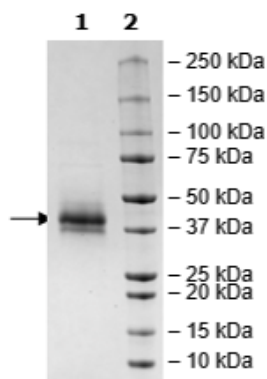


## Product Information

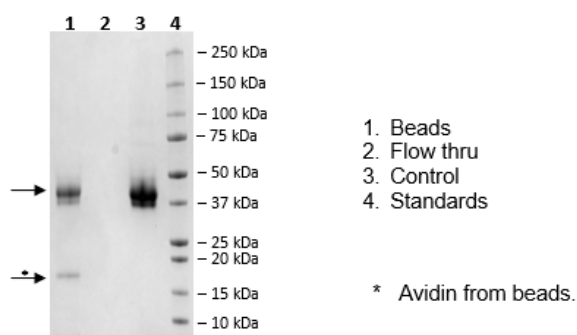
<b>Construct:</b>	IL2RB (27-239-Avi-FLAG)-(Biotin)
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
<b>Concentration:</b>	1.13 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% 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_000878
<b>MW:</b>	28 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Purity:</b>	≥90%
<b>Aggregation:</b>	<10%

## Quality Control Data

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



### Biotin-Avidin Pulldown



Gel Filtration Trace

