

Product Information

Description:	Biotinylated anti-CD20 IgG antibody is a purified recombinant antibody which recognizes human CD20 antigen. This antibody has been tested for specific binding to purified human CD20 protein.
Concentration:	1.24 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol More than 90% biotinylation
Expression System:	Heavy chain (HC) and Light chain (LC) co-expressed in HEK293
Purification:	Protein A affinity
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	Heavy Chain: 53 kDa; Light Chain 23 kDa
Purity:	≥90%
Assay Conditions:	Functional validation: The antibody was validated by ELISA using a purified CD20 antigen. The CD20 protein was coated onto a 96-well plate overnight at 4°C (50 µl/well at a concentration of 4 µg/ml in PBS). The plate was washed 3 times with Immuno Buffer 1 (BPS Bioscience #79311) and blocked using 100 µl of Blocking Buffer 2 (BPS Bioscience #79728) for 1 hour at room temperature. After removing the blocking buffer, 50 µl/well of purified biotinylated anti-CD20 antibody (BPS Bioscience #101207), serially diluted in Blocking Buffer 2, was added for 30 minutes at room temperature. After 3 more washes, the plate was incubated with Streptavidin-HRP, washed, and incubated with the Colorimetric HRP substrate. The reaction was stopped, and absorbance was read at 450 nm. The Blank value was subtracted from all values.
Applications:	This product is for research use only. It is not suitable for human diagnostic or therapeutic use.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

CD20:anti CD20-Biotin Binding Assay

