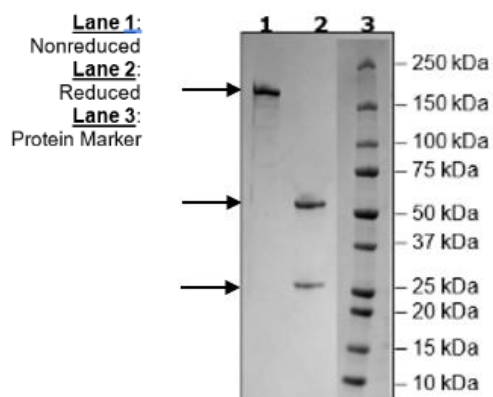


Product Information

Description:	Anti-CD2 antibody is a human recombinant antibody which recognizes human CD2 antigen. This antibody has been tested for specific binding to human CD2 (BPS Bioscience #101239).
Concentration:	0.12 mg/ml
Species:	Human
Isotype:	IgG
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
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:	150 kDa
Purity:	≥90%
Assay Conditions:	Functional validation: The antibody was validated by measuring anti-CD2 binding to CD2 antigen in ELISA assay. The anti-CD2 antibody 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) for 1 hour at room temperature. After removing the blocking buffer, 50 µl/well of purified biotinylated CD2-Avi-His protein (BPS Bioscience #101239), serially diluted in Blocking Buffer 2, was added for 30 minutes at room temperature. After 3 more washes, the plate was incubated with Streptavidin-HR, 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



Anti-CD2:CD2-Biotin Binding Assay

