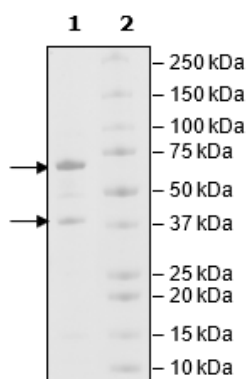


Product Information

Construct:	PRMT5 (FLAG-2-637(end)) / MEP50 (His-2-342(end))
Concentration:	1.27 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT, and 100 µg/ml FLAG peptide
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	PRMT5: NM_006109; MEP50: NM_024102
MW:	PRMT5: 74 kDa; MEP50: 38 kDa
Purity:	≥90%
Assay Conditions:	Assay was performed according to the PRMT5 Chemiluminescent Assay Kit (BPS Bioscience #52002L) with PRMT5 enzyme titrated from 800 ng/rxn – 0.78 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



PRMT5/MEP50 Activity

