

## Data Sheet

### SUV39H1, GST-tag

Human, recombinant, N-terminal GST tag

**Catalog #:** 51070

**Lot #:** 140817-G1

**Conc.:** 0.04 mg/ml

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

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

1. Kang, M.Y. *et al.*, *Int. J. Cancer* **121** (10), 2192-2197 (2007).
2. Spensberger, D. and Delwel, R. *FEBS Lett.* **582** (18), 2761-2767 (2008).

**Description:** Human SUV39H1 (GenBank Accession No. NM\_003173), amino acids 82-412, with N-terminal GST tag, MW= 64 kDa, expressed in *E. coli*.

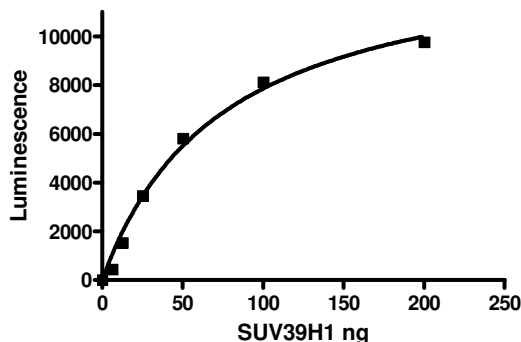
**Specific Activity:** 2.4 pmol/min/ $\mu\text{g}$ .

Assay conditions: 50  $\mu\text{l}$  reaction mix (50 mM TRIS pH 8.8, 5mM MgCl<sub>2</sub>, 4mM DTT, 0.05% Tween-20, 2.5  $\mu\text{M}$  S-adenosylmethionine, and SUV39H1 (0 – 200 ng) add to the wells coated with the substrate on G bioscience Neutravidin black plate. Incubate for 1 hr. Add antibody against methylated K9 residue of histone H3, 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

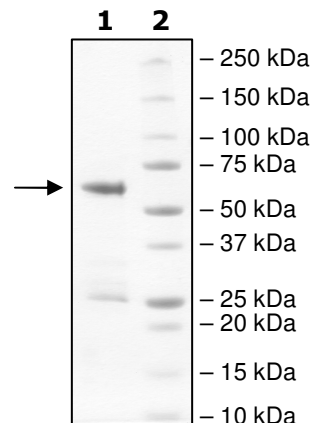
**Specific Activity**



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

**Lane 1:**  
 2 $\mu\text{g}$  SUV39H1  
**Lane 2:**  
 Protein Marker

**MW:** 65 kDa  
**Purity:**  $\geq 80\%$



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)