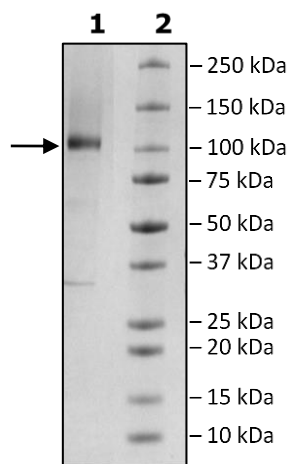


Product Information

Construct:	PSMA (His-Avi-44-750(end))-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
Concentration:	0.40 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM 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 MW due to glycosylation.
Purity:	83%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown

