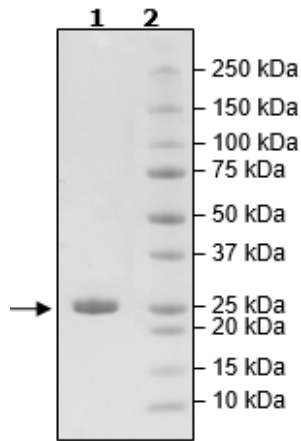


Product Information

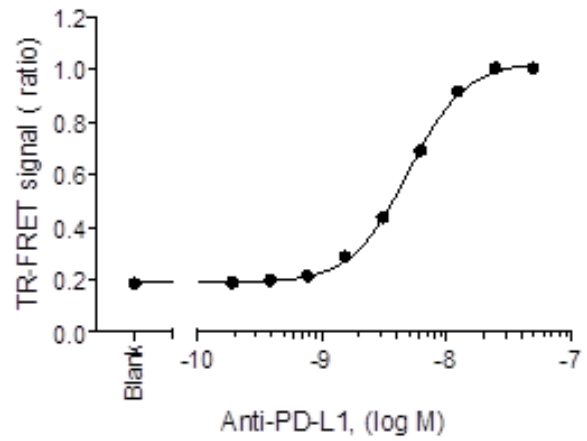
Description:	Recombinant humanized monoclonal anti-PD-L1 (Programmed death-ligand 1) antibody Fab (fragment antigen-binding) able to recognize the extracellular domain of human PD-L1 and block the binding of PD-L1 to PD-1. This antibody has an Avi-Tag™ at the N-terminus of the heavy chain.
Species:	Human
Isotype:	IgG1
Clonality:	Monoclonal
Concentration:	1.64 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Purification:	Protein A affinity chromatography from HEK293 supernatants.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	Heavy Chain: 25 kDa; Light Chain: 23 kDa
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation was confirmed to be ≥90%.
Stability:	At least 12 months at -80°C. Protect from light.
Storage:	-80°C
Cross Reactivity:	This antibody Fab has been shown to cross-react with mouse PD-L1. No other species have been tested.
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:	Anti-PD-L1-Biotin (BPS Bioscience #11032) was titrated from 0 to 50 nM in a 2-fold dilution series in presence of 100 nM PD-L1-FLAG-tag (BPS Bioscience #71183) and incubated for 1 hour at RT. Binding was detected with Anti-FLAG-Eu and Streptavidin-APC.
Applications:	Useful as an antagonist to block PD-L1 signaling and binding to PD-1 in biochemical assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Anti-PD-L1: PD-L1 TR-FRET Binding Assay



Biotin-Avidin Pulldown

