

## Data Sheet

### CTLA4 (Woodchuck), Fc-Fusion (Human), Avi-tag

Woodchuck, Recombinant, N-terminal  
 Fc-Avi-Tag

**Catalog #:** 100037

**Lot #:** 180404      **Conc.:** 0.74 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. Dilek, N., *et al.*, *PLOS ONE*. 2013 Dec. 23: DOI: 10.1317.
2. Leach, D.R., *et al.*, *Science*. 1996; **271(5256)**:1734-1736.

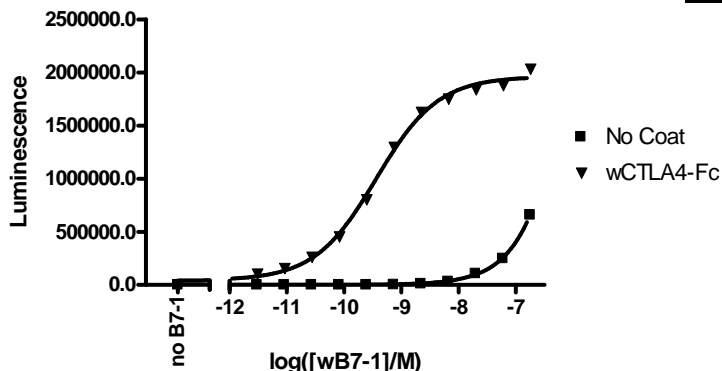
**Description:** Woodchuck secreted CTLA4, also known as Cytotoxic T-lymphocyte-associated protein 4 and CD152, with N-terminal Avi-Tag™ fused to the Fc region of Human IgG1, a.a. 36-162 expressed in a HEK293 cell expression system. MW = 42 kDa (monomer). This protein runs at a higher MW by SDS-PAGE due to glycosylation.

**Assay Conditions:** Reaction done similar to human CTLA4:B7-1[biotin] assay kit (BPS Catalog # 72009) protocol. Buffer is PBS with 0.1% BSA. Blocking buffer is Superblock + 0.05% Tween-20. The coat protein, woodchuck CTLA4 (wCTLA4), was added ON at 4 deg (various concentrations as indicated). Binding reaction initiated with addition of woodchuck B7-1-biotin (wB7-1-biotin) for 2h at RT. Binding was detected using Strep-HRP generated luminescence.

**Applications:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance

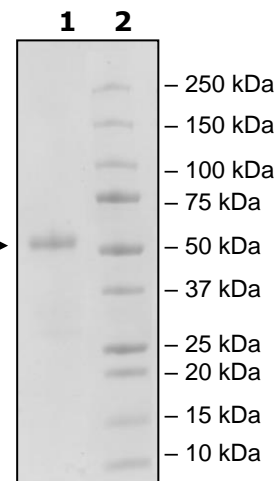
**Woodchuck CTLA4: B7-1 Interaction**



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

**Lane 1:**  
 4μg CTLA4  
 (Woodchuck)  
**Lane 2:**  
 Protein Marker

**MW:**  
 42 kDa + glycans  
**Purity:**  $\geq 90\%$



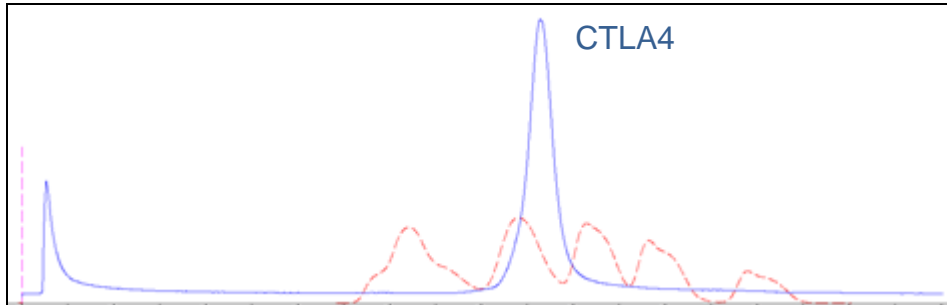
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto: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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)