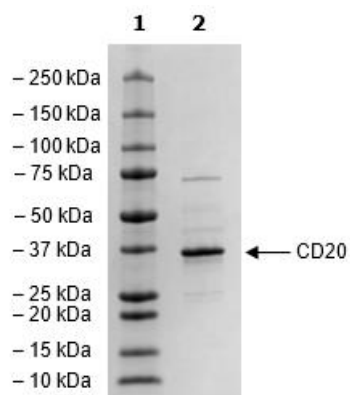


## Product Information

|                              |   |
|------------------------------|---|
| <b>Description:</b>          | Recombinant human CD20, also known as membrane spanning 4-domainsA1 (MS4A1), full length, encompassing amino acids 1-297(end). This construct contains a C-terminal FLAG-tag. This recombinant protein was affinity purified.   |
| <b>Species</b>               | Human   |
| <b>Construct:</b>            | CD20 (1-297(end)-FLAG)  |
| <b>Concentration:</b>        | 0.42 mg/ml  |
| <b>Expression System:</b>    | HEK293  |
| <b>Purity:</b>               | 85%   |
| <b>Format:</b>               | Aqueous buffer solution.  |
| <b>Formulated In:</b>        | 50 mM HEPES (pH 8), 150 mM NaCl, 10% (vol/vol) glycerol, 0.1% DDM, 0.01% CHS, and 100 µg/ml FLAG peptide  |
| <b>MW:</b>                   | 34 kDa  |
| <b>Genbank Accession:</b>    | NM_021950.3   |
| <b>Stability:</b>            | At least 6 months at -80°C.   |
| <b>Storage:</b>              | -80°C   |
| <b>Instructions for Use:</b> | Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.  |
| <b>Assay Conditions:</b>     | The protein was validated by measuring its binding to Anti-CD20 antibody, Biotin-labeled (BPS Bioscience #101207) in biochemical binding (ELISA) assay. The CD20 protein was coated onto a 96-well plate overnight at 4°C (50 µl/well at a concentration of 2 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) with DDM/CHS detergent (dodecyl-β-D-maltoside (DDM) 0.05%, cholesteryl hemisuccinate (CHS) 0.01%) for 1 hour at room temperature. After removing the blocking buffer, 50 µl/well of purified anti-CD20 antibody, biotin-labeled, serially diluted in Blocking Buffer 2 with DDM/CHS detergent, was added for 60 minutes at room temperature. After 3 more washes, the plate was incubated with HRP Streptavidin (BioLegend #405210), washed, and incubated with Colorimetric HRP substrate (BPS Bioscience #79651). The reaction was stopped, and absorbance was read at 450 nm. The Blank value was subtracted from all values. |
| <b>Applications:</b>         | Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.   |

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining



### CD20: Anti-CD20 Binding Assay

