

## Data Sheet

### PD-L1 (Monkey), Fc fusion (Human IgG1), Biotin-labeled

Macaca fascicularis, Recombinant, Fc fusion protein, C-terminal Avitag™

**Catalog #:** 71154

**Lot #:** 150529

**Conc.:** 1.37 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; **192(7)**:1027-34.
2. Dong, H., *et al.*, *Nat. Med.* 1999; **5(12)**:1365-9.

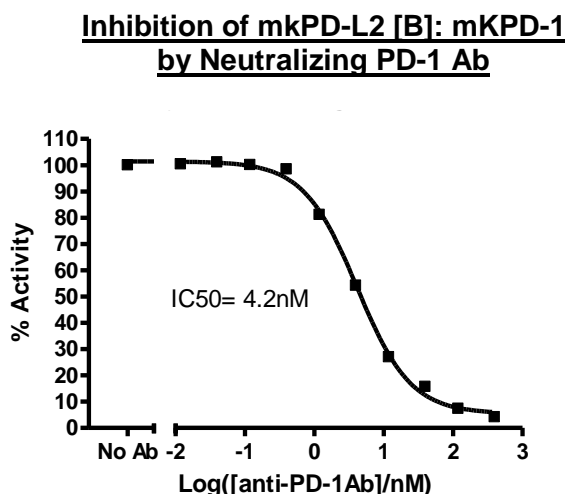
**Description:**

Monkey 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. XP\_005581836.1, a.a. 25-167, fused at the C-terminus to the Fc portion of human IgG1 with an Avitag™. This expressed in a HEK293 cell expression system and is enzymatically biotin-labeled using Avitag™ technology. MW = 54 kDa. Note: this protein runs at an apparent higher molecular weight due to glycosylation.

**Application:**

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

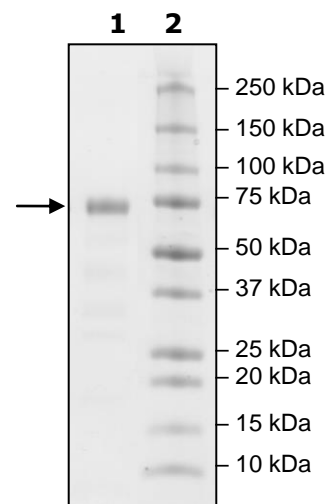
### Quality Assurance



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

**Lane 1:** 2µgPD-L1  
**Lane 2:** Protein Marker

**MW:** 54 kDa + glycans  
**Purity:** ≥80%



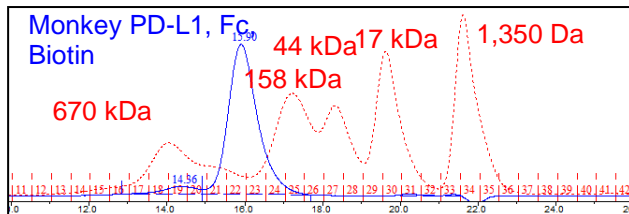
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** info@bpsbioscience.com

### Gel Filtration Curve



<10 % Aggregation

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)