

Product Information

Construct:	PTPRQ (His-2003-2259)
Concentration:	1.30 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT, 200 mM imidazole
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_001145026
MW:	31 kDa
Purity:	90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

