

Data Sheet

PD-1 (P130A), Fc fusion, Biotin-labeled

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

Catalog #: 71246

Lot #: 160303-G **Conc.:** 1.61 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, a.a. 25-167 with an P130A 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

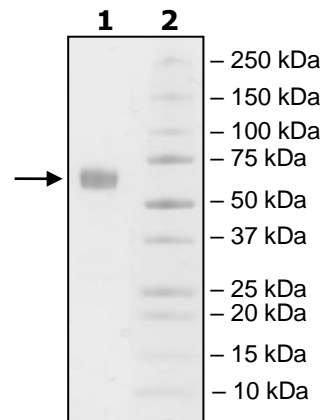
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 µg PD-1 (P130A)

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**

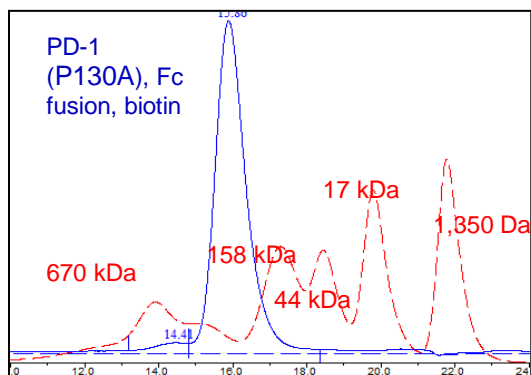
Or you can Email us at: support@bpsbioscience.com

Please visit our website at: 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