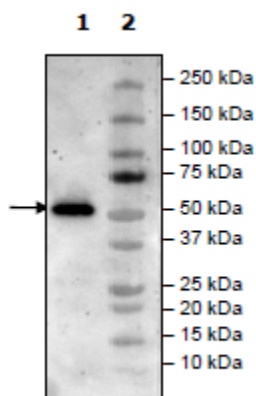


## Product Information

<b>Description:</b>	This antibody binds and blocks the activation of the PD-1 receptor. The antibody is humanized, containing a human IgG4 Fc region with an engineered hinge region mutation (S228P) designed to prevent the exchange of IgG4 molecules.
<b>Concentration:</b>	2.22 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>	HEK293
<b>Clonality:</b>	Monoclonal
<b>Isotype:</b>	IgG4
<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>	Heavy Chain: 50 kDa, Light Chain: 25 kDa
<b>Purity:</b>	≥90%
<b>Purification:</b>	Protein A affinity chromatography from HEK293 supernatants.
<b>Cross reactivity:</b>	This antibody has not been tested with other species
<b>Assay Conditions:</b>	Reaction done according to PD-1 [Biotinylated]: PD-L1 Inhibitor Screening Assay Kit (BPS Bioscience, #72005) protocol. Buffer is PBS + 0.1% BSA. Blocking buffer is Superblock + 0.05% Tween-20. Coat protein was added ON at 4 deg 100 ng/well. Various concentrations of PD-1 neutralizing antibody (0.02-300 nm) added for 2 hours at room temperature. Strep-HRP added for 1 hour at room temperature.
<b>Applications:</b>	Useful for inhibition of PD-1:PD-L1 and PD-1:PD-L2 interactions and for western blot analysis.

## Quality Control Data

### Western Blot



### Inhibition of PD-1:PD-L1 by Neutralizing PD-1 (S228P) Antibody

