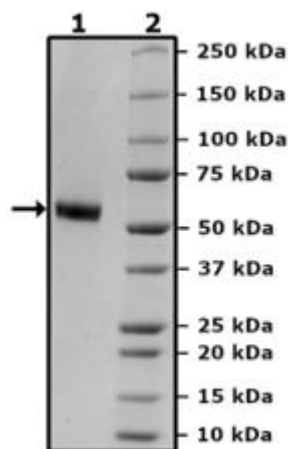


## Product Information

<b>Description:</b>	Recombinant human CD8a (cluster of differentiation 8 a) transcript 1 encompassing amino acids 22-182. This construct contains a C-terminal Fc domain followed by an Avi-Tag™. The recombinant protein was affinity purified. HiP™ indicates a high purity protein (≥90% pure) with less than 10% aggregation as measured by gel filtration.
<b>Background:</b>	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.
<b>Species:</b>	Human
<b>Construct:</b>	CD8a (22-182-Fc(IgG1)-Avi)
<b>Concentration:</b>	2.08 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>	47 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Aggregation:</b>	<10%
<b>Genbank Accession:</b>	NM_001768.7
<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.

## Quality Control Data

## 4-20% SDS-PAGE Coomassie Staining



## Quality Control Data

### Gel Filtration Curve

