

## Data Sheet

### PD-L1 (A121Q), Fc fusion, Biotin-labeled

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

**Catalog #:** 71320

**Lot #:** 150722

**Conc.:** 1.44 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  $-80^{\circ}\text{C}$ . Avoid  
freeze/thaw cycles. Protein may be diluted  
to  $\geq 100 \mu\text{g/ml}$  in PBS + glycerol and stored  
at  $-80^{\circ}\text{C}$ .

**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 A121Q 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:**

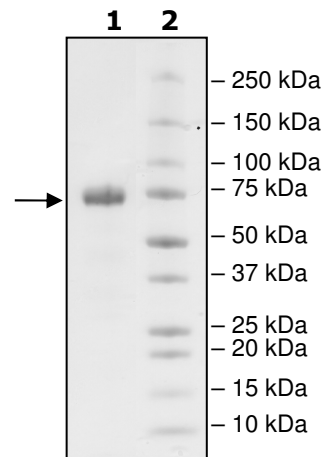
3  $\mu\text{g}$  PD-L1

**Lane 2:**

Protein Marker

**MW:** 54kDa+glycans

**Purity:**  $\geq 90\%$



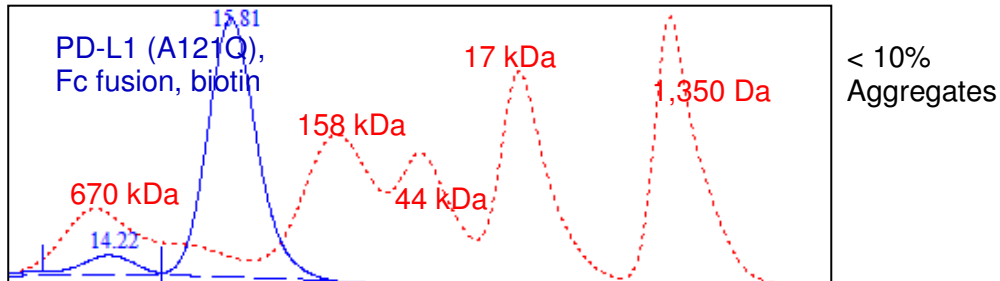
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

**Gel Filtration Curve**



**Sequence:**

FTVTVPKDL YVVEYGSNMTIECKFPVEKQLDLAALIVQWEMEDKNI IQFVHG EEDLKVQHSSYRQ  
 RARLLK DQLSLGNAALQITDVKLQDAGVYRCMISYGGQDYKRITVKVNAPYNKINQRILVVD PVT  
 SEHELTCQAEGYPKAEVIWTSSDHQVLSGKTTTTNSKREEKLFNVTSTLRINTTNEIFYCTFRR  
 LDPEENHTAELVIPELPLAHPNERTDIEGRMDPKSCDKTHTCPPCPAPELLGGPSVFLFPPKPK  
 DTLMI SRTPEVTCVVVDVSHEDPEVKFNWYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLHQDW  
 LNGKEYKCKVSNKALPAPIEKTI SKAKGQPREPQVYTLPPSREEMTKNQVSLTCLVKGFYPSDIA  
 VEWESNGQPENNYKTPPVLDSDGSFFLYSKLTVDKSRWQQGNV FSCSVMHEALHNHYTQKLSLSL  
 SPGKGGGLNDIFEAQKIEWHE

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

Green: Linker

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

Blue: Avi-tag™

Purple: A121Q 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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)