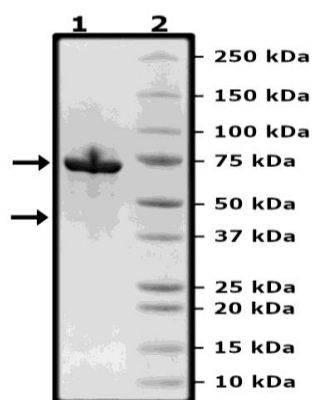


## Product Information

<b>Construct:</b>	PRMT5 (FLAG-2-637(end)) / MEP50 (His-2-342(end))
<b>Concentration:</b>	0.32 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution.
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	PRMT5: NM_006109; MEP50: NM_024102
<b>MW:</b>	PRMT5: 73 kDa; MEP50: 38 kDa
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	Assay was performed according to PRMT5 Chemiluminescent Assay Kit (BPS Bioscience #52002L) with PRMT5 (BPS Bioscience #51048) enzyme titrated from at various concentrations.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PRMT5 Activity

