

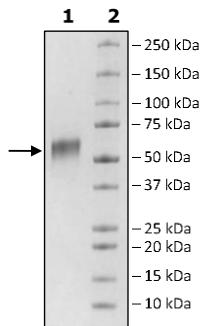
Product Information

Construct:	CD123 (20-305(end)-Avi-His)
Concentration:	0.43 mg/ml
Species:	Human
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM002183.4
MW:	36 kDa + glycans
Glycosylation:	This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.
Purity:	≥90%
Assay Conditions:	Experimental design and assay protocol for measuring anti-CD123 (BPS Bioscience, #101036) 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 (BPS Bioscience, #101036) (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 an equal volume of 1N HCl. The blue color will change to yellow. Measure the absorbance at 450 nm.
- Applications:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Binding assay of anti-CD123 and CD123 in ELISA

