

Product Information

Construct:	SOS1 (FLAG-594-1049-Avi)
Concentration:	0.55 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, 80 µg/ml FLAG peptide
Expression System:	Sf9
Format:	Aqueous buffer solution.
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession No.:	NM_005633
MW:	57 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

