

## Data Sheet

### **RANK, Fc fusion (IgG1), Biotin Labeled**

Human, Recombinant, C-terminal Avi-tag

**Catalog #:** 70822

**Lot #:** 160811

**Conc.:** 0.71 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. Annis, A., J. *American Chemical Society.*, 2004 May; **126(4)**: 15495-15503.
2. Yan, T., J. *Cellular Biochemistry.* 2001 Aug; **83(2)**:320-325.

**Description:**

Human Receptor Activator of Nuclear Factor  $\kappa$  B (RANK), also known as TRANCE Receptor and TNFRSF11A, (Genbank Accession No. NM\_003839), a.a. 30-212, fused at the C-terminus to the Fc portion of human IgG1, with a C-terminal Avi-tag, expressed in a HEK293 cell expression system. This protein is enzymatically biotinylated using AviTag™ technology. MW = 49 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation

**Application:**

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Assurance

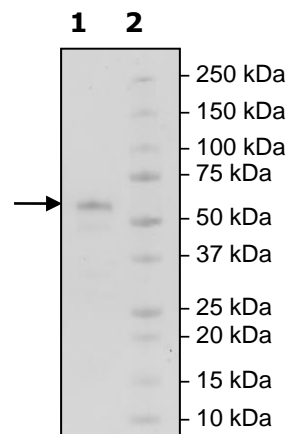
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
2 $\mu\text{g}$  RANK

**Lane 2:**  
Protein Marker

**MW:** 49 kDa + glycans

**Purity:**  $\geq 71\%$



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**

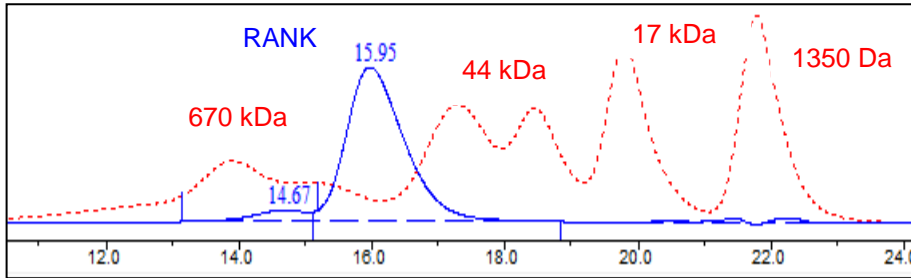
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



6042 Cornerstone Court West, Suite 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)