

Data Sheet

PRMT4/CARM1 (HEK293)

Human, recombinant, N-terminal FLAG tag
Catalog #: 51047
Lot#: 130204 **Conc.:** 0.2 mg/ml

Formulated in: 40 mm Tris-HCl, pH 8.0,
 110 mm NaCl, 2.2 mm KCl, 80 ng/μl flag
 peptide, 20% glycerol, and 3 mM DTT.

Stability: >6 months at -80 °C

References:

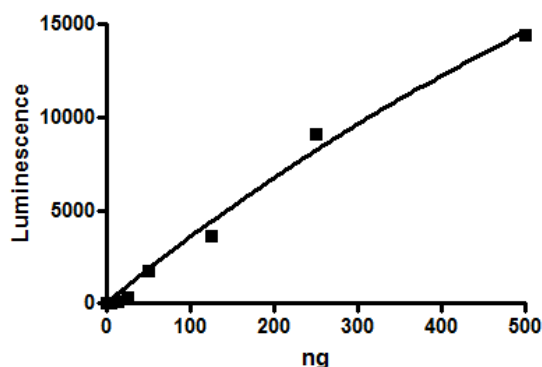
1. Li H. *et al.*, *J. Biol. Chem.* **277 (47)**, 44623-44630 (2002).
2. Frieze S. *et al.*, *Cancer Res.* **68 (1)**, 301-306 (2008).

Description: Human PRMT4/CARM1 (GenBank Accession No. NM_199141), amino-acids 2-end, with N-terminal FLAG tag, MW=66 kDa, expressed in a FreeStyle 293-F cells.

Assay conditions: 50 μl reaction mix (1x HMT Assay Buffer 5, BPS Cat# 52191 and 5-1000 ng methyltransferase PRMT4) add to the wells coated with the substrate on a Thermo HBC Neutravidin black plate. Incubate for 1 hr. Add antibody against methylated substrate. Incubate for 1 hr. Then, add secondary HRP-labeled antibody and incubate for 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

Lane 1:
 2 μg PRMT4/CARM1
Lane 2:
 Protein Marker

MW: 66 kDa
Purity: ≥65%

