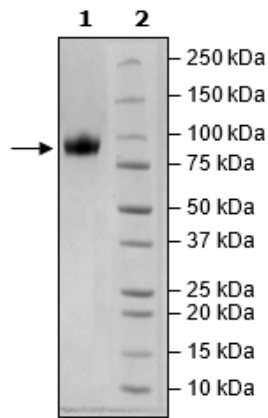


Product Information

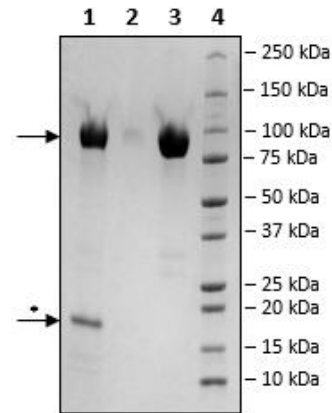
Description:	Recombinant human CD155, encompassing amino acids 27-343. This construct contains a C-terminal human IgG1 Fc domain followed by an Avi-Tag™. This protein was affinity purified.
Background:	CD155 (also known as PVR cell adhesion molecule) is a transmembrane glycoprotein receptor of the immunoglobulin superfamily of proteins. It is highly expressed in dendritic cells, fibroblasts, endothelial cells, and multiple tumor cells including ovarian carcinoma, non-small cell lung, glioblastoma, and colorectal carcinoma. It has a high affinity for a T cell regulatory transmembrane surface protein called TIGIT (T cell immunoglobulin and ITIM domains) and plays a key role in tumor cell invasion and migration. Interaction between CD155 and TIGIT causes immunosuppressive effects on CD4 ⁺ and CD8 ⁺ T cells, which hinders T cell proliferation and function. CD155 also interacts with the regulatory receptors CD226 (expressed on natural killer (NK) cells, monocytes and CD4 ⁺ T cells) and CD96. While CD155-CD226 engagement activates NK cell cytotoxicity and T cell response, CD155-CD96 interaction inhibits NK cell function. Due to these opposing regulatory effects, blockade of CD155-TIGIT and CD155-CD226 interactions are attractive therapeutic strategies in cancer immunotherapy and autoimmune diseases, respectively.
Species:	Human
Construct:	CD155 (27-343-Fc(IgG1)-Avi)-(Biotin)
Concentration:	1.40 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:	63 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_006505
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 pull down assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown



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

* Avidin from beads.