

## Data Sheet

### PD-1(PCD1)-FLAG-Avi-His-tag

Human, recombinant, C-terminal FLAG-Avi-His-tag

**Catalog #:** 71198

**Lot #:** 150702-G      **Conc.:** 0.89 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  $-80^{\circ}\text{C}$ . 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:**

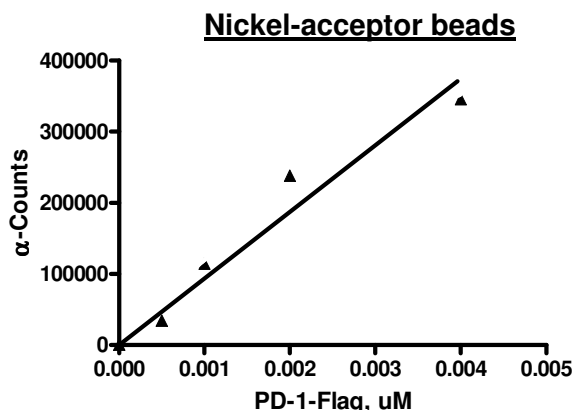
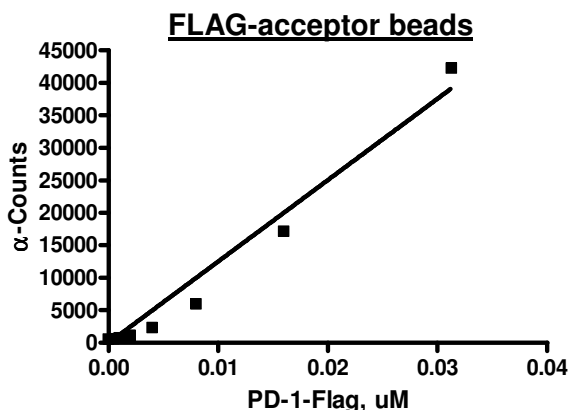
1. Shinohara, T., *et al.*, *Genomics*. 1994 Oct;**23(3)**:704-6.
2. James, E.S., *et al.*, *Genes Immun*. 2005 Aug;**6(5)**:430-7.

**Description:** Human secreted programmed cell death 1 (PD-1), also known as PDCD1, SLEB2, CD279 and HPD-L, GenBank Accession No. NM\_005018, a.a. 25-167, with C-terminal FLAG-Avi-His-tags, MW=20 kDa, expressed in a HEK293 cell expression system. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

**Assay Conditions:** 10 nM PD-L1, Biotin-labeled (BPS Cat. #71105) was mixed with varying amounts of PD-1, FLAG-tag (BPS Cat. #71198) in PD-1 Assay Buffer for 1 hr at RT. FLAG-acceptor beads (left) or Nickel-acceptor beads (right) were added and incubated for 30 min at RT followed by streptavidin-conjugated donor beads for 30 min at RT. Alpha counts were read.

**Application:** Useful for studying protein binding and screening small molecular inhibitors.

## Quality Assurance



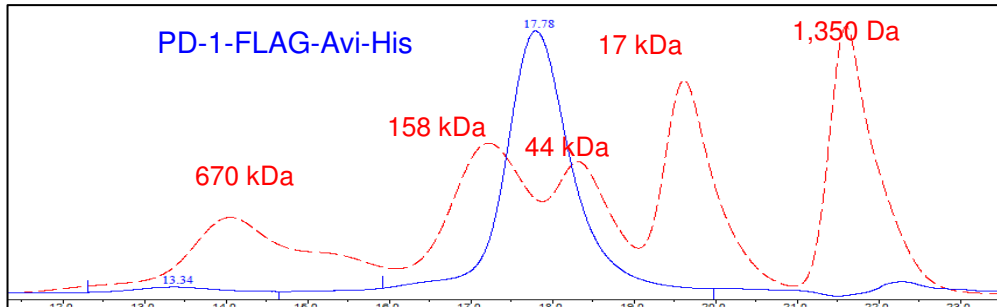
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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)

**Gel Filtration Curve**



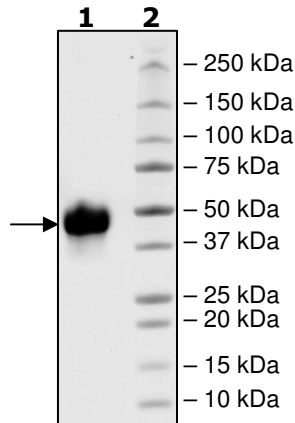
<5%  
 Aggregation

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

**Lane 1:**  
 3 µg PD-1

**Lane 2:**  
 Protein Marker

**MW:** 20 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)