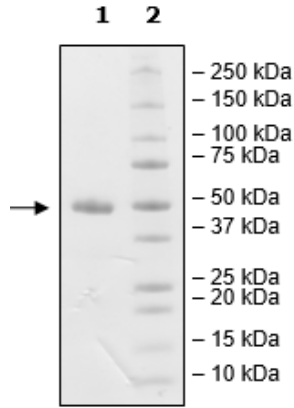


Product Information

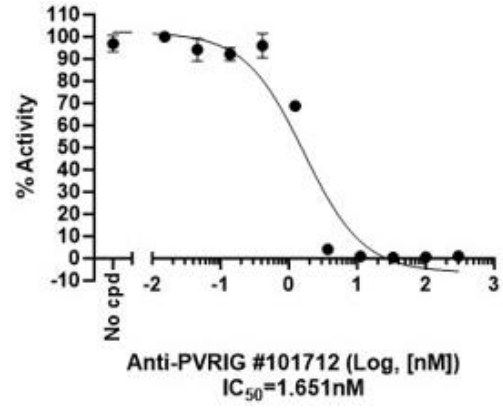
Description:	Human secreted CD112, encompassing amino acids 32-360. This construct has a C-terminal His-tag followed by an Avi-Tag™. This protein was enzymatically biotinylated using Avi-tag™ technology and was affinity purified. HiP™ indicates a high purity protein with ≥90% purity as measured by gel filtration.
Background:	CD112 (Poliovirus receptor-related 2, PVRL2, or nectin-2), is widely expressed on antigen presenting cells and tumor cells and is the high affinity ligand of CD112R (Poliovirus receptor related immunoglobulin domain containing, PVRIG). CD112R is found in NK and T cells, and in T cells it is found particularly in CD8 ⁺ T cells. High levels of this protein are found in NK, CD8 ⁺ and CD4 ⁺ T cells of patients with kidney, ovary, lung, prostate and endometrium cancer and acute myeloid leukemia (AML). CD112 has also been linked to tumor angiogenesis, growth and metastasis. It has a wider pattern of expression, being also present in endothelial cells, neurons and fibroblasts. CD112-CD112R interaction is a positive immune checkpoint that enhances human T cell response and has emerged as an attractive therapeutic target for oncology. TIGIT (T cell immunoreceptor with Ig and ITIM domains) and CD226 (also called DNAM-1, DNAX Accessory Molecule-1) also bind CD112. The development of inhibitors able to prevent the interaction between CD112R-CD112 alone or in combination with blockage of TIGIT-CD155/CD112 interaction may prove beneficial in cancer therapy.
Species:	Human
Construct:	CD112 (32-360-His-Avi)-(Biotin)
Concentration:	0.58 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:	38 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Aggregation:	<10%
Genbank Accession:	NM_002856
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation 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.
Assay Conditions:	Assay run according to CD112R:CD112 [Biotinylated] Inhibitor Screening Assay Kit (BPS Bioscience #79732). 100 ng CD112R were used to coat plates overnight at 4°C. Various amounts of CD112 [Biotinylated] incubated for 2 hours at room temperature with the coated plates.
Applications:	Useful for neutralizing of CD112 in ELISA and in cellular assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PVRIG:PVRL2-Biotin Neutralizing Activity



Biotin-Avidin Pulldown



Quality Control Data

Gel Filtration Curve

