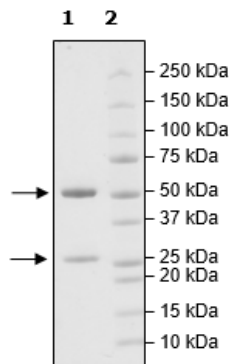


Product Information

Description:	Recombinant monoclonal antibody recognizing human CD28. This antibody binds to CD28 on T cells and promotes T cell response.
Background:	The T cell-specific receptor CD28 is an important regulator of the immune system. The ligation of CD28 and its ligands CD80/CD86 on the APC provides a critical costimulatory signal alongside T cell receptor (TCR) for T cell activation. The costimulatory signal can also be delivered by the interaction between CD28 and its agonist antibody.
Concentration:	1.05 mg/ml
Host Species:	Human
Isotype:	IgG1
Clonality:	Monoclonal
Cross Reactivity:	This antibody has not been tested for cross reactivity with any other species.
Formulated In:	8 mM Na Phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl and 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	Heavy Chain: 51 kDa, Light Chain: 27 kDa
Purity:	≥90%
Purification:	Protein A affinity chromatography from HEK293 supernatants.
Assay Conditions:	Serial dilutions of anti-CD3 (BPS Bioscience, #71274) and CD28 antibodies were added to the IL-2 reporter-Jurkat cells, then incubated at 37°C incubator for 16 hrs. After the treatment, perform Luciferase assay using One-Step Luciferase assay system (BPS Bioscience, #60690)
Applications:	This antibody is useful for activation of CD28 costimulatory pathway in T cells.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Costimulatory Activity of Anti-CD28 Antibody on IL-2 reporter/Jurkat Cells

