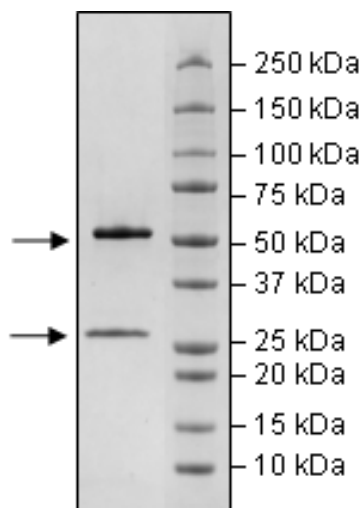


Product Information

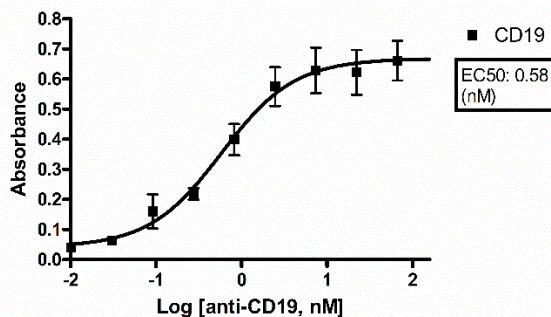
Description:	This purified recombinant monoclonal antibody is a human Anti-CD19 IgG which recognizes CD19 antigen. The variable regions for this antibody are based on an FDA-approved drug. This antibody has been tested for specific binding affinity to purified human CD19 protein (BPS Bioscience, #101015) in a colorimetric ELISA binding assay.
Concentration:	1.53 mg/ml
Species:	Human
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	Heavy chain (HC) and Light chain (LC) co-expressed in HEK293 cells
Purification:	Protein A affinity purification from HEK293 cells.
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	Anti-CD19 HC: ~49.9 kDa; Anti-CD19 LC: ~23.9 kDa
Purity:	≥90%
Assay Conditions:	<p><i>Experimental design and assay protocol for measuring anti CD19 specific binding to CD19 antigen in ELISA assay:</i></p> <p>Purified human his-tagged CD19 was bound to clear 96-well nickel plate overnight in 4°C (1ug/mL). Upon washing and blocking wells for 1 hour at room temp, serial dilutions of anti-CD19 (200 nM to 0nM in 3-fold dilutions) were incubated in each well for 1 hr at room temp (slow shaking). Next, wells were washed and incubated with anti-human IgG-HRP for 1 hour at room temp (slow shaking) then washed again. For detection, the wells were incubated with Colorimetric HRP Substrate (BPS Bioscience #79651) for 1-10 minutes until a blue color developed in the positive control. The reaction was then quenched with equal volume of 1N HCL and absorbance read at 450 nm.</p>
Applications:	The anti-CD19 IgG format antibody can be used for labeling cells expressing CD19 for flow cytometry and immunofluorescence microscopy.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Binding affinity of Anti-CD19 IgG to target protein in ELISA assay



Purified CD19-His-Tag (1 $\mu\text{g/ml}$) was bound to a clear, 96-well, nickel-coated plate. Anti-CD19 (0 - 200 nM) was added and following incubation, the binding affinity was determined using a colorimetric detection assay.