

Data Sheet

PD-L1 (Monkey), Fc fusion (Human IgG1)

Macaca fascicularis, Recombinant, Fc fusion protein

Catalog #: 71153

Lot #: 150421

Conc.: 1.63 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:

Monkey 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. XP_005581836.1, a.a. 25-167, fused at the C-terminus to the Fc portion of human IgG1 expressed in a HEK293 cell expression system. MW = 52 kDa. Note: this protein runs at an apparent higher molecular weight due to glycosylation.

Application:

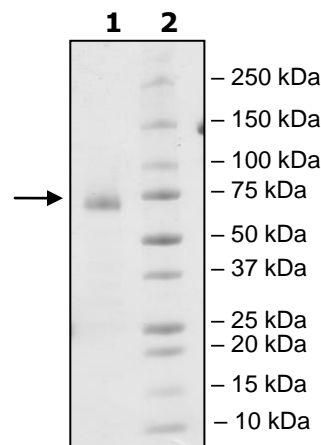
Useful for the study of protein-protein interaction and *in vitro* and *in vivo* protein function studies.

Quality Assurance

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

Lane 1:
3 μg PD-L1-Fc
Lane 2:
Protein Marker

MW: 52 kDa + glycans
Purity: $\geq 90\%$



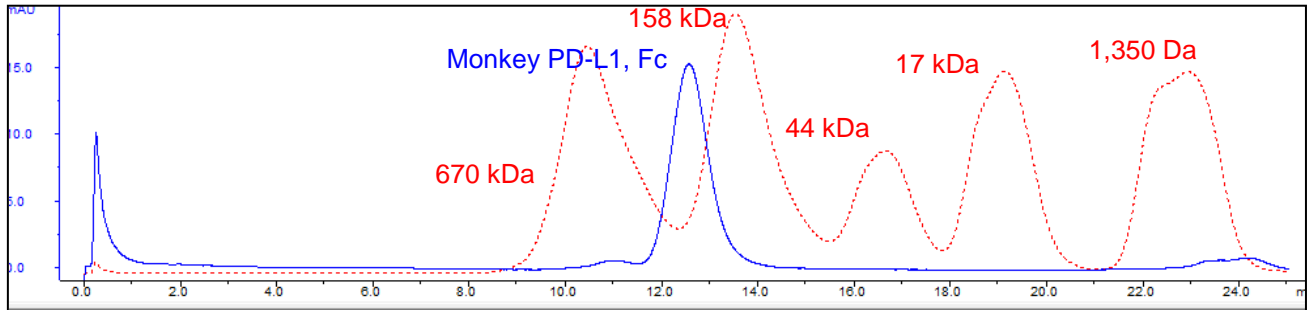
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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



6042 Cornerstone Court W, Ste B
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