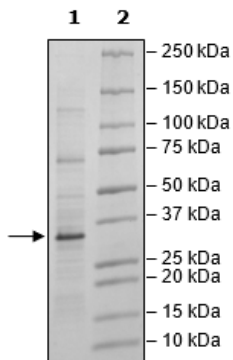


Product Information

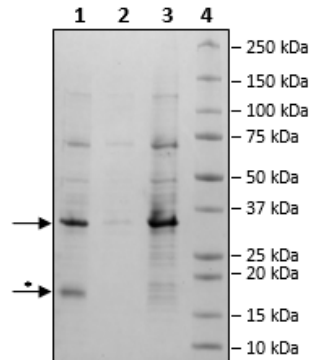
Construct:	LIN28A (FLAG-Avi-2-209(end))-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag technology. Biotinylation is confirmed to be $\geq 90\%$.
Concentration:	0.21 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT, and 100 $\mu\text{g}/\text{mL}$ FLAG peptide
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_024674
MW:	26 kDa
Purity:	53%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown



1. Beads
2. Flow Thru
3. Control
4. Standards

* Avidin from beads.