

Data Sheet

RANK, Fc fusion (IgG1), Avi-tag

Human, Recombinant, C-terminal Avi-tag

Catalog #: 70823

Lot #: 160809

Conc.: 0.8 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 $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°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 κ 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. 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:

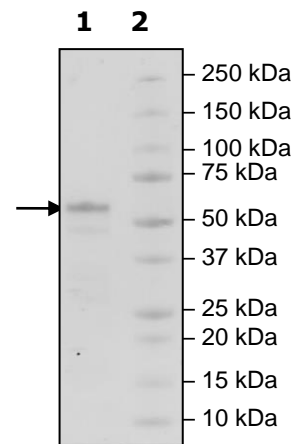
2 μg RANK

Lane 2:

Protein Marker

MW: 49 kDa + glycans

Purity: $\geq 73\%$



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**

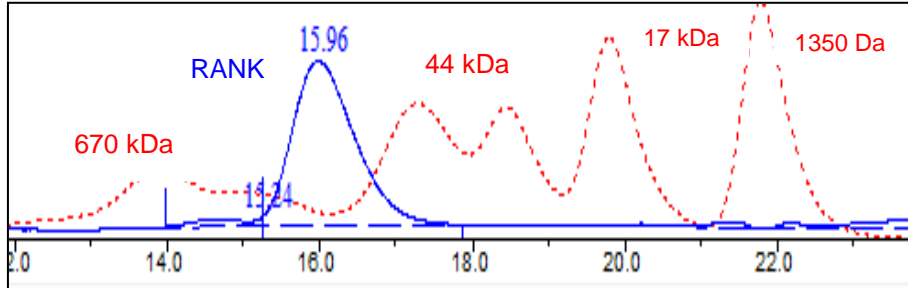
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: 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

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
Please visit our website at: www.bpsbioscience.com