

Data Sheet

Anti-RANKL Neutralizing Antibody

Recombinant, Human, Monoclonal,
Anti-RANKL

Catalog #: 100874

Lot #: 200923 **Conc.:** 3.45 mg/ml

Host Species | Isotype:

Human | IgG1

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

Stability: Stable for at least 12 months at -80°C. Avoid freeze/thaw cycles.

References:

1. Ominsky, M., *et al.*, *Bone*. 2011 Aug; **49(2)**: 162-173.
2. Kostenuik, P. J., *et al.*, *Bone*. 2011 Aug; **49(2)**: 151-161

Description: Human, recombinant monoclonal neutralizing antibody recognizing the RANKL protein, also known as TNFSF11 antibody and CD254 antibody. This antibody recognizes human RANKL. It has not been tested with other species. MW= 50 kDa (Heavy chain); 24 kDa (Light chain).

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: Useful for studying the inhibition of RANKL in cellular assays

Quality Assurance

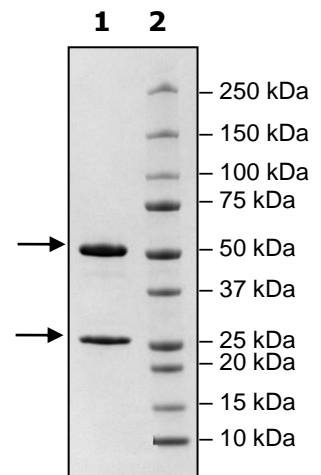
4-20% SDS-PAGE
Coomassie staining

Lane 1:
5µg Anti-RANKL

Lane 2:
Protein Marker

MW:
Heavy chain: 50 kDa
Light chain: 24 kDa

Purity: ≥90%

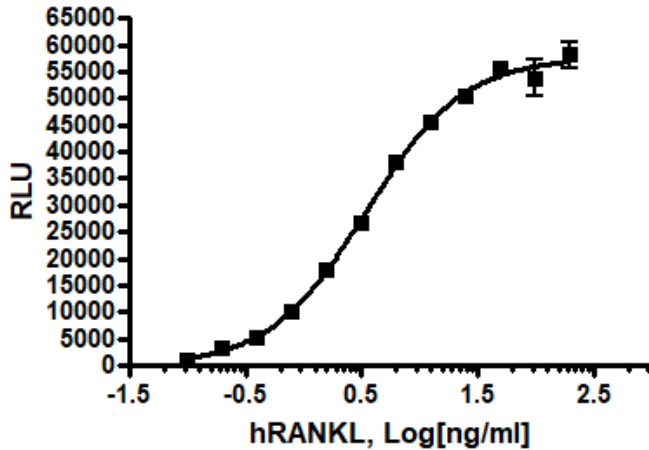
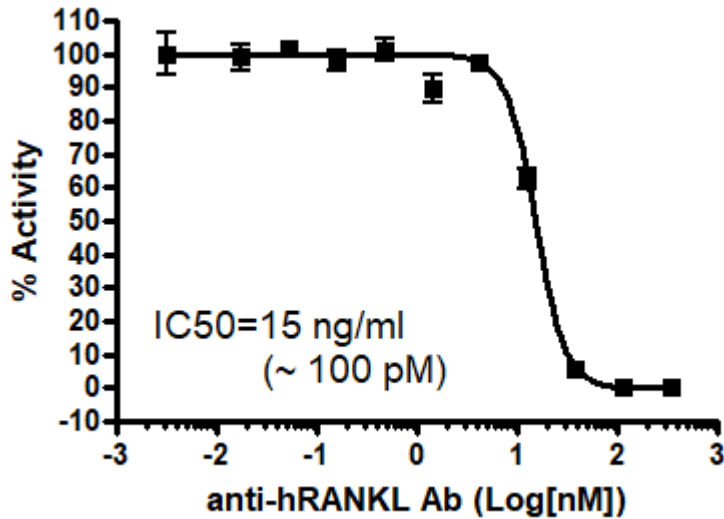


OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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

Figure 1. Human RANKL dose response in NF- κ B reporter (Luc)-HEK293**Figure 2. NFB Inhibition by RANKL**

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