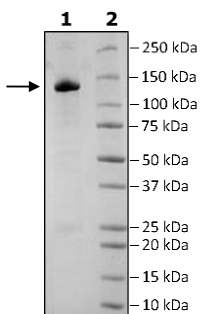


Product Information

Construct:	DNMT1 (His-FLAG-501-1632(end))
Concentration:	0.81 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 ng/μl FLAG peptide, 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_001130823
MW:	130 kDa
Purity:	≥90%
Assay Conditions:	Assay was performed using the DNMT1 Chemiluminescent Assay Kit (BPS Bioscience #52050L) with DNMT1 titrated from 800 ng/rxn – 12.5 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



DNMT1 Activity

