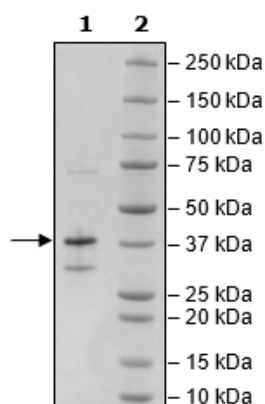


Product Information

Construct:	BCMA (1-54-Fc(IgG1)-Avi)-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%
Concentration:	0.20 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
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_001192
MW:	34 kDa
Purity:	74%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown

