

## Product Information

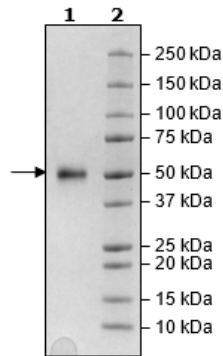
<b>Description:</b>	Recombinant human CD112R, also known as PVRIG (Poliovirus receptor-related Immunoglobulin domain-containing protein), encompassing amino acids 41-172. This construct contains a C-terminal Human IgG1, Fc region followed by an Avi-tag™. This recombinant protein was affinity purified.
<b>Background:</b>	PVRIG, also known as Poliovirus receptor related immunoglobulin domain containing protein or CD112R, is a natural killer (NK) cell immune checkpoint receptor, and it is also found in CD8 <sup>+</sup> T cells. PVRIG works as an inhibitory receptor and recognizes CD112 (also known as poliovirus receptor-related 2, PVRL2 or nectin-2) as its ligand. This binding results in lower responses of NK and T cells towards cancer cells. In T cells it inhibits the nuclear factor of activated T cells (NFAT). Blocking the interaction between PVRIG and its ligand, for instance with an antibody directed towards PVRIG, represents a novel strategy in checkpoint blockade immunotherapy.
<b>Species:</b>	Human
<b>Construct:</b>	CD112R (41-172-Fc(IgG1)-Avi)
<b>Concentration:</b>	0.33 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	42 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Aggregation:</b>	<10%
<b>Genbank Accession:</b>	NM_024070
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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.
<b>Assay Conditions:</b>	Assay run according to CD112R:CD112 [Biotinylated] Inhibitor Screening Assay Kit (BPS Bioscience #79732). 100 ng of CD112R was used to coat plates overnight at 4°C. Various amounts of CD112 [Biotinylated] were incubated for 2 hours at room temperature.
<b>Applications:</b>	Useful for binding and neutralizing of PVRIG in ELISA and in cellular assays.

# CD112R (PVRIG), Fc-Fusion, Avi-Tag (Human) HiP™ Recombinant

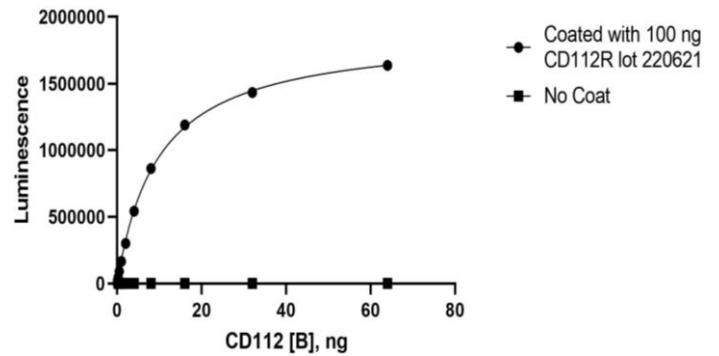
Catalog: 79116  
Lot: 220621

## Quality Control Data

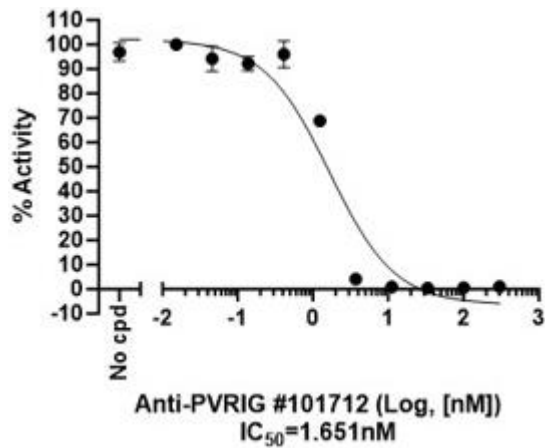
### 4-20% SDS-PAGE Coomassie Staining



### CD112R:CD112 [Biotinylated] Binding Activity



### PVRIG:PVRL2-Biotin Neutralizing Assay



Quality Control Data

Gel Filtration Curve

