

## Data Sheet

### IL-17RA, Fc fusion, Avi-Tag

Human, Recombinant, C-terminal Fc-Avi-tag

**Catalog #:** 11264

**Lot #:** 161102

**Conc.:** 0.80 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.

**References:**

1. Wright, J.F., *et al.* *J. Immun.* 2008; **181(4)**: 2799-2805.
2. Rickel, E.A., *et al.*, *J. Immun.* 2008; **181(6)**: 4299-4310.

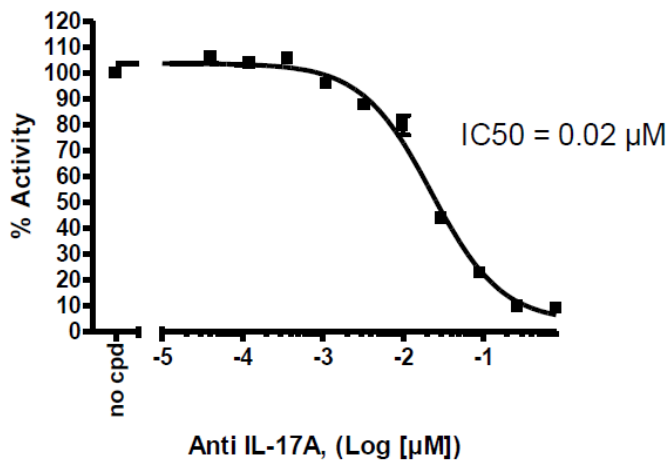
**Description:** Human Interleukin-17RA (IL-17RA), also known as CD217 and CANDF5, GenBank Accession No. NM\_014339, a.a 33-320, fused at the C-terminus of the Fc portion of human IgG1, with C-terminal Avi-tag<sup>TM</sup>, expressed in a HEK293 cell expression system, MW=62 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

**Assay Conditions:** IL-17RA-IL-17A[B] binding activity was measured using the IL-17RA:IL-17A[Biotin] Binding Assay Kit, BPS Bioscience #79891 and Anti-IL-17A Antibody, BPS Bioscience #91015.

**Applications:** Useful for studying protein binding and screening small molecules for drug discovery.

## Quality Assurance

### IL-17RA:IL17[Biotin] Interaction



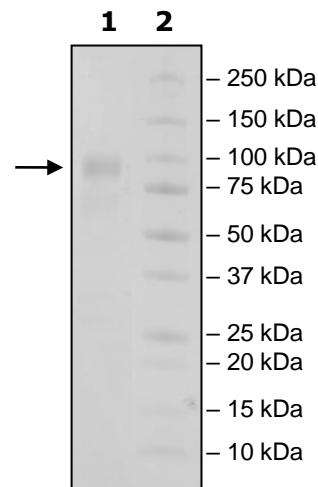
### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
3µg IL17 RA

**Lane 2:**  
Protein Marker

**MW:** 62 kDa

**Purity:** ≥80%



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)