

## Data Sheet

### PD-L1 (Y123A), Fc fusion, Avi-Tag

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

**Catalog #:** 71992

**Lot #:** 160926

**Conc.:** 1.24 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 Y123A 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.

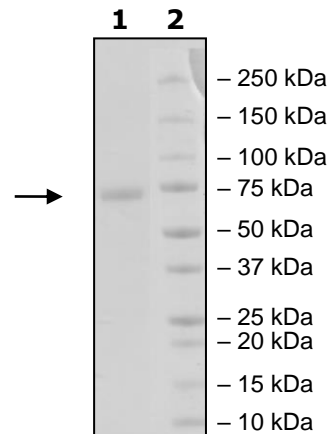
**Application:**

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(Y123A)  
**Lane 2:**  
Protein Marker  
**MW:** 54 kDa + 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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)