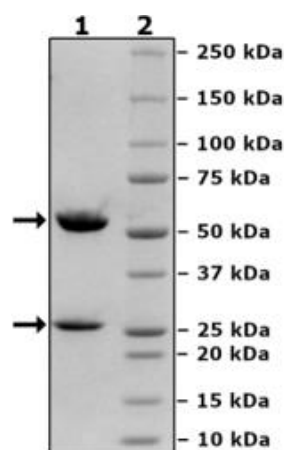


Product Information

Description:	This purified recombinant antibody is a humanized mouse anti-CD8 a (cluster of differentiation 8 a) antibody. This antibody has a C-terminal Avi-Tag™ on its heavy chain. This antibody recognized human CD8.
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 CTLs (cytotoxic T-lymphocytes). It is thus critical for the lysis of cancer cells. Antibodies against CD8 are currently used in the clinic for the diagnostic of cancer, as in the case of crefmirlimab.
Species:	Mouse
Clonality:	Monoclonal
Concentration:	1.76 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:	Heavy Chain: 51 kDa; Light Chain: 23 kDa + glycans
Glycosylation:	This antibody runs at a higher MW by SDS-PAGE due to glycosylation.
Label:	This antibody 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

