

Data Sheet

PD-L1 (I54A), Fc fusion, Biotin-labeled

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

Catalog #: 71314

Lot #: 150722

Conc.: 1.30 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) with a I54A mutation, 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.

Application:

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

Quality Assurance

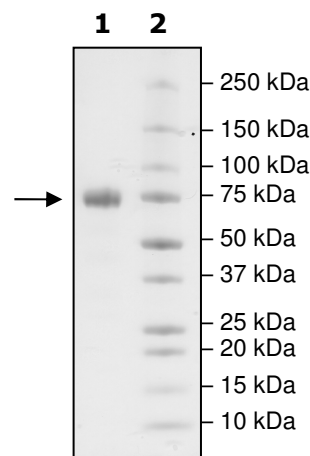
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
3 µg PD-L1

Lane 2:
Protein Marker

MW: 54kDa+glycans

Purity: ≥90%



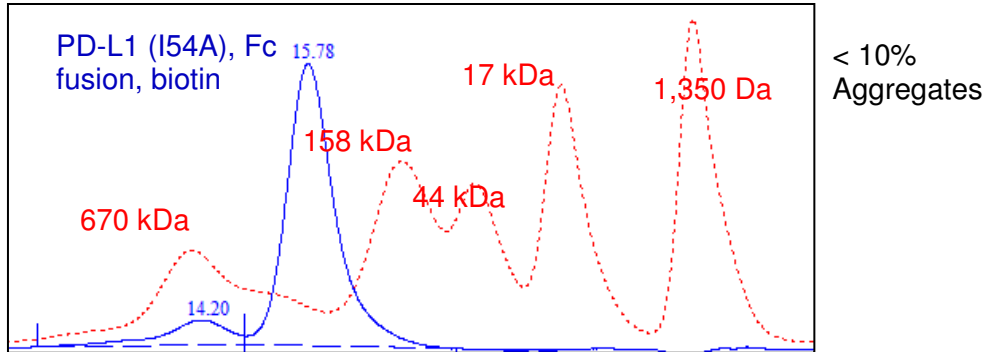
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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

Gel Filtration Curve



Sequence:

FTVTVPKDLYVVEYGSNMTIECKFPVEKQLDLAALAVQWEMEDKNI IQFVHGEEEDLKVQHSSYRQ
 RARLLKDQLSLGNAALQITDVKLQDAGVYRCMISYGGADYKRITVKVNAPYNKINQRILVVDPVT
 SEHELTCQAEQYPAEVIWTSSDHQVLSGKTTTTNSKREEKLFNVTSTLRINTTTNEIFYCTFRR
 LDPEENHTAELVIPELPLAHPNERTDIEGRMDPKSCDKTHTCPPCPAPELLGGPSVFLFPPKPK
 DTLMI SRTPEVTCVVVDVSHEDPEVKFNWYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLHQDW
 LNGKEYKCKVSNKALPAPIEKTI SKAKGQPREPQVYTLPPSREEMTKNQVSLTCLVKGFYPSDIA
 VEWESNGQPENNYKTPPVLDSDGSFFLYSKLTVDKSRWQQGNVFSQVMSHEALHNHYTQKLSLSL
 SPGKGGGLNDIFEAQKIEWHE

Black: PD-L1 (a.a. 19-239)

Green: Linker

Red: Fc of human IgG1 (a.a.100-330)

Blue: Avi-tag™

Purple: I54A mutation

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