

Data Sheet

PD-L1 (CD274), Fc fusion, Biotin-labeled

Human, Recombinant, Fc fusion protein
 with C-terminal Avi-tag™

Catalog #: 71105-2

Lot #: 170323

Conc.: 0.99 mg/ml

Formulated in: 8 mM phosphate, pH 7.4,
 110 mM NaCl, 2.2 mM KCl, and 20%
 glycerol

Stability: At least 6 months at -80°C. Avoid
 freeze/thaw cycles. Protein may be diluted
 to ≥ 100 µg/ml in PBS + glycerol and stored
 at -80°C.

References:

1. Freeman, G.J., *et al.*, *J Exp Med.* 2000
 Oct 2;**192(7)**:1027-34.
2. Dong, H., *et al.*, *Nat Med.* 1999
 Dec;**5(12)**:1365-9.

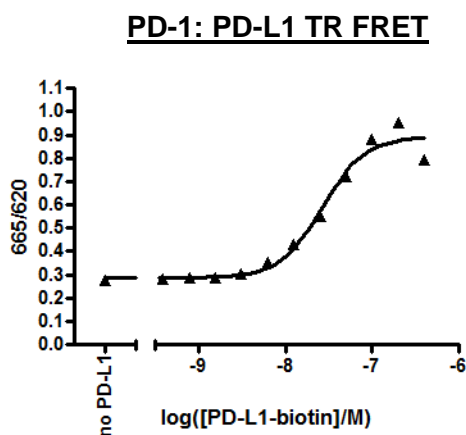
Description: Human secreted Programmed
 Cell Death 1 Ligand 1 (PD-L1)-Fc fusion
 protein, also known as CD274 and B7
 homolog 1 (B7-H1), GenBank Accession
 No. NM_014143, amino acids 19-239, fused
 at the C-terminus with the Fc portion of
 human IgG1, with C-terminal Avi-tag™
 expressed in a HEK293 cell expression
 system and enzymatically biotin-labeled
 using Avi-tag™ technology. MW = 54 kDa.

Assay Conditions: PD-1-PD-L1 binding
 activity, measured using the PD-1:PD-
 L1[Biotinylated] Inhibitor Screening Assay
 Kit, BPS Bioscience, Catalog #72003.

Application:

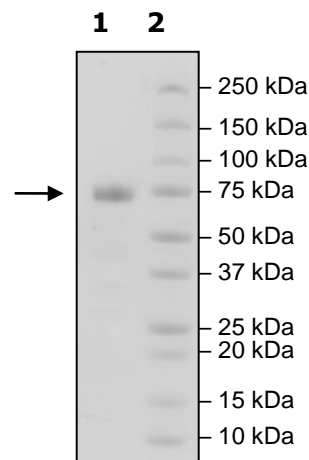
Useful for the study of protein-protein
 interaction and *in vitro* and *in vivo* protein
 function.

Quality Assurance



**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 4µg PD-L1
Lane 2:
 Protein Marker
MW: 54kDa+glycans
Purity: ≥90%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com