

Data Sheet

Mouse PD-1, Fc fusion, Biotin-labeled

Mouse, recombinant, Fc fusion protein,
 C-terminal biotin tag

Catalog #: 71118

Lot #: 150724-G2 **Conc.:** 0.39 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 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 $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

1. Agata, Y., *et al.*, *Internat. Immunol.* 1996; **8(5)**: 765-772.
2. Freeman, G.J. *et al.*, *J. Exp. Med.* 2000; **192(7)**: 1027-1034.

Description:

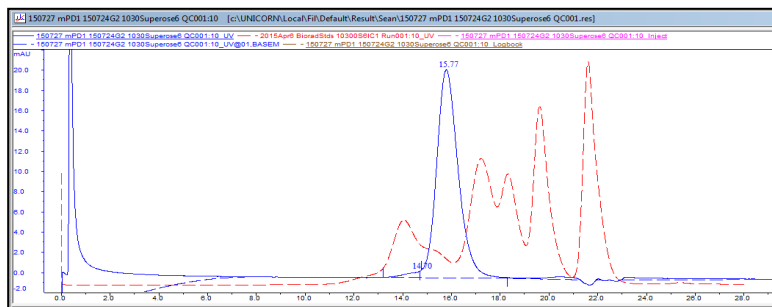
Mouse secreted Programmed Cell Death 1 (PD-1)-Fc fusion protein, also known as CD279, PDCD1, and SLEB2, GenBank Accession No. NM_008798, a.a. 25-167, fused at the C-terminus to the Fc portion of mouse IgG2a and with a C-terminal Avi-tag, expressed in a HEK293 cell expression system. This protein was biotinylated using AviTag™ technology. MW = 45 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

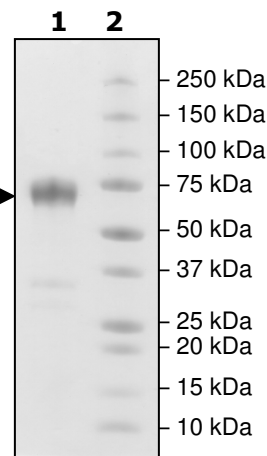
Gel Filtration Trace



4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg PD-1
Lane 2:
Protein Marker

MW: 45 kDa + glycans
Purity: $\geq 90\%$



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