

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_014143, 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.

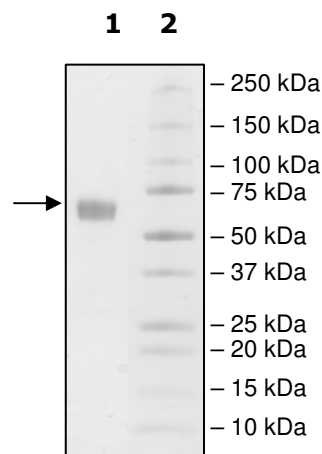
Application:

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 **1.858.829.3082** Fax **1.858.481.8694**

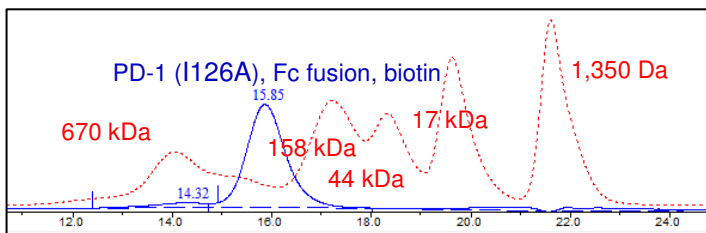
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@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 **1.858.829.3082** 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)