

Product Information

Construct:	WRN (GST-517-1093)
Concentration:	0.58 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 200 mM NaCl, 2.2 mM KCl, 0.05% Tween-20, 10 mM glutathione, 1 mM TCEP, 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_000553
MW:	92 kDa
Purity:	89%

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

