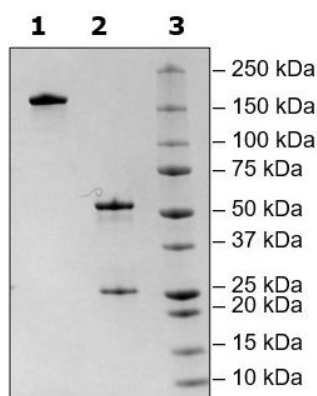


Product Information

Description:	Anti-PD-1 IgG antibody is a purified recombinant antibody which recognizes human PD1 antigen. This antibody has been functionally tested in two co-culture assays.
Concentration:	0.50 mg/ml
Species:	Human
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	Heavy chain (HC) and Light chain (LC) co-expressed in HEK293
Purification:	Protein A affinity purification from HEK293 cells
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	150 kDa (non-reduced); Heavy chain: 50 kDa; Light chain: 24 kDa
Purity:	≥90%
Assay Conditions:	The antibody acted as a PD-1 agonist when added to FcGR2B/TCR activator CHO cells co-cultured with PD-1/NFAT Jurkat cells, as indicated by PD-1-mediated inhibition of TCR activation (co-culture assay described in BPS Bioscience #78436). In addition, this anti-PD-1 antibody did not antagonize PD-1/PD-L1 binding since the TCR activation signal was still inhibited by PD-1/PD-L1 signaling in the presence of the antibody (co-culture assay described in BPS Bioscience #60535).
Applications:	This product is for research use only. It is not suitable for human diagnostic or therapeutic use.

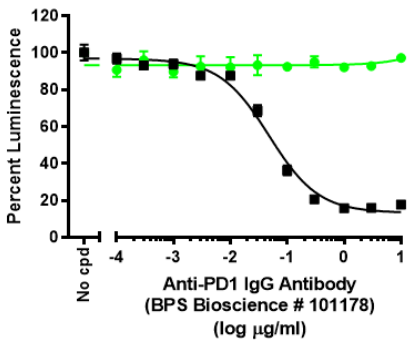
Quality Control Data

4-20% SDS-Page Coomassie Staining



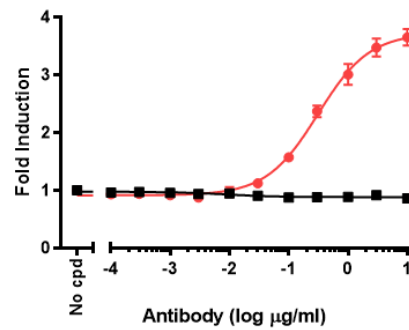
Lane 1: 5 µg Anti-PD1 (Non-reduced)
Lane 2: 5 µg Anti-PD1 (Reduced)
Lane 3: Protein Marker

Agonistic Effect in Co-Culture Assay



- FcGR2B/TCR Activator-CHO Recombinant Cell Line (BPS Bioscience # 78436)
EC₅₀ = 0.05 µg/ml
- TCR Activator-CHO Recombinant Cell Line (BPS Bioscience # 60539)
EC₅₀ = not determined

Lack of Effect on PD-1/PD-L1 Interaction



- Anti-PD1 IgG Antibody (BPS Bioscience # 101178)
EC₅₀ = not determined
- Anti-PD1 Neutralizing Antibody (BPS Bioscience # 71120)
EC₅₀ = 0.31 µg/ml