

Product Information

Construct:	PSMA(His-Avi-44-750(end))
Concentration:	0.44 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110mM NaCl, 2.2mM KCl, 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_004476
MW:	84 kDa + glycans
Glycosylation:	This protein runs at a higher molecular weight due to glycosylation.
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

