

Product Information

Construct:	DNMT3a (GST-2-912(end))
Concentration:	0.20 mg/ml
Species:	Human
Formulated In:	45 mM Tris-HCl pH 8.0, 124 mM NaCl, 2.4 mM KCl, 0.04% Tween-20, 3 mM DTT, 10% glycerol, and 2 mM glutathione
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_175629
MW:	128 kDa
Purity:	10%

Quality Control Data

4-20% SDS-Page Coomassie Staining

