

## Data Sheet

### **B7-2 (CD86), Fc fusion, Biotin labeled**

Human, recombinant, C-terminal Fc fusion  
Protein, C-terminal Avi-tag™

**Catalog #:** 71159

**Lot #:** 150417

**Conc.:** 0.38 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  
≥ 100 µg/ml in PBS + glycerol and stored at  
-80°C.*

**References:**

1. Ertelt, J.M., *et al.* *J. Leukocyte Biol.* 2013; **94**(2): 367-376.
2. Kuchroo, V.K., *et al.*, *Cell.* 1995; **80**(5):707-718.

**Description:**

Human secreted B7-2, Fc fusion protein, also known as T-lymphocyte activation antigen CD86, B7.2, FUN-1, B70, BU63, and CD86. GenBank Accession No. NM\_006889, a.a. 20-239 expressed in a HEK293 cell expression system. This protein is fused at the C-terminus to the Fc region of human IgG1 and is enzymatically biotinylated using Avi-tag™ technology. MW = 53.8 kDa (monomer). This protein runs at a higher MW by SDS-PAGE due to glycosylation.

**Application:**

Useful for studying protein binding and screening small molecules.

## **Quality Assurance**

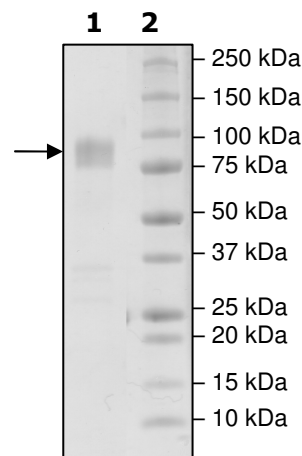
**4-20% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
4 µg B7-2-biotin

**Lane 2:**  
Protein Marker

**MW:** 53.8 kDa+glycans

**Purity:** ≥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**

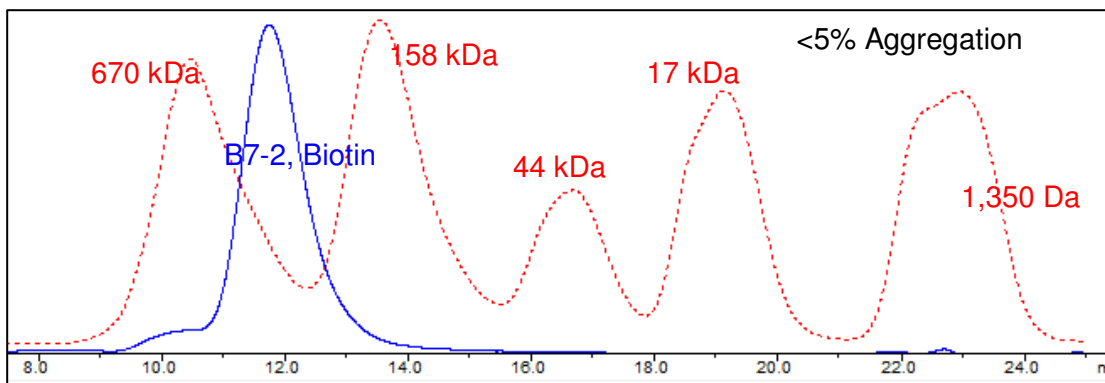
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E  
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



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)