

Product Information

Description:	Recombinant KIR3DL3 (Killer Cell Immunoglobulin-like Receptor 3DL3) encompassing amino acids 26-322. This construct contains a C-terminal Avi-Tag™, followed by a His-Tag (6xHis). The recombinant protein was affinity purified.
Background:	KIR3DL3, also known as Killer Cell Immunoglobulin-like Receptor 3DL3, is an inhibitory receptor present in NK cells. It is a framework gene, and present in all human haplotypes. In healthy individuals it is expressed at very low levels in CD8 ⁺ T _{EMRA} and CD56 ^{dim} NK cells, but in certain pathologies its levels are increased. It was recently found that B7-H7 is a ligand for KIR3DL3, an interaction that leads to NK cell suppression. Its role as immune checkpoint makes it an attractive target in cancer therapy.
Species:	Human
Construct:	KIR3DL3 (26-322-Avi-His)
Concentration:	0.51 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:	36 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_153443.5
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

