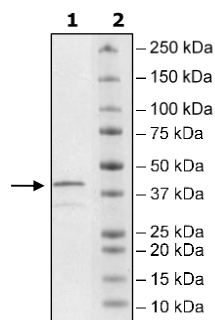


Product Information

<b>Construct:</b>	BCMA (1-54-Fc (IgG1)-Avi)
<b>Concentration:</b>	1.11 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_001192
<b>MW:</b>	34 kDa +glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Purity:</b>	87%
<b>Assay Conditions:</b>	Assay was performed according to BCMA:APRIL[Biotinylated] Inhibitor Screening Assay Kit (BPS Bioscience #79722) with APRIL[Biotinylated] enzyme titrated from 40 ng/rxn – 0.03 ng/rxn.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



BCMA:APRIL[Biotinylated] Binding Interaction

