

## Data Sheet

### **PD-1 (I126A), Fc fusion, Biotin-labeled**

Human, recombinant, Fc fusion protein  
with C-terminal AviTag™

**Catalog #:** 71238

**Lot #:** 150827

**Conc.:** 1.16 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 (PD-1)-Fc fusion protein, also known as CD279, PDCD1, and SLEB2, GenBank Accession No. NM\_005018, amino acids 25-167 with an I126A mutation, fused at the C-terminus with the Fc portion of human IgG1, and C-terminal AviTag™, expressed in a HEK293 cell expression system and enzymatically biotin-labeled using AviTag™ technology. MW = 44.5 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

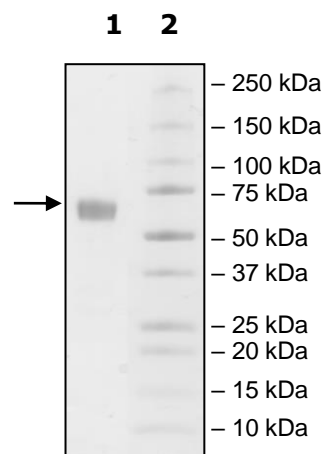
**Application:**

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:**  
3 µg PD-1(I126A)  
**Lane 2:**  
Protein Marker  
**MW:** 44.5 kDa+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 **1.858.202.1401** Fax **1.858.481.8694**

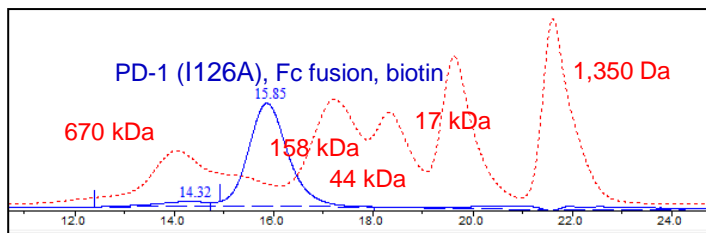
Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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



6405 Mira Mesa Blvd Ste 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

### Gel Filtration Curve



<10% Aggregates

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**  
Or you can Email us at: **support@bpsbioscience.com**  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)