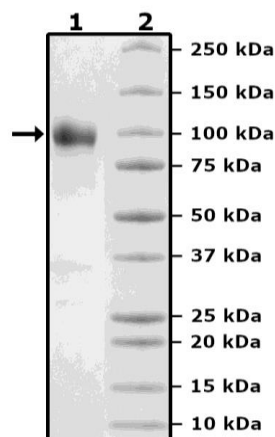


## Product Information

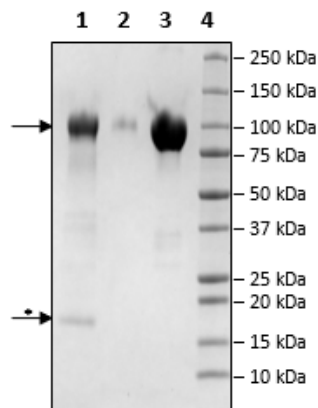
<b>Construct:</b>	TPBG (32-355-Fc(IgG1)-Avi)-(Biotin)
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
<b>Concentration:</b>	1.62 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>Expression System:</b>	HEK293
<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_006670
<b>MW:</b>	64 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<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.