

Product Information

Construct:	MMP13 (20-471(end)-His)
Concentration:	0.19 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 200 mM imidazole
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_002427
MW:	53 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

