

Product Information

Construct:	DPY30 (His-2-99(end))
Concentration:	3.07 mg/ml
Species:	Human
Formulated In:	45 mM Tris-HCl pH 8.0, 124 mM NaCl, 2.4 mM KCl, 225 mM Imidazole, 1 mM TCEP, 1 μM ZnCl ₂ , 10% Glycerol.
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_032574
MW:	12 kDa
Purity:	90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

