

## Data Sheet

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

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

**Catalog #:** 71114

**Lot #:** 160602

**Conc.:** 0.69 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:

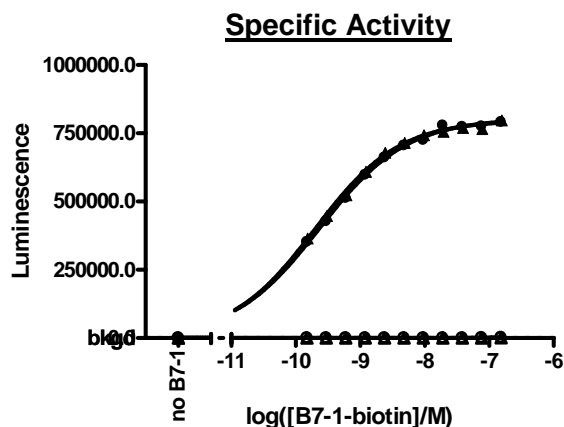
1. Byrum, J.N., Van Komen, J.S.,  
Rodgers, W., *J. Immunol.* 2013 Sep  
15;191(6):3073-3081.
2. Legat, A., et al., *Front Immunol.* 2013  
Dec 19;4:455.

**Description:** Human secreted B7-1-Fc  
fusion protein, also known as CD80, with C-  
terminal Avi-Tag. GenBank Accession No.  
NM\_005191, a.a 35-242 expressed in a  
HEK293 cell expression system and  
enzymatically biotin-labeled using Avi-tag™  
technology. MW = 53 kDa (monomer). This  
protein runs at a higher M.W. by SDS-  
PAGE due to glycosylation.

**Assay Description:** Reaction done  
according 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 was added  
ON at 4 deg 100 ng/well. Various  
concentrations of biotinylated ligand (0-  
8ng/ul, 0.5 ng/ul is the recommended kit  
condition) added for 2h at RT. Strep-HRP  
added for 1h at RT.

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

## Quality Assurance



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

**Lane 1:**

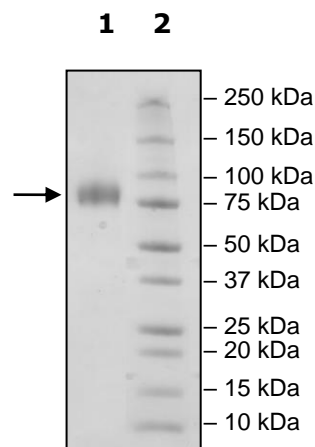
3μg B7-1

**Lane 2:**

Protein Marker

**MW:** 53 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**

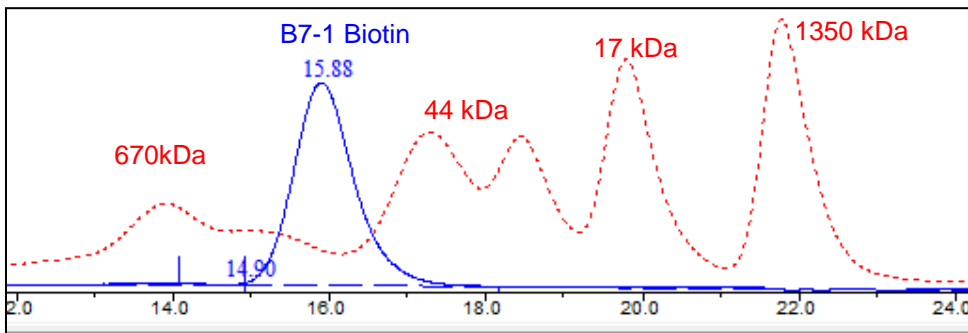
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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)