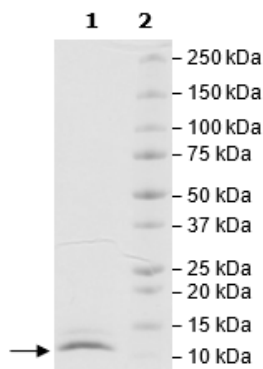


Product Information

Construct:	Ubiquitin (His-Avi-2-76)-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag technology. Biotinylation is confirmed to be $\geq 90\%$.
Concentration:	0.91 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 100 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_021009
MW:	12 kDa
Purity:	90%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown

