

## Data Sheet

### PRMT6, His-Tag

Human, recombinant, N-terminal His-tag

**Catalog #:** 51049

**Lot#:** 180320-2

**Conc.:** 0.32 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM imidazole, 20% glycerol, and 3 mM DTT.

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

**References:**

- Hyllus D. *et al.*, *J. Genes Dev.* **21** (24), 3369-3380 (2007).
- Guccione E. *et al.*, *Nature* **449** (7164), 933-937 (2007).

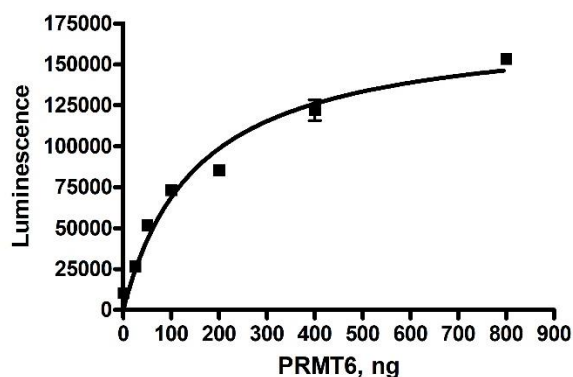
**Description:** Human PRMT6 also known as HRMT1L6 and protein arginine N-methyltransferase 6, GenBank Accession No. NM\_018137, a.a. 2-375(end), with N-terminal His-tag, expressed in an *E. coli* expression system. MW=43 kDa.

**Specific Activity:** 0.036 pmol/min/ $\mu\text{g}$   
**Assay Conditions:** For standard, asymmetric H4 R3(2Me)-Biotin was coated onto a neutravidin plate at varying concentrations. Biotin was coated onto a neutravidin plate at 0.4  $\mu\text{g}/\text{ml}$ . Reaction mixture contained 50 mM Tris buffer pH 8, 10 mM NaCl, 1 mM DTT, 0.05% Tween-20, 40  $\mu\text{M}$  S-adenosylmethionine, and various amounts of PRMT6—incubated for 1 hour at room temperature. Followed by 1-hour incubation with primary antibody against asymmetric H4 R3(2Me), and 30-minute incubation with secondary antibody. Luminescence was developed with HRP substrate A and B.

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

### Quality Assurance

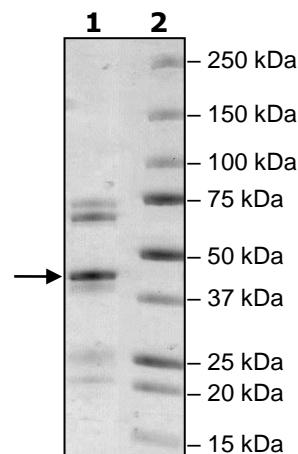
**Specific Activity**



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

**Lane 1:**  
2  $\mu\text{g}$  PRMT6  
**Lane 2:**  
Protein Marker

**MW:** 43 kDa  
**Purity:**  $\geq 41\%$



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)