

## Data Sheet

### PD-1, Avi-His-Tag, Biotin-Labeled (Mouse)

Mouse, Recombinant, C-terminal Avi-His-Tag,  
Biotin-Labeled

**Catalog #:** 79189

**Lot #:** 170921-2      **Conc.:** 1.62 mg/ml

**Formulated in:** 8mM phosphate, pH7.4,  
110mM NaCl, 2.2mM KCl, and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid  
freeze/thaw cycles. Protein may be diluted to  
 $\geq 100 \mu\text{g/ml}$  in PBS + glycerol and stored at  $-80^{\circ}\text{C}$ .

**References:**

1. Agata, Y., *et al.*, *Internat. Immunol.* 1996;  
**8(5):** 765-772.
2. Freeman, G.J. *et al.*, *J. Exp. Med.* 2000;  
**192(7):** 1027-1034.

**Description:**

Mouse secreted Programmed Cell Death 1 (PD-1)-Fc fusion protein, also known as CD279, PDCD1, and SLEB2, GenBank Accession No. NM\_008798, a.a. 25-167, with a C-terminal Avi-His-tag, expressed in a HEK293 cell expression system. This protein is enzymatically biotinylated using Avi-tag™ technology. MW = 19 kDa. This protein runs at an apparent higher molecular weight due to glycosylation.

---

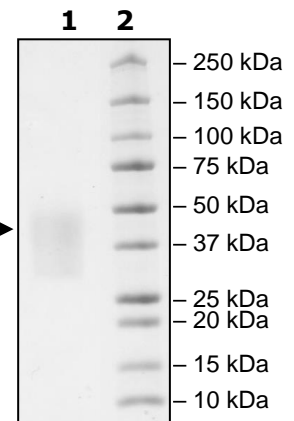
### **Quality Assurance**

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

**Lane 1:**  
2 $\mu\text{g}$  PD-1

**Lane 2:**  
Protein Marker →

**MW:** 19 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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)