

Data Sheet

Mouse PD-L1, Fc fusion, Biotin-labeled

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

Catalog #: 71119

Lot #: 150205-D

Conc.: 2.47 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
 $\geq 100 \mu\text{g/ml}$ 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:

Mouse 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. NM_021893,
a.a. 18-238, 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
= 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

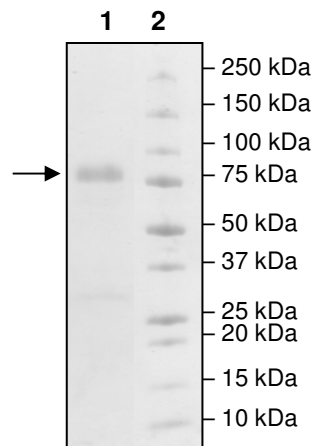
4-20% SDS-PAGE
Coomassie staining

Lane 1:
4 μg PD-L1

Lane 2:
Protein Marker

MW: 53.8 kDa

Purity: $\geq 80\%$



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**

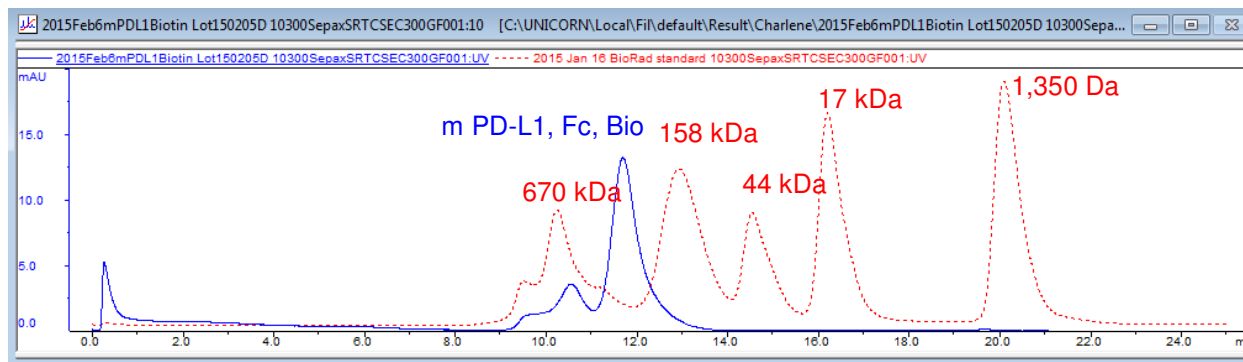
Or you can Email us at: info@bpsbioscience.com

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



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