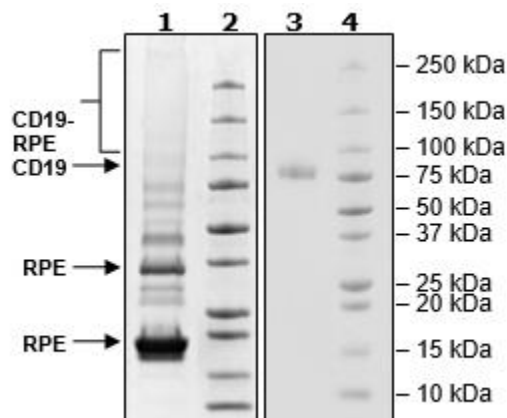


Product Information

Construct:	CD19 (21-291-Fc (IgG1)-Avi)-(PE)
Label:	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, α subunit (18 kDa) and β -subunit (20 kDa), and an additional γ -subunit (34 kDa). PE is covalently attached randomly through lysines on the target protein.
Concentration:	1.25 mg/ml
Species:	Human
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles. Protect from light.
Storage:	-80°C
Genbank Accession:	NM_001770
MW:	59 kDa + glycans + PE
Glycosylation:	This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.
Purity:	≥90%
Application:	Useful for labeling cells expressing Anti-CD19 antibodies for flow cytometry and immunofluorescence microscopy.

Quality Control Data

4-20% SDS-Page Coomassie Staining

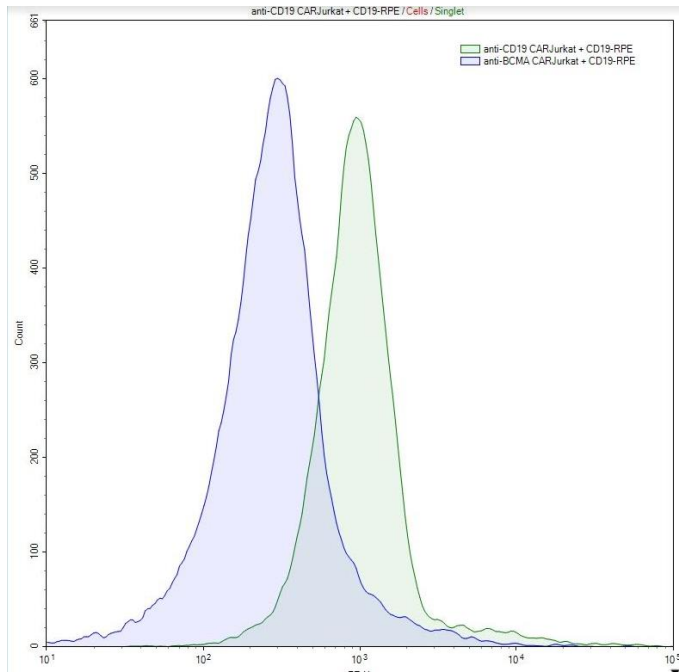


Lane 1: 8 μ g CD19-PE

Lane 3: 4 μ g CD19

Lanes 2 & 4: Protein Marker

FACS Assay



CD19, Fc-Fusion (IgG1), Avi-Tag, PE-Labeled (BPS Bioscience #100732) 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-CD3 ζ) (BPS Bioscience #79853).