

## Product Information

<b>Description:</b>	Anti-CD123 IgG antibody is a purified recombinant antibody which recognizes CD123 antigen. This antibody has been tested for specific binding to purified human CD123 protein (BPS Bioscience, #101035) in a colorimetric ELISA binding assay.
<b>Concentration:</b>	0.50 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
<b>Expression System:</b>	Heavy chain (HC) and Light chain (LC) co-expressed in HEK293
<b>Purification:</b>	Protein A affinity purification from HEK293 cells
<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 (non-reduced); Heavy chain: 50 kDa; Light chain: 24 kDa
<b>Purity:</b>	≥90%

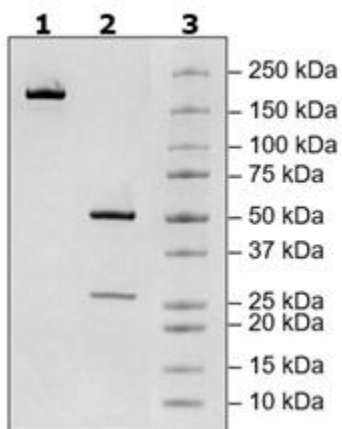
**Assay Conditions:** Experimental design and assay protocol for measuring anti-CD123 specific binding to CD123 antigen in ELISA assay:

1. Purified human His-tagged CD123 (BPS Bioscience, #101035) is bound to a clear 96-well nickel plate overnight at 4°C (1 µg/ml in PBS, 50 µl per well).
2. Wash each well 3x with Blocking Buffer 2, excess liquid is removed by tapping onto paper towels.
3. Wells are blocked by adding 100 µl of blocking buffer to each well. Incubate for 1 hour at room temperature with slow shaking.
4. Serial dilutions of anti-CD123 (200 nM to 0 nM in 3-fold dilutions, 50 µl per well) are incubated with bound CD56 for 1 hour at room temperature with slow shaking. Exclude wells designated "blank" and add 50 µl blocking buffer instead.
5. Next, wells are washed and incubated with 50 µl of anti-human IgG-HRP (1 µg/ml in Blocking Buffer 2) for 1 hour at room temperature with slow shaking. Then wells are washed again.
6. For detection, the wells are incubated with 100 µl Colorimetric HRP Substrate (BPS Bioscience #79651) for 1-10 minutes until a blue color develops in the positive control.
7. Quench the reaction using equal volume of 1N HCl. The blue color will change to yellow. Measure the absorbance at 450 nm.

**Applications:** This product is for research use only. It is not suitable for human diagnostic or therapeutic use. The anti-CD123 IgG format antibody can be used for flow cytometry and immunofluorescence microscopy.

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining



Lane 1: 5 µg Anti-CD123 (Non-reduced)  
Lane 2: 5 µg Anti-CD123 (Reduced)  
Lane 3: Protein Marker

### Binding assay of anti-CD123 and CD123 in ELISA

