

Product Information

Construct:	GRP78 (FLAG-19-654(end))
Concentration:	0.26 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 100 µg/ml FLAG peptide, 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_005347
MW:	72 kDa
Purity:	87%

Quality Control Data

4-20% SDS-Page Coomassie Staining

