

Data Sheet

PD-L1(CD274), FLAG-Tag

Human, Recombinant, C-terminal FLAG-tag

Catalog #: 71183

Lot #: 180607 **Conc.:** 0.86 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:

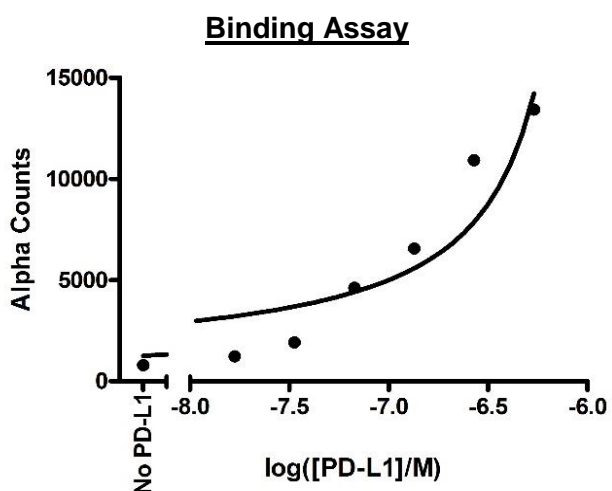
- Freeman, G.J., *et al. J. Exp. Med.* 2000; **192(7)**: 1027-1034
- Dong, H., *et al., Nat. Med.* 1999; **5(12)**: 1365-1369.

Description: Human Programmed Cell Death 1 Ligand 1 (PD-L1), also known as CD274 and B7 homolog 1 (B7-H1), PDCD1 ligand 1, programmed death ligand 1, PDCD1L1, GenBank Accession No. NM_014143, a.a. 19-239, with a C-terminal FLAG-tag, expressed in a HEK293 cell expression system. MW = 26 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Assay Conditions: 10 nM PD-1, Biotin-labeled (BPS Cat. #71109) was mixed with varying amounts of PD-L1, FLAG-tag (BPS Cat. #71183) in PD-1 Assay Buffer for 1 hr. at RT. FLAG-acceptor beads were added and incubated for 30 min at RT followed by strep-conjugated donor beads for 30 min at RT. Alpha counts were read.

Application: Useful for studying protein binding and for screening small molecules and antibodies.

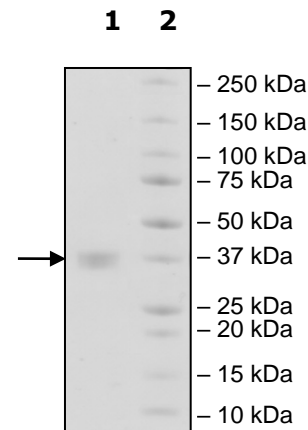
Quality Assurance



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

Lane 1:
4 μg PD-L1
Lane 2:
Protein Marker

MW: 26 kDa+glycans
Purity: $\geq 90\%$



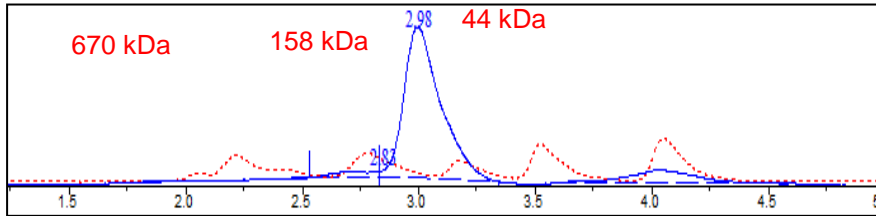
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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



<10% Aggregation

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