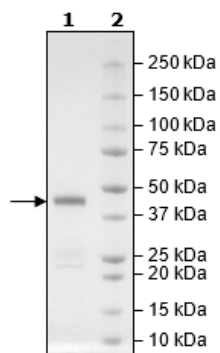


## Product Information

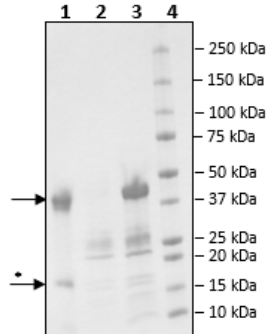
<b>Construct:</b>	CBL-B (His-Avi-39-426)-(Biotin)
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
<b>Concentration:</b>	0.4 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	45 mM Tris-HCl pH 8.0, 124 mM NaCl, 2.4 mM KCl, 10% glycerol, 3 mM DTT
<b>Expression System:</b>	<i>E. coli</i>
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_170662
<b>MW:</b>	48 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining



### Biotin-Avidin Pulldown



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

\* Avidin from beads.