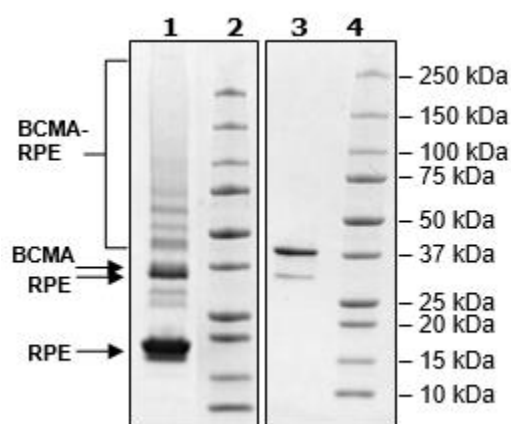


## Product Information

<b>Construct:</b>	BCMA (1-54-Fc(IgG1)-Avi)-(PE)
<b>Label:</b>	R-Phycoerythrin (PE) is an oligomeric protein complex (270 kDa) from red algae that exhibits intensely bright red-orange fluorescence with high quantum yields. The complex consists of six heterodimers, $\alpha$ subunit (18 kDa) and $\beta$ -subunit (20 kDa), and an additional $\gamma$ -subunit (34 kDa). PE is covalently attached randomly through lysines on the target protein.
<b>Concentration:</b>	1.97 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. Protect from light.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_001192
<b>MW:</b>	34 kDa + glycans + PE
<b>Glycosylation:</b>	This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.
<b>Purity:</b>	$\geq 90\%$
<b>Applications:</b>	Useful for labeling cells expressing Anti-BCMA antibodies for flow cytometry and immunofluorescence microscopy.

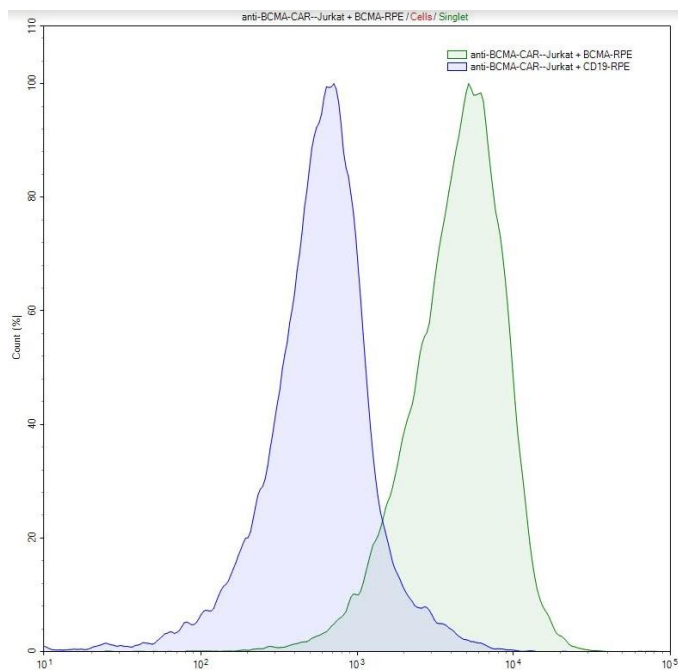
## Quality Control Data

## 4-20% SDS-Page Coomassie Staining



**Lane 1:** 8 µg BCMA-PE  
**Lane 3:** 4 µg BCMA  
**Lane 2 & 4:** Protein Marker

## FACS Assay



BCMA, Fc-Fusion, Avi-Tag, PE-Labeled (BPS Bioscience #100733) was tested using FACS with Anti-BCMA CAR /NFAT (Luciferase) Reporter Jurkat Cell Line (BPS Bioscience #79694) and Anti-CD19 CAR / NFAT (Luciferase) Reporter Jurkat Cell Line (CD19 SCFV-CD28-4-1BB-CD3ζ) (BPS Bioscience #79853).