

Data Sheet

BAFF-R, Fc-fusion (IgG1), Avi-Tag

Human, recombinant, C-terminal Fc-Avi-Tag
Catalog #: 100286
Lot #: 190220 **Conc.:** 0.99 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.

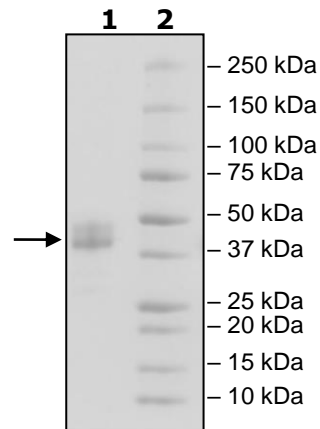
Description: Human BAFF-R, also known as TNF Receptor Superfamily Member 13C, Tumor Necrosis Factor Receptor Superfamily Member 13C, BlyS Receptor 3, BAFFR, and B Cell-Activating Factor Receptor, GenBank Accession No. NM_052945, a.a. 23-71 fused at the C terminus to the Fc portion of human IgG1 followed by a C-terminal Avi-tag™ and expressed in a HEK293 cell expression system. MW=33.5 kDa.

Quality Assurance

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

Lane 1:
4µg 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
Please visit our website at: www.bpsbioscience.com