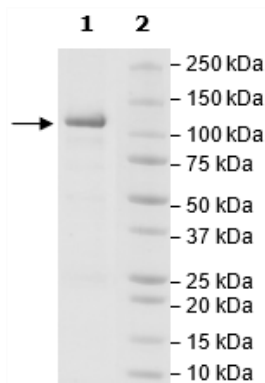


Product Information

Construct:	SetDB1 (His-GST-510-1290(end))
Concentration:	0.34 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 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_012432
MW:	114 kDa
Purity:	85%
Assay Conditions:	Assay run according to SetDB1 Chemiluminescent Assay Kit (BPS Bioscience #51056).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



SetDB1 Activity

