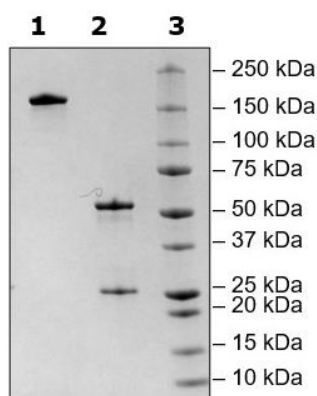


## Product Information

<b>Description:</b>	Anti-PD-1 agonist antibody is a purified recombinant antibody which recognizes human PD1 antigen. This antibody has been functionally tested in two co-culture assays.
<b>Concentration:</b>	0.98 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>Expression System:</b>	Heavy chain (HC) and Light chain (LC) co-expressed in HEK293
<b>Purification:</b>	Protein A affinity purification from HEK293 cells
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 12 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>MW:</b>	150 kDa (non-reduced); Heavy chain: 50 kDa; Light chain: 24 kDa
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	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).
<b>Applications:</b>	This product is for research use only. It is not suitable for human diagnostic or therapeutic use.

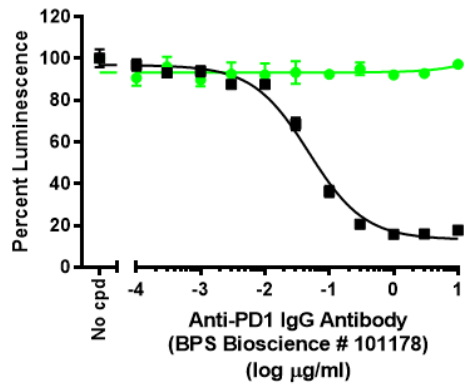
## Quality Control Data

## 4-20% SDS-PAGE Coomassie Staining



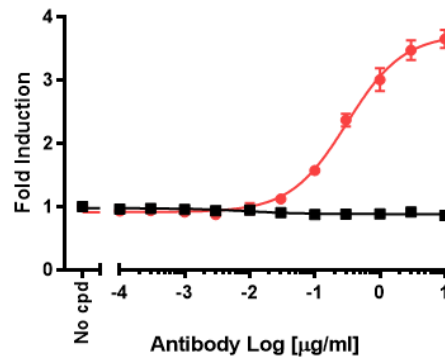
Lane 1: 5 µg Anti-PD1 (Non-reduced)  
Lane 2: 5 µg Anti-PD1 (Reduced)  
Lane 3: Protein Marker

## Agonist Effect in Co-Culture Assay



- FcGR2B/TCR Activator-CHO Recombinant Cell Line (BPS Bioscience # 78436)  
EC<sub>50</sub> = 0.05 µg/ml
- TCR Activator-CHO Recombinant Cell Line (BPS Bioscience # 60539)  
EC<sub>50</sub> = not determined

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



- Anti-PD-1 Agonistic Antibody (BPS Bioscience # 101178)  
EC<sub>50</sub> = not determined
- Anti-PD1 Neutralizing Antibody (BPS Bioscience # 71120)  
EC<sub>50</sub> = 0.31 µg/ml