

## Data Sheet

### **PD-L1 (D26A), Fc fusion, Biotin-labeled**

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

**Catalog #:** Custom

**Lot #:** 150803-G

**Conc.:** 0.54 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 D26A 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 and *in vitro* and *in vivo* protein  
function.

## **Quality Assurance**

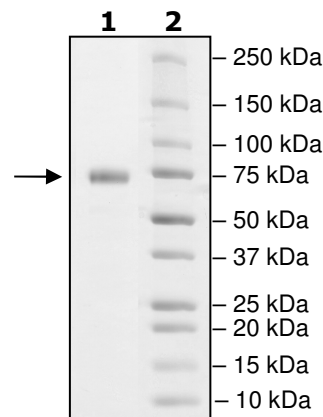
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
2 µg PD-L1(D26A)

**Lane 2:**  
Protein Marker

**MW:** 53.8kDa+glycans

**Purity:** ≥90%



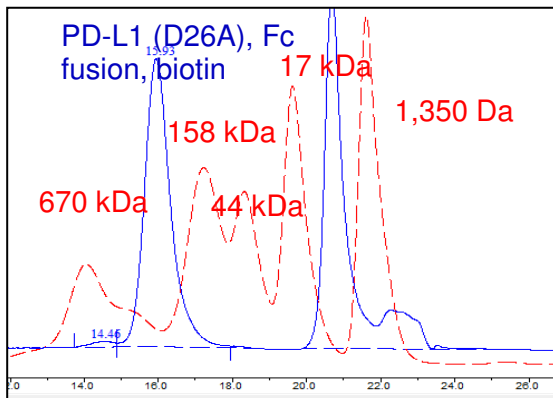
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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](mailto:info@bpsbioscience.com)

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

**Gel Filtration Curve**



< 10%  
 Aggregates

**Sequence:**

FTVTVPK<sup>A</sup>LYVVEYGSNMTIECKFPVEKQLDLAALIVQWEMEDKNI IQFVHGEE<sup>L</sup>DKVQHSSYRQ  
 RARLLKDQLSLGNAALQITDVKLQDAGVYRCMISYGGADYKRITVKVNAPYNKINQRILVVD<sup>P</sup>VT  
 SEHELTCQAEGYPKAEVIWTSSDHQVLSGKTTTTNSKREEKLFNVTSTLRINTTTNEIFYCTFRR  
 LDPEENHTAELVIPELPLAHPPNERT<sup>D</sup><sup>I</sup><sup>E</sup><sup>G</sup><sup>R</sup><sup>M</sup><sup>D</sup><sup>P</sup><sup>K</sup><sup>S</sup><sup>C</sup><sup>D</sup><sup>K</sup><sup>T</sup><sup>H</sup><sup>T</sup><sup>C</sup><sup>P</sup><sup>P</sup><sup>C</sup><sup>P</sup><sup>A</sup><sup>P</sup><sup>E</sup><sup>L</sup><sup>L</sup><sup>G</sup><sup>G</sup><sup>P</sup><sup>S</sup><sup>V</sup><sup>F</sup><sup>L</sup><sup>F</sup><sup>P</sup><sup>P</sup><sup>K</sup><sup>P</sup><sup>K</sup>  
<sup>D</sup><sup>T</sup><sup>L</sup><sup>M</sup><sup>I</sup><sup>S</sup><sup>R</sup><sup>T</sup><sup>P</sup><sup>E</sup><sup>V</sup><sup>T</sup><sup>C</sup><sup>V</sup><sup>V</sup><sup>D</sup><sup>V</sup><sup>S</sup><sup>H</sup><sup>E</sup><sup>D</sup><sup>P</sup><sup>E</sup><sup>V</sup><sup>K</sup><sup>F</sup><sup>N</sup><sup>W</sup><sup>Y</sup><sup>V</sup><sup>D</sup><sup>G</sup><sup>V</sup><sup>E</sup><sup>V</sup><sup>H</sup><sup>N</sup><sup>A</sup><sup>K</sup><sup>T</sup><sup>K</sup><sup>P</sup><sup>R</sup><sup>E</sup><sup>E</sup><sup>Q</sup><sup>N</sup><sup>S</sup><sup>T</sup><sup>Y</sup><sup>R</sup><sup>V</sup><sup>V</sup><sup>S</sup><sup>V</sup><sup>L</sup><sup>T</sup><sup>V</sup><sup>L</sup><sup>H</sup><sup>Q</sup><sup>D</sup><sup>W</sup>  
 LNGKEYKCKVSNKALPAIEKTI SKAKGQPREPQVYTLPPSREEMTKNQVSLTCLVKGFYPSDIA  
 VEWESNGQPENNYKTPPVLDSDGSFFLYSKLTVDKSRWQQGNVFC<sup>S</sup><sup>S</sup><sup>V</sup><sup>M</sup><sup>H</sup><sup>E</sup><sup>A</sup><sup>L</sup><sup>H</sup><sup>N</sup><sup>H</sup><sup>Y</sup><sup>T</sup><sup>Q</sup><sup>K</sup><sup>S</sup><sup>L</sup><sup>S</sup><sup>L</sup>  
 SPGKGGGLNDIFE<sup>A</sup><sup>Q</sup><sup>K</sup><sup>I</sup><sup>E</sup><sup>W</sup><sup>H</sup><sup>E</sup>

Black: PD-L1 (a.a. 19-239)  
 Green: Linker  
 Red: Fc of human IgG1 (a.a.100-330)  
 Blue: Avi-tag™  
 Purple: D26A 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)