

Data Sheet

PD-L1 (K124A), Fc fusion, Avi-Tag, HiP™

Human, Recombinant, Fc fusion protein with C-terminal Avi-tag™

Catalog #: 71993

Lot #: 160927

Conc.: 1.02 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 ≥ 100 µg/ml in PBS + glycerol and stored at -80°C.

References:

1. Freeman, G.J., *et al.*, *J Exp Med.* 2000 Oct 2;**192(7)**:1027-34.
2. Dong, H., *et al.*, *Nat Med.* 1999 Dec;**5(12)**:1365-9.

Description: Human secreted Programmed Cell Death 1 Ligand 1 (PD-L1)-Fc fusion protein, also known as CD274 and B7 homolog 1 (B7-H1) with a K124A mutation, GenBank Accession No. NM_014143, amino acids 19-239, fused at the C-terminus with the Fc portion of human IgG1, with C-terminal Avi-tag™ expressed in a HEK293 cell expression system. MW = 54 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Application:

Useful for the study of protein-protein interaction, epitope mapping, and *in vitro* and *in vivo* protein function.

Quality Assurance

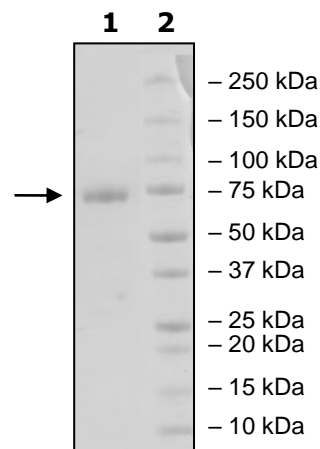
**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 3µg PD-L1(K124a)

Lane 2:
 Protein Marker

MW: 54k Da+glycans

Purity: ≥90%



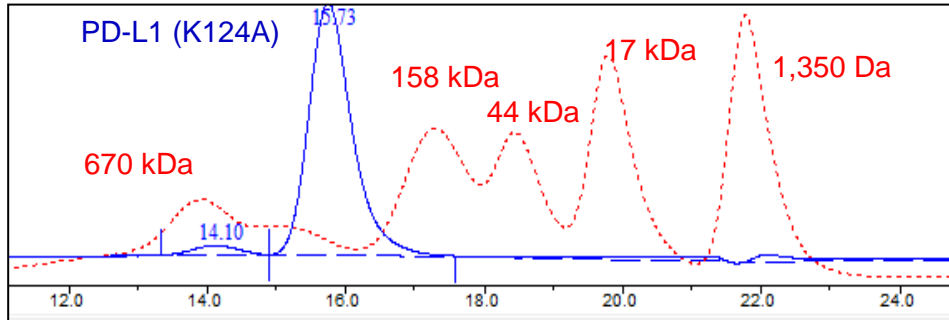
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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