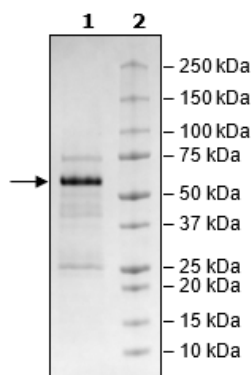


Product Information

Construct:	SUV39H2 (GST-26-350(end))
Concentration:	0.22 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 20 mM glutathione
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_024670
MW:	63 kDa
Purity:	62%
Assay Conditions:	Assay was done according to SUV39H2 Activity Assay Kit (BPS Bioscience #52008L) with various amounts of SUV39H2.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



SUV39H2 Activity

