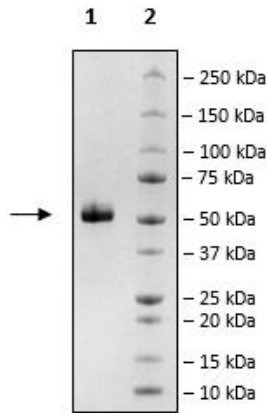


Product Information

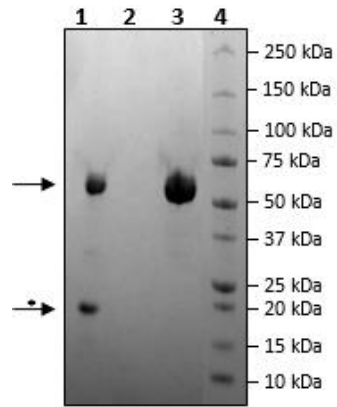
Description:	Recombinant human CD4, encompassing amino acids 26-396. This construct contains a C-terminal Avi-Tag™ followed by an His-tag (6xHis). This protein was affinity purified.
Background:	CD4 (cluster of differentiation 4) is part of the immunoglobulin superfamily, and it can be found in T-helper cells, monocytes, macrophages, and dendritic cells. It is a co-receptor in the TCR (T-cell receptor), binding to MHC (major histocompatibility complex) class II molecules. CD4 binds to the tyrosine kinase Lck (lymphocyte-specific protein tyrosine kinase), which phosphorylates the ITAM (immunoreceptor tyrosine activation motifs) domain of the CD3, activating CD3 related signaling. CD4 is a typical T cell marker of T helper cells, and it has been linked to cancer, autoimmune diseases such as vitiligo and type I diabetes. In addition, HIV-1 makes use of CD4 to trigger viral envelope protein conformational changes that allow cell infection. Ibalizumab, an anti-CD4 antibody, is currently used in the treatment of HIV, being considered a first-in-class medication. Further studies into CD4 and potential strategies around it may benefit patients and CD4 related diseases.
Species:	Human
Construct:	CD4 (26-396-Avi-His)-(Biotin)
Concentration:	0.62 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:	45 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_000616.5
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is 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.
Applications:	Useful for avidin binding studies.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown



1. Beads
2. Flow Thru
3. Control
4. Standards

* Avidin from beads.