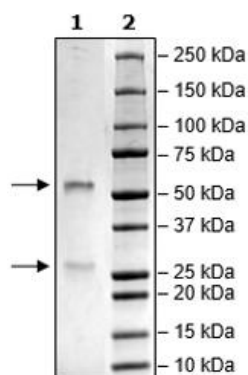


## Product Information

<b>Description:</b>	Biotinylated anti-CD123 IgG antibody is a purified recombinant antibody that recognizes human CD123 antigen. This antibody has been tested for specific binding affinity to purified human CD123 protein.
<b>Label:</b>	This antibody is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
<b>Concentration:</b>	1.41 mg/ml
<b>Species:</b>	Human
<b>Isotype:</b>	IgG1
<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>	Total: 150 kDa; HC: 51 kDa; LC: 24 kDa
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	Functional validation: The antibody was validated by measuring anti-CD123 binding to CD123 antigen in ELISA assay. The CD123 protein (BPS Bioscience #101035) was coated onto a 96-well plate overnight at 4°C (50 µl/well at a concentration of 4 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 anti-CD123 antibody (BPS Bioscience #101140), serially diluted in Blocking Buffer 2, was added for 30 minutes at room temperature. After 3 more washes, the plate was incubated with Streptavidin-HRP, 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



CD123: Anti-CD123-Biotin Binding Assay

