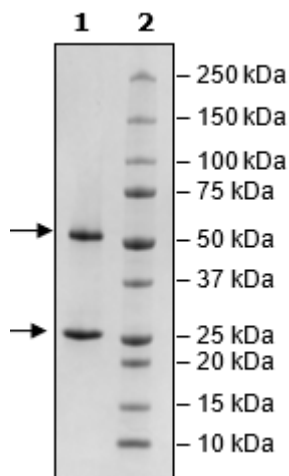


Product Information

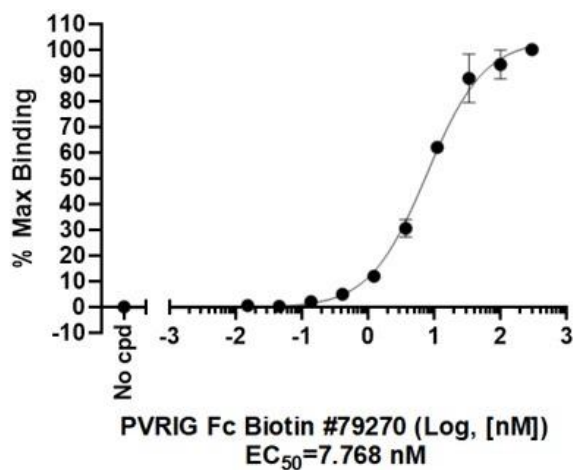
Description:	This anti-PVRIG (Poliovirus receptor related immunoglobulin domain containing) antibody is a purified, recombinant human monoclonal antibody which recognizes human PVRIG. This recombinant antibody has been tested for specific binding to purified recombinant human PVRIG in an ELISA binding assay.
Background:	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 ⁺ 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.
Species:	Human
Isotype:	IgG1
Clonality:	Monoclonal
Concentration:	1.47 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:	Heavy Chain: 50 kDa; Light Chain: 24 kDa
Stability:	At least 12 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:	The antibody was validated by measuring its binding to a PVRIG antigen in ELISA. PVRIG protein (BPS Bioscience #101712) was coated onto a 96-well plate overnight at 4°C (50 µl/well at a concentration of 2 µg/ml in PBS). The plate was washed 3 times with Immuno Buffer 1 (BPS Bioscience #79311) and blocked using 100 µl of Blocking Buffer 2 (BPS Bioscience #79728) for 1 hour at Room Temperature (RT). After removing the blocking buffer, 50 µl/well of serially diluted CD112R (PVRIG), Fc-Fusion, Avi-Tag, Biotin-Labeled (Human) Recombinant (BPS Bioscience #79270) in Blocking Buffer 2, was added for 1 hour at RT. After 3 washes, the plate was incubated with Streptavidin-HRP (BPS Bioscience #79742), washed again, and colorimetric HRP substrate was added. The reaction was stopped, and absorbance was read at λ=450 nm. The Blank value was subtracted from all values.
Applications:	Useful for binding and neutralization of PVRIG in ELISA and in cellular assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Anti-PVRIG: PVRIG Biotin Binding Assay



PVRIG: PVRL2-Biotin Neutralizing Assay

