

## Data Sheet

### PRMT1

Human, recombinant, N-terminal GST tag

**Catalog #:** 51040

**Lot#:** 110915      **Conc.:** 0.13 mg/ml

**Formulated in:** 40mM Tris pH 8.0, 110mM NaCl, 2.2mM KCl, 0.04% Tween, 20% Glycerol, 3mM DTT.

**Stability:** >6 months at -80°C

#### References:

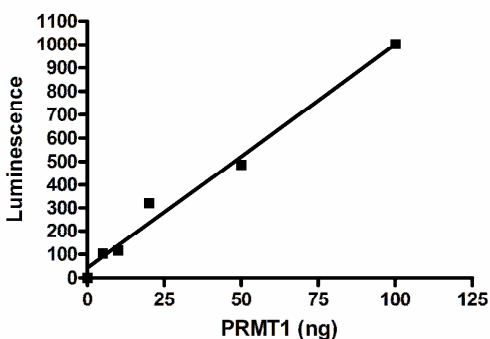
1. Le Romancer, M. *et al.*, *Mol. Cell* **31** (2), 212-221 (2008)
2. Fronz, K. *et al.*, *J. Biol. Chem.* **283** (29), 20408-20420 (2008)

**Description:** Human PRMT1 (GenBank Accession No. NM\_001536), amino acids 2-end, with N-terminal GST tag, MW= 68 kDa, expressed in an *E. coli* expression system.

**Specific Activity:** 0.52 pmol/min/ug.  
 Assay conditions: 50 µl reaction mix (20 mM phosphate buffer pH7.4, 20 µM S-adenosyl methionine, and 1-5 ng methyltransferase PRMT1) add to the wells coated with the substrate. Incubate at room temperature for 1 hr. Add antibody against methylated R3 residue of histone H4, incubate 1 hr. Then, add secondary HRP-labeled antibody and incubate 30 min. Finally, add HRP chemiluminescent substrates and read luminescence.

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

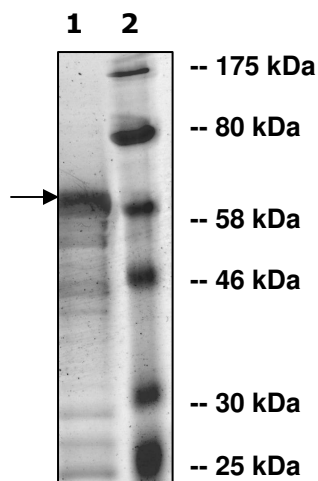
### Quality Assurance



### 10% SDS-PAGE Coomassie staining

Lot 110316G  
**Lane 1:**  
 1.5 µg PRMT1  
**Lane 2:**  
 Protein Marker  
 BioLabs (#P7708L)

**MW:** 68 kDa  
**Purity:** ≥49%



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

To place your order, please contact us by Phone **1.858.829.3082** 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)