

Product Information

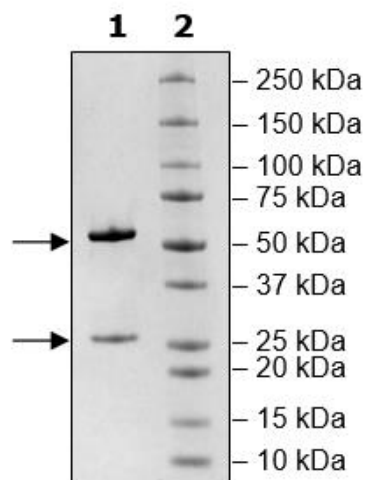
Description:	This anti-CD19 IgG antibody is a purified recombinant antibody, which was enzymatically biotinylated using Avi-tag™ technology. This antibody has been tested for specific binding to purified human CD19 protein (BPS Bioscience #101015).
Construct:	Anti-CD19 LC (1-240(end)) / Anti-CD19 HC (1-490(end))-Fc (Human IgG1)-Avi
Concentration:	1.72 mg/ml
Species:	Human
Isotype:	IgG1
Formulated In:	8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
Expression System:	HEK293
Clonality:	Monoclonal
Purification:	Protein A affinity chromatography from HEK293 supernatants
Cross Reactivity:	This antibody recognizes human CD19. It has not been tested with other species.
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	Anti-CD19 HC: 52 kDa; Anti-CD19 LC: 24 kDa
Purity:	≥90%
Assay Conditions:	<i>Experimental design and assay protocol for measuring biotinylated anti-CD19 specific binding to His-tagged CD19 recombinant protein in ELISA:</i>

1. Coat a clear 96-well nickel-coated plate with purified human his-tagged CD19 (BPS Bioscience #101015) at 1 µg/ml in PBS. Incubate overnight at 4°C to bind protein to surface
2. Dilute 3x BPS Immuno Buffer 1 (BPS Bioscience #79311) with sterile distilled water 1:3 to make 1x Immuno Buffer 1
3. Wash each well with 100 µl of 1x BPS Immuno Buffer 1 three times. Tap upside down on paper towels or absorbent pads.
4. Block each well with 100 µl of Blocking Buffer 2 (BPS Bioscience #79728) for 1 hour at room temperature (slow shaking).
5. After blocking, remove blocking buffer and tap to dry.
6. Add serial dilutions of the antibody (300 nM to 0 nM in 3-fold dilutions recommended for initial analysis). Incubate for 1 hour at room temperature (with slow shaking).
7. Wash wells three times and incubate with HRP-Streptavidin (1:1000, BPS Bioscience #79742) for 30 minutes at room temperature (with slow shaking).
8. Wash again three times. Tap to dry.
9. For detection, add 100 µl of Colorimetric HRP Substrate (BPS Bioscience #79651) for 1-10 minutes until a blue color develops in the positive control.
10. Once color develops in the positive control and samples, the reaction must be immediately quenched with an equal volume (100 µl) of 1N HCL
11. Absorbance is measured at 450 nm.

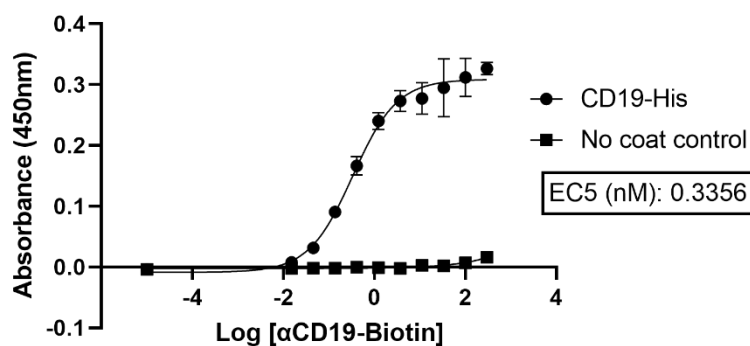
Applications: Useful for studying the inhibition of CD19 in ELISA and in cellular assays.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Binding Assay



Biotinylated anti-CD19 (BPS Bioscience #101093) binding assay to recombinant CD19 (BPS Bioscience #101015). The ELISA was performed in replicates following the assay conditions described and the EC50 of this interaction determined using increasing concentrations of antibody

Related Products

Products	Catalog #	Size
Anti-CD19 IgG Antibody	100981	50 µg/100 µg
Anti-CD19-Anti-CD3 Bispecific Antibody	100441	50 µg/100 µg
Anti-BCMA-Anti-CD19-Anti-CD3-His Trispecific Antibody	100761	50 µg/100 µg
CD19, Avi-His-Tag	101015	100 µg
CD19, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	79475	25 µg/50 µg