

Product Information

Description:	Recombinant human IL-2RB (Interleukin-2 receptor subunit beta), encompassing amino acids 27-239. This construct contains a C-terminal Avi-Tag™ followed by a His-tag (6xHis). The recombinant protein was affinity purified.
Background:	Interleukin-2 (IL-2) is a cytokine secreted by activated T lymphocytes to stimulate the proliferation of both B and T cells. IL-2 is involved in the extreme inflammation response termed “cytokine storm” observed following viral infection in susceptible patients or following adaptive cell therapy. Thus, it is considered a therapeutic target for the treatment of severe COVID-19. IL-2 acts by binding to a membrane heterotrimeric receptor consisting of IL-2 receptor alpha (IL2RA, also known as CD25) and beta (IL2RB, also known as CD122) chains, which together with the common gamma chain (IL2RG) form a high-affinity receptor, whereas IL-2RA homodimers form a low-affinity receptor. Polymorphisms in the genes coding for IL2RA and IL2RB may be associated with risk of lung cancer.
Species:	Human
Construct:	IL2RB (27-239-Avi-His)
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:	28 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_000878
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 SDS-PAGE.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

