

Product Information

Construct:	HSP40 (1-340(end)-His)
Concentration:	1.49 mg/ml
Species:	Human
Formulated In:	25 mM Tris-HCl, pH 7.5, 400 mM NaCl, 0.05% Tween-20, 10% glycerol and 3 mM DTT.
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_006145
MW:	40 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

