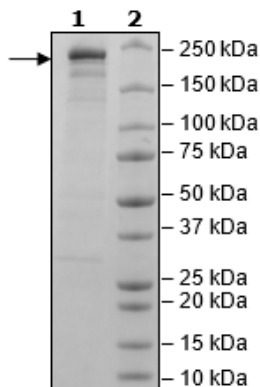


Product Information

Construct:	DNMT1 (GST-2-1632(end))
Concentration:	0.62 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 3 mM DTT
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:	211 kDa
Purity:	78%
Assay Conditions:	Assay was performed according to the protocol in DNMT1 Chemiluminescent Assay Kit (BPS Bioscience #52050L) with DNMT1 titrated at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



DNMT1 Activity

