

# CD123, Avi-Tag, His-Tag, Biotin-Labeled Recombinant

Catalog: 101817

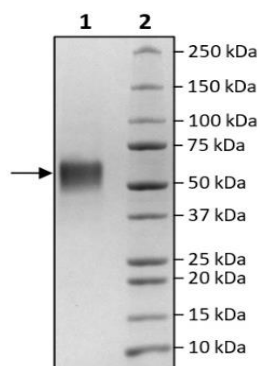
Lot: 240514

## Product Information

<b>Description:</b>	Recombinant human CD123 (also known as IL3Ra, interleukin 3 receptor alpha) encompassing amino acids 20-305 (end). This construct contains a C-terminal Avi-Tag™ followed by a C-terminal His-tag (6xHis). The recombinant protein was affinity purified.
<b>Background:</b>	CD123, also known as interleukin 3 receptor alpha (IL3Ra), is a cell surface receptor found in pluripotent progenitor cells and involved in hematopoietic lineage differentiation. It can also be found in basophils, plasmacytoid dendritic cells and dendritic cells. CD123 is overexpressed on leukemic hematopoietic stem cells (HSCs), but not normal HSCs. It has been found in acute myeloid and B-lymphoid leukemias, blastic plasmacytoid dendritic neoplasms (BPDCN), and hairy cell leukemia. This expression profile makes CD123 an attractive target in cancer therapy. CAR-Ts and other therapies targeting CD123 are being evaluated in clinical trials.
<b>Species:</b>	Human
<b>Construct:</b>	CD123 (20-305-Avi-His)-(Biotin)
<b>Concentration:</b>	0.35 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	36 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_002183.4
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Applications:</b>	Useful for SDS-PAGE and avidin pull-down assays.

## Quality Control Data

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



### Biotin-Avidin Pulldown

