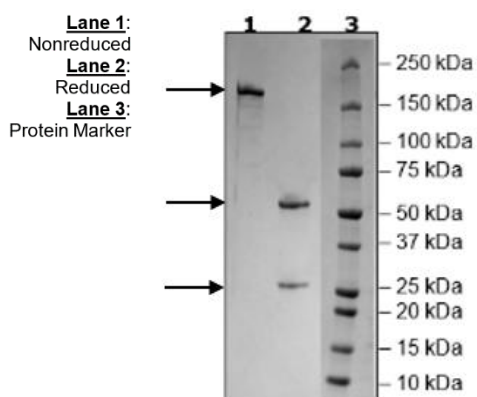


## Product Information

<b>Description:</b>	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).
<b>Concentration:</b>	0.12 mg/ml
<b>Species:</b>	Human
<b>Isotype:</b>	IgG
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>Expression System:</b>	Heavy chain (HC) and Light chain (LC) co-expressed in HEK293
<b>Purification:</b>	Protein A affinity
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 12 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>MW:</b>	150 kDa
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	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.
<b>Applications:</b>	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

