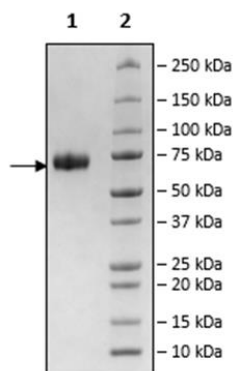


Product Information

Description:	Recombinant human KIR2DL3 (Killer cell immunoglobulin-like receptor 2DL3), encompassing amino acids 22-245. This construct contains at the C-terminus the Fc domain of a human IgG1, followed by an Avi-Tag™. This protein was affinity purified. This protein is enzymatically biotinylated using Avi-Tag™ technology.
Background:	KIR2DL3 (Killer cell immunoglobulin-like receptor 2DL3), also known as CD158b, regulates the immune response by binding MHC class I molecules and inhibiting natural killer (NK) cells. This protein belongs to the KIR (killer cell immunoglobulin-like receptors) family, which are immune-regulating glycoproteins expressed on NK cells and some T cells. KIR2DL3 is associated with graft-versus-host disease and prostate cancer. Further research will highlight the role of KIR2DL3 in immune response pathways.
Species:	Human
Construct:	KIR2DL3 (22-245-Fc(IgG1)-Avi)-(Biotin)
Concentration:	1.2 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:	53 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Genbank Accession:	NM_015868
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 and avidin pulldown assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

