

## Data Sheet

### NNMT, His-Tag

Human, recombinant, C-terminal His-tag

**Catalog #:** 11357

**Lot#:** 140717

**Conc.:** 1.94 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 3 mM DTT, 145 mM Imidazole.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**References:**

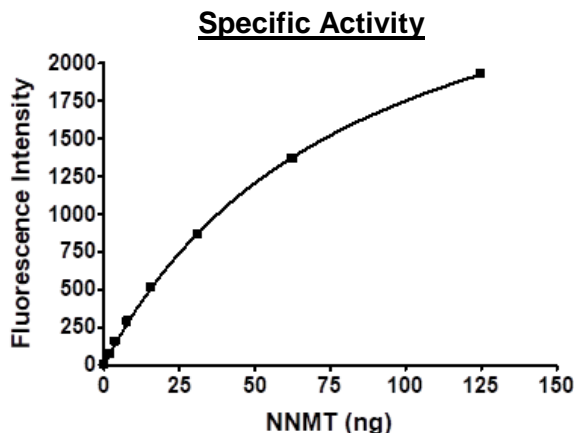
1. Tan, H.H. and Porter, A.G. *Biochem. Biophys. Res. Commun.* **382 (2)**, 462-467 (2009).
2. Fan, H. *et al.*, *World J. Gastroenterol.* **15 (16)**, 2020-2026 (2009).

**Description:** Human NNMT1 also known as Nicotinamide N-Methyltransferase, GenBank Accession No. NM\_006169, a.a. 1-264(full length), with C-terminal His-tag, expressed in an *E. coli* expression system. MW=30 kDa.

**Specific Activity:** 57 pmol/min/ $\mu\text{g}$ .  
 Assay Conditions: 10 mM Tris-HCl, pH 7.4, 0.05% Tween 20, 1 mM DTT, 100  $\mu\text{M}$  Quinoline, 100  $\mu\text{M}$  SAM, 30 min, room temperature.

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Assurance



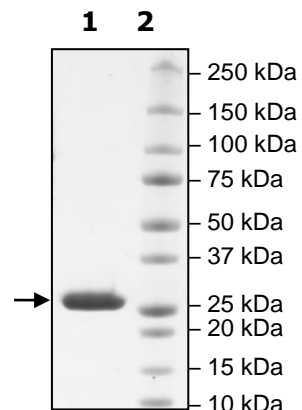
**4-20% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
 NNMT  
 (5.8  $\mu\text{g}$  loaded)

**Lane 2:**  
 Protein Marker

**Purity:**  $\geq 90\%$

**MW:** 30 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)