

Data Sheet

Mouse PD-L1 (CD274), Fc fusion

Mouse, recombinant, C-terminal Fc fusion

Catalog #: 71117

Lot #: 150211

Conc.: 1.0 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. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

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. 19-238, fused at the C-terminus to the Fc portion of mouse IgG2a, expressed in a HEK293 cell expression system. MW = 51.8 kDa.

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:

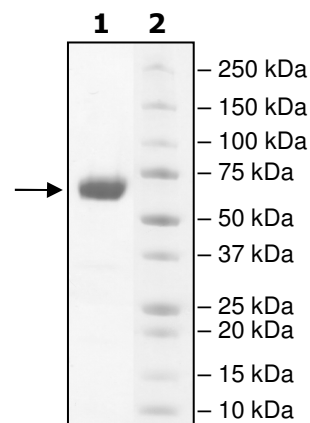
3 μg PD-L1-Fc

Lane 2:

Protein Marker

MW: 51.8 kDa

Purity: $\geq 90\%$



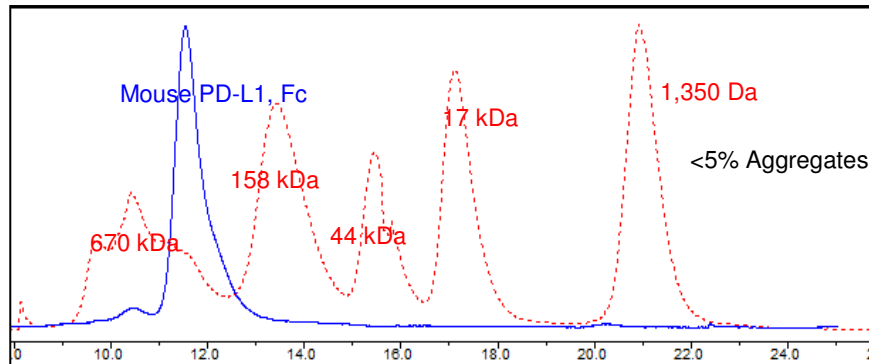
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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

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