

Data Sheet

B7-1 (Woodchuck), Fc-fusion (Human), Avi-Tag, Biotin-labeled

Woodchuck, Recombinant, Fc fusion protein,
 C-terminal Avi-Tag, Biotin labeled

Catalog #: 100040

Lot #: 180405 **Conc.:** 0.86 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 $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .*

References:

1. Byrum, J.N., Van Komen, J.S., Rodgers, W., *J. Immunol.* 2013 Sep 15; **191(6)**:3073-3081.
2. Legat, A., *et al.*, *Front Immunol.* 2013 Dec 19; **4**:455.

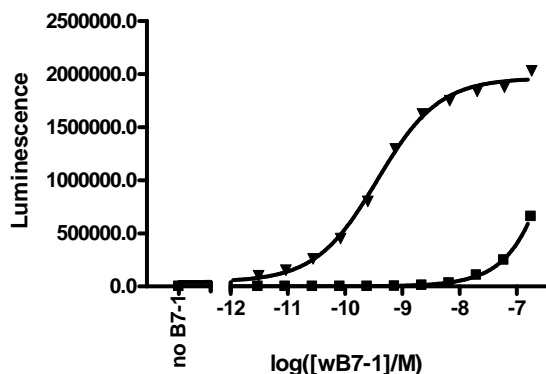
Description: Woodchuck B7-1, also known as CD80, with C-terminal Avi-Tag fused to the Fc-region of Human IgG1. a.a 30-237, expressed in a HEK293 cell expression system and enzymatically biotin-labeled using Avi-tag™ technology. MW = 52 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

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.

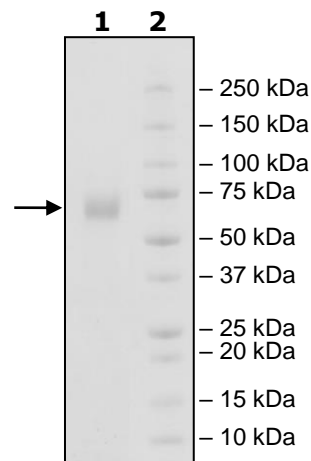
Quality Assurance

Woodchuck CTLA4: B7-1 Interaction



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

Lane 1:
 B7-1 (Woodchuck)
Lane 2:
 Protein Marker
MW:
 52kDa + 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
 Please visit our website at: www.bpsbioscience.com