

## Data Sheet

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

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

**Catalog #:** 100287

**Lot #:** 190307 **Conc.:** 3.62 mg/ml

**Formulation:** 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. Thompson, J.S., *et al.*, *Science*. 2001 Sep; **293(5537)**: 2108-2111.
2. Castigli, E. *et al.*, *JEM*. 2005 Jan; **201(1)**: 35.

**Description:** Human BAFF-R, also known as TNF Receptor Superfamily Member 13C, Tumor Necrosis Factor Receptor Superfamily Member 13C, BLyS Receptor 3, BAFFR, and B Cell-Activating Factor Receptor, GenBank Accession No. NM\_052945, a.a. 23-71 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=33.5 kDa. This protein is enzymatically biotinylated using Avi-Tag technology.

---

## **Quality Assurance**

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

**Lane 1:**

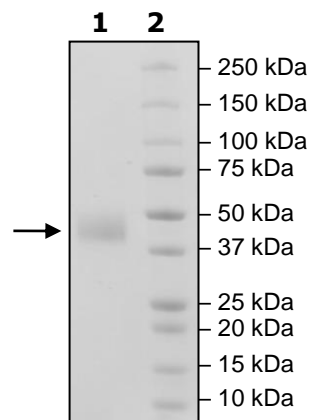
11µg BAFF-R

**Lane 2:**

Protein Marker

**MW:** 33.5 kDa

**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)