

## Data Sheet

### PD-1, Fc Fusion (Woodchuck)

Woodchuck, Recombinant, C-terminal  
Fc-Fusion Protein

**Catalog #:** 79314

**Lot #:** 171010      **Conc.:** 1.01 mg/ml

**Formulated in:** 8 mM Phosphate, pH 7.4,  
400 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:

Woodchuck secreted programmed cell death 1 (PD-1) also known as, PDCD1, SLEB2, CD279 and HPD-L, GenBank Accession No. HQ403652.1, a.a 25-172, fused at the C-terminus to the Fc portion of Human IgG1, expressed in a HEK293 cell expression system. MW = 43 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

### Quality Assurance

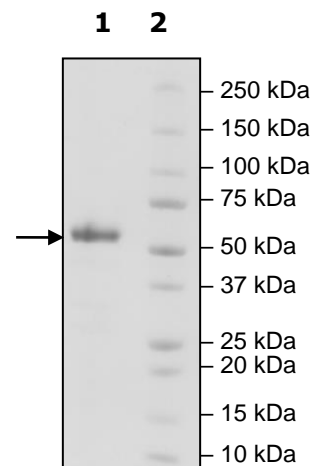
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4 $\mu\text{g}$  PD-1(Woodchuck)

**Lane 2:**  
Protein Marker

**MW:**  
43kDa + 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)