catalyzed methylation of nicotinamide is quantified using SAHH and Thioglo IV. 20 µL assays containing variable amounts of NNMT were incubated with 25 µM nicotinamide, 50 µM SAM and 20 µM TG IV (EMD Millipore). CLARIOstar plate reader (BMG Labtech) was used for real time monitoring with Ex/Em 400/465nm.

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