

Product Information

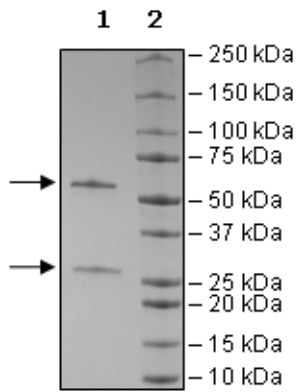
Description:	This anti-NCAM1 (CD56) IgG antibody is a purified recombinant antibody, which was enzymatically biotinylated using Avi-tag™ technology. This antibody has been tested for specific binding affinity to purified human NCAM1 (aka CD56) protein (BPS Bioscience #101043).
Concentration:	0.55 mg/ml
Species:	Human
Formulated In:	8mM phosphate pH 7.4, 110mM NaCl, 2.2mM KCl, and 20% glycerol
Expression System:	Heavy chain (HC) and Light chain (LC) co-expressed in HEK293
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:	150 kDa
Purity:	≥99%
Assay Conditions:	<i>Experimental design and assay protocol for measuring anti-CD56/NCAM1 specific binding to CD56/NCAM1 antigen in ELISA assay:</i>

- 1. Purified human His-tagged NCAM1 (CD56) (BPS Bioscience #101043) was bound to a clear 96-well nickel plate overnight at 4°C (1 µg/ml in PBS).*
- 2. Wells were washed with BPS Immuno Buffer 1 (BPS Bioscience #79311) three times and tapped upside down on paper towels or absorbent pads.*
- 3. Each well was blocked with 100 µl of blocking buffer (BPS Bioscience #79728) for 1 hour at room temperature (slow shaking).*
- 4. Serial dilutions of the antibody (600 nM to 0nM in 3-fold dilutions) were incubated in each well for 1 hour at room temperature (slow shaking).*
- 5. Next, wells were washed three times and incubated with HRP-Streptavidin (1:1000, BPS Bioscience #79742) for 30 minutes at room temp (slow shaking)*
- 6. Wells were washed again three times with BPS Immuno Buffer 1 and tapped to dry.*
- 7. 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.*
- 8. The reaction was then immediately quenched with an equal volume of 1N HCl and absorbance was measured at 450 nm.*

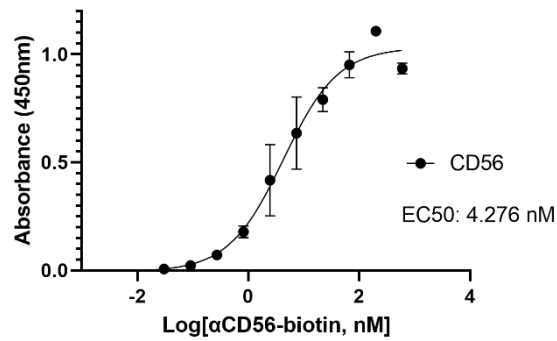
Applications: This product is for research use only. It is not suitable for human diagnostic or therapeutic use. The anti-NCAM1 (CD56) IgG format antibody can be used for flow cytometry and immunofluorescence microscopy.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Binding Assay



Purified NCAM1 (CD56) His-tagged protein (BPS Bioscience #101043) was immobilized on the surface of a microtiter plate, then tested for binding using increasing amounts of biotin-labeled NCAM1 (CD56) antibody (BPS Bioscience #101112) in a colorimetric assay, as described in Assay Conditions.

Related Products:

Products	Catalog #	Size
NCAM1, Avi-His-Tag, HiP™	101043	Multiple Sizes
NCAM1, Avi-His-Tag, Biotin-Labeled	101044	Multiple Sizes