

CD8a, Avi-Tag, His-Tag, Biotin-Labeled Recombinant

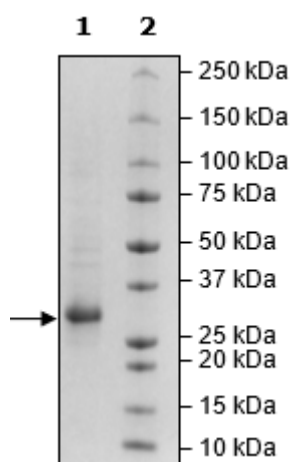
Catalog: 101873
Lot: 230627

Product Information

Description:	Recombinant human CD8a (cluster of differentiation 8 a) transcript 1 encompassing amino acids 22-182. This construct contains a C-terminal Avi-Tag™ followed by a His-Tag (6xHis). This protein was affinity purified.
Background:	CD8 is a co-receptor for the TCR (T cell receptor) in T cells, binding to MHC (major histocompatibility complex) class I proteins. CD8 is a typical marker of cytotoxic T cells and is involved in signaling. CD8 has two isoforms, a and b. CD8b recruits Lck (lymphocyte-specific protein tyrosine kinase) to the TCR-CD3 complex, and Lck phosphorylates multiple proteins involved in activation of cytotoxic T-lymphocytes. It is thus critical for the lysis of cancer cells.
Species:	Human
Construct:	CD8a (22-182-Avi-His)-(Biotin)
Concentration:	0.50 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	21 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_001768.7
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation was confirmed to be ≥90%.
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	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.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

