

## Data Sheet

### **B7-1(CD80) (Mouse), Fc fusion, Avi-Tag, Biotin-labeled**

Mouse, Recombinant,  
 C-terminal Fc-Avi-tag, Biotin labeled

**Catalog #:** 100478

**Lot #:** 190815

**Conc.:** 1.01 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.*

**References:**

1. Pentcheva-Hoang, T., *et al. Immunity.*  
 2004 Sep; **21(3)**: 401-413.
2. Hirano, F., *et al., Cancer Research.*  
 2005 Feb; **65(3)**: 1089-1096.

**Description:** Mouse secreted B7-1, also known as CD80, GenBank Accession No. NM\_009855, a.a 38-246, fused at the C-terminus of the Fc portion of human IgG1, with C-terminal Avi-tag, expressed in a HEK293 cell expression system and enzymatically biotinylated using Avi-tag™ technology. MW=52 kDa. Biotinylation is confirmed to be  $\geq 90\%$ .

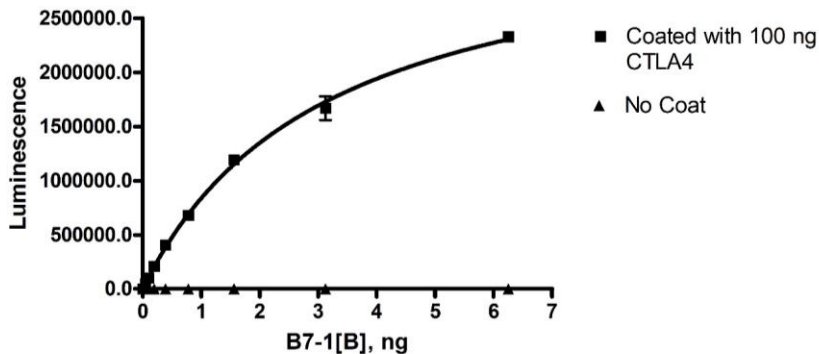
**Assay Conditions:** Reaction done similar to CTLA4:B7-1[biotin] assay kit (Catalog # 72009) protocol. Buffer is PBS + 0.1% BSA. Blocking buffer is Superblock + 0.05% Tween-20.

Coat protein (mouse CTLA4) was added overnight at 4 deg 100 ng/well. Various amounts of biotinylated ligand (mouse B7-1[B]) (0-50 ng) added for 2h at RT. Strep-HRP added for 1h at RT.

**Applications:** Useful for the study of protein-protein interaction and in vitro and in vivo protein function studies.

### Quality Assurance

**Mouse CTLA4: B7-1 [Biotin] Interaction**

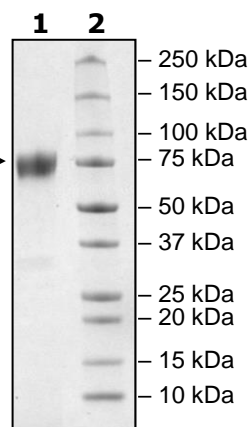


**4-20% SDS-PAGE  
 Coomassie staining**

**Lane 1:** → 4µg B7-1  
**Lane 2:** Protein Marker

**MW:** 52  
 kDa+glycans

**Purity:**  $\geq 90\%$



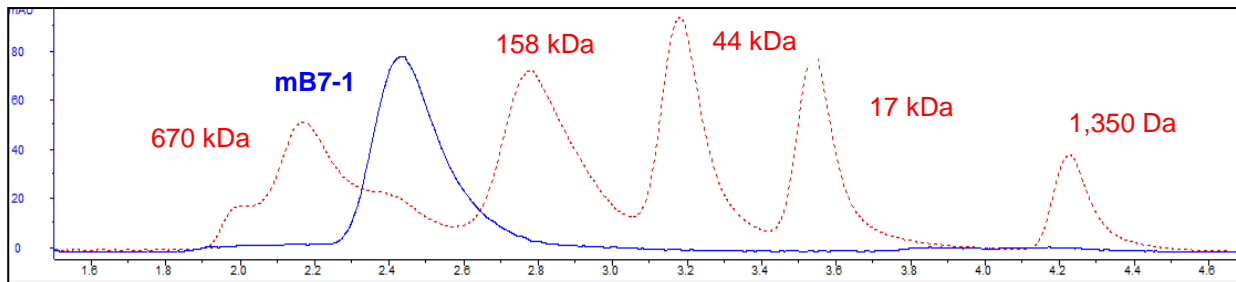
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** 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:** 1.858.202.1401  
**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 **1.858.202.1401** 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)