

# PD-L1 (CD274), Fc Fusion, Biotin-Labeled (Human) HiP™ Recombinant

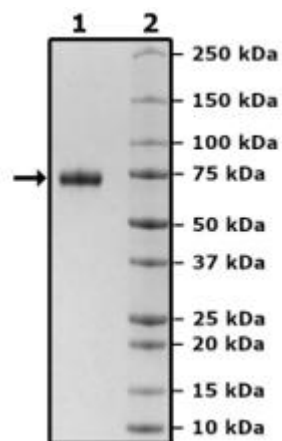
Catalog: 71105  
Lot: 230801-G

## Product Information

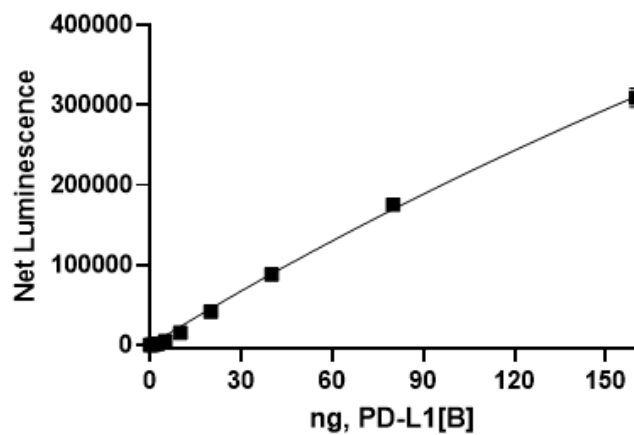
<b>Description:</b>	Human secreted Programmed Cell Death 1 Ligand 1 (PD-L1)-Fc fusion protein, encompassing amino acids 19-239, fused at the C-terminus with the Fc portion of human IgG1 followed by an Avi-tag™. HiP™ indicates a high purity protein (≥90% pure) and less than 10% aggregation as measured by gel filtration.
<b>Background:</b>	Programmed Cell Death 1 Ligand 1 (PD-L1), also known as CD274 and B7 homolog 1 (B7-H1), is an immune checkpoint acting as a negative regulator of immune cells. It helps maintain immune homeostasis and prevent excessive immune responses. Normally expressed at the surface of antigen-presenting cells, it is also found in tumor cells and contributes to cancer-associated immune suppression through its interaction with PD-1 expressed in T cells. Checkpoint inhibitors targeting the interaction between PD-L1 and PD-1 (such as pembrolizumab and nivolumab) have demonstrated significant clinical activity in various types of cancer and proved transformative for cancer therapy.
<b>Species:</b>	Human
<b>Construct:</b>	PD-L1 (19-239-Fc(IgG1)-Avi)-(Biotin)
<b>Concentration:</b>	1.16 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	54 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Aggregation:</b>	<10%
<b>Genbank Accession:</b>	NM_014143
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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.
<b>Assay Conditions:</b>	Assay was performed according to PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit (BPS Bioscience #72003) with PD-L1 (CD274), Fc Fusion, Biotin-labeled (Human) Recombinant titrated at various concentrations.
<b>Applications:</b>	Useful for binding studies and screening inhibitors of PD1 binding to PD-L1.

Quality Control Data

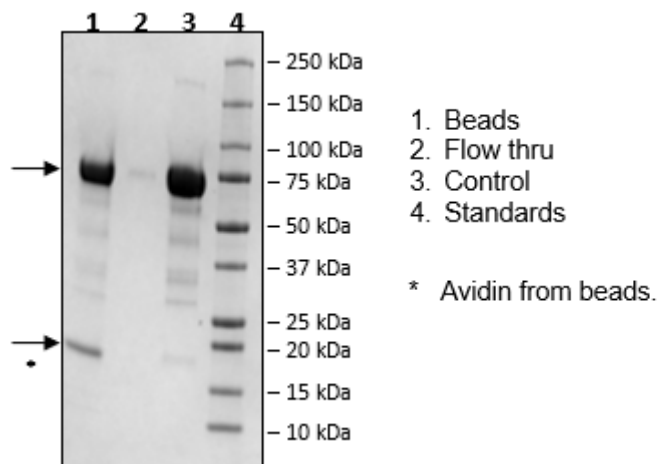
4-20% SDS-PAGE Coomassie Staining



PD-1:PD-L1[Biotinylated] Binding



Biotin-Avidin Pulldown



Quality Control Data

Gel Filtration Curve

