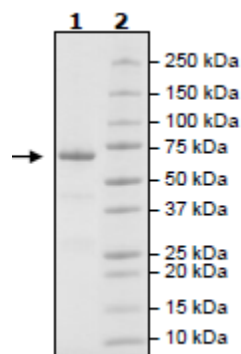


Product Information

Construct:	PRMT1(His-GST-2-371(end))
Concentration:	0.54 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20 and 20% glycerol
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_001536
MW:	70 kDa
Purity:	86%
Assay Conditions:	Enzyme was incubated with 0.25 μ M of biotinylated H4 1-21 peptide with 1 μ M AdoMet in HMT buffer 2A in a 10 μ l reaction. 5 μ l of Anti-Rabbit AlphaLISA acceptor beads (1:250) and 5 μ l of Primary Antibody 4 (1:20) were added and incubation continued for 30 min. Streptavidin Donor beads (1:125) were added (10 μ l) and mixture was further incubated for 15 min at room temperature before A-counts were measured.
Applications	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

