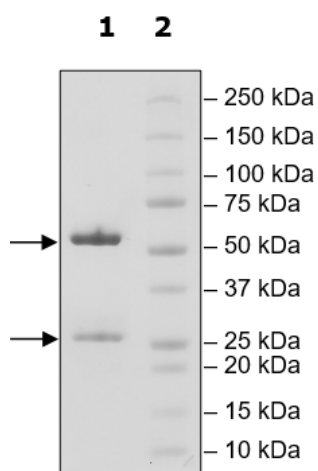


Product Information

Description:	Purified, recombinant, humanized anti-PD-L1 (IgG1) antibody with C-terminal Avi-tag, expressed in HEK293 cells. Purified antibody was enzymatically biotinylated using Avitag™ technology. This antibody reacts with human PD-L1 and cross-reacts with mouse PD-L1. It has not been tested for cross-reactivity with other species.
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Species	Human
Isotype:	IgG1
Clonality:	Monoclonal
Concentration:	0.60 mg/ml
Expression System:	HEK293
Purity:	≥90%
Purification:	Protein A affinity chromatography from HEK293 supernatants
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, 110 mM NaCl, 2.2 mM KCl, pH 7.4, and 20% glycerol
MW:	Heavy Chain: 53 kDa; Light Chain: 25 kDa
Stability:	At least 12 months at -80°C.
Storage:	-80°C
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.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

