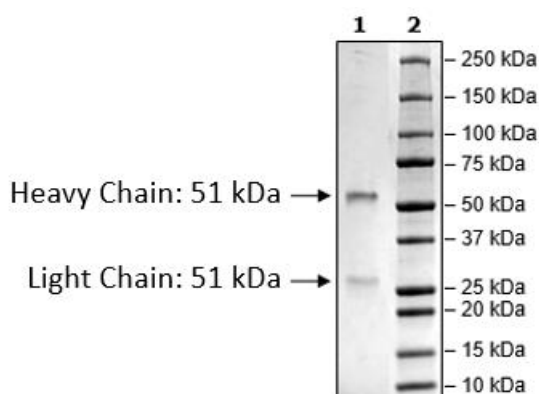


Product Information

Description:	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.
Label:	This antibody is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Concentration:	1.41 mg/ml
Species:	Human
Isotype:	IgG1
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:	Total: 150 kDa; HC: 51 kDa; LC: 24 kDa
Purity:	≥90%
Assay Conditions:	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.
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



CD123: anti-CD123-Biotin Binding Assay

