

Product Information

Construct:	Axl (GST-473-894(end))
Concentration:	0.50 mg/ml
Species:	Human
Formulated In:	25 mM Tris-HCl, pH 8.0, 69 mM NaCl, 1.3 mM KCl, 0.02% Tween-20, 50% 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_021913
MW:	74 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

