

HNMT (Histamine N-methyltransferase)

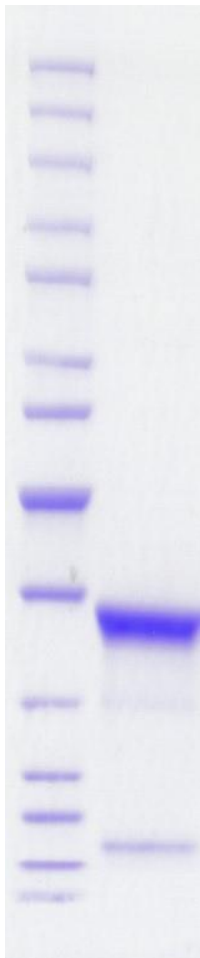
CATALOG NO.: MT-11-670

DESCRIPTION: Human recombinant HNMT (residues 2-292; Genbank Accession # NP_008826.1; MW = 38.1 kDa) expressed with an N-terminal His tag and C-terminal StrepII tag in *E. coli*.

PURITY: > 85% by SDS-PAGE.

SUPPLIED AS: ___µg/µL in 50mM Tris-HCl, pH 7.5, 500mM NaCl, 10% glycerol, 0.25mM TCEP as determined by OD₂₈₀.

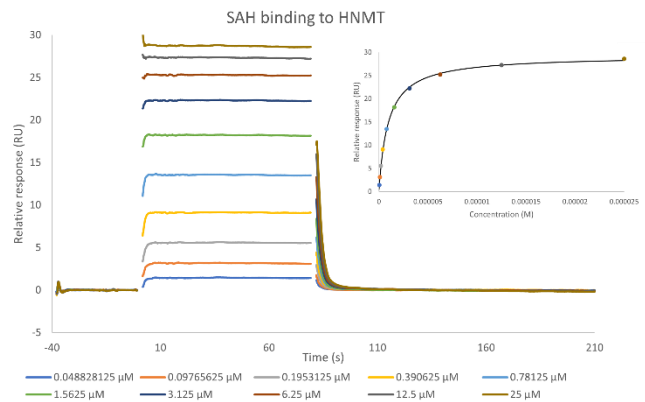
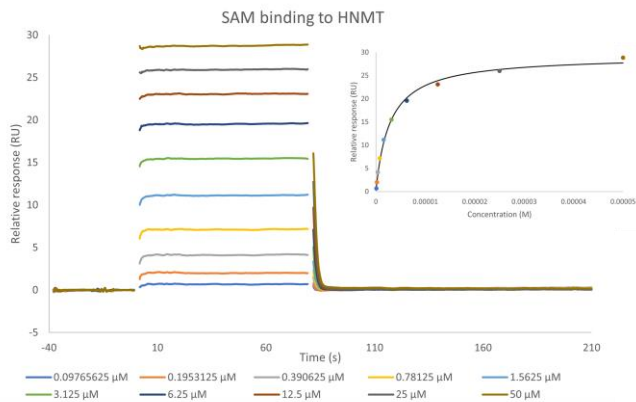
STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice/ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



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

*Enzymatic activity has not been tested.
Protein was active in SPR assay (see page 2)*

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



Binding Model	Target	Analyte	Chi ² (RU ²)	K _D (M)	R _{max} (RU)	Offset (RU)	Binding Model	Target	Analyte	Chi ² (RU ²)	K _D (M)	R _{max} (RU)	Offset (RU)
Steady state affinity	HNMT	SAM	5.91e-1	2.85e-6	28.9	0.4	Steady state affinity	HNMT	SAH	8.70e-2	9.53e-7	28.9	0.4

Titration of SAM (left) and SAH (right) against immobilized HNMT measured on a Biacore 8K+ in multi-cycle kinetics mode.

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