

## Data Sheet

### **CD83, Fc-fusion (IgG1), Avi-Tag, Biotin labeled**

Human, recombinant, C-terminal Fc-Avi-Tag,  
Biotin labeled

**Catalog #:** 100348

**Lot #:** 190411

**Conc.:** 1.82 mg/ml

**Formulated In:** 8 mM Phosphate, pH 7.4,  
110 mM NaCl, 2.2 mM KCl, and 20%  
glycerol.

**Stability:** Stable for at least 6 months at  
-80°C. Avoid freeze/thaw cycles.

#### **References:**

1. Zhou, L.J, *et al.*, *The Journal of Immunology*. 1995 Apr; **154(8)**: 3821-3835.
2. Lechmann, M. *et al.*, *Trends in Immunology*. 2002 Jun; **23(6)**: 273-275.

**Description:** Human CD83, also known as Cell Surface Protein HB15, B-Cell Activation Protein, CD83 Antigen, and HCD83, GenBank Accession No. NM\_004233, a.a. 20-144 fused at the C-terminus to the Fc portion of human IgG1 followed by a C-terminal Avi-tag™ and expressed in a HEK293 cell expression system. MW=43 kDa. This protein is enzymatically biotinylated using Avi-Tag technology. Biotinylation is confirmed to be ≥90%.

---

### **Quality Assurance**

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

**Lane 1:**

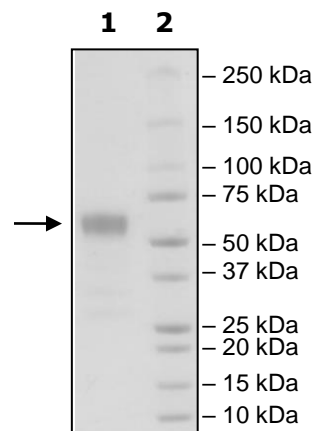
4µg CD83

**Lane 2:**

Protein Marker

**MW:** 43 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 **1.858.202.1401** 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)