

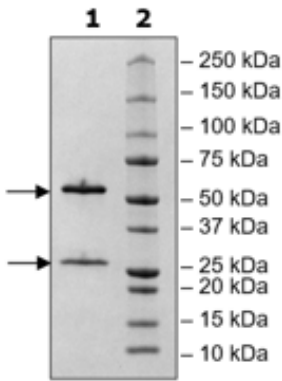
Product Information

Description:	Anti-CD235a antibody is a purified biotinylated recombinant human antibody which recognizes the CD235a/GYPA protein. Glycophorin A is the major intrinsic membrane protein of the erythrocyte. The N-terminal sialoglycosylated segment, which lies outside the erythrocyte membrane, has MN and Ss blood group receptors. This antibody has been tested for specific binding to purified human CD235a protein (BPS Bioscience #101196) in a colorimetric ELISA binding assay.
Concentration:	1.07 mg/ml
Isotype:	Human IgG1
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2mM KCl, 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
Tag:	C-terminal Avi-Tag™
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	~150 kDa (HC: 51 kDa + glycans; LC: 23 kDa)
Purity:	≥90%
Assay Conditions:	<p><i>Experimental design and assay protocol for measuring anti-CD235a specific binding to CD235a antigen in ELISA assay:</i></p> <ol style="list-style-type: none"><i>1. Purified human CD235a protein (cat#101196) was coated onto a clear 96-well nickel plate overnight at 4°C (1 µg/ml in PBS, 50 µl per well).</i><i>2. The next day, the wells were washed three times with 100 µl of BPS Immuno Buffer 1 (BPS Bioscience, #79311). The plate was tapped upside down on paper towels to remove excess buffer.</i><i>3. The wells were blocked with 100 µl of Blocking Buffer 2 (BPS Bioscience, #79728) per well for 1 hour at room temperature with slow shaking.</i><i>4. Serial dilutions of anti-CD235a diluted in Blocking Buffer 2 were incubated in each well for 1 hour at room temperature (with slow shaking). (suggested range: 300 nM to 0 nM in replicates)</i><i>5. Wells were washed as in step 2 and incubated with Streptavidin-HRP for 1 hour at room temperature (with slow shaking), then washed again.</i><i>6. For detection, the wells were incubated with 50 µl of Colorimetric HRP Substrate (BPS Bioscience, #79651) for 1-10 minutes until a blue color developed in the positive control.</i><i>7. The reaction was then immediately quenched with an equal volume of 1N HCl and absorbance was measured at 450 nm.</i>

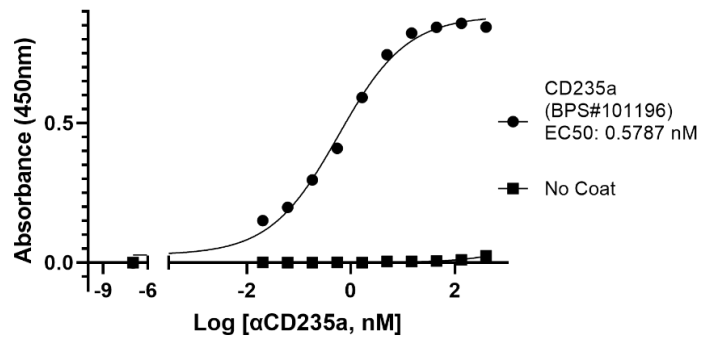
Applications: The anti-CD235a IgG1 antibody can be used for flow cytometry and immunofluorescence microscopy.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Binding assay of anti-CD235a and CD235 in ELISA



The plate was coated with CD235a protein (1 µg/ml) overnight. A "No Coat" control was included without CD235a protein. The binding assay was performed following the assay conditions described above, using a serial dilution of anti-CD235a. The absorbance of the "No Coat" control (blank) was subtracted from all other values.

Related Products

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
CD235a, Fc Fusion (IgG1), Avi-Tag HiP™	101196-1	100 µg
Blocking Buffer 2	79728	50 ml
Immuno Buffer 1	79311	50 ml