

Data Sheet

IL-17A Neutralizing Antibody

Recombinant, Human Monoclonal
Anti-IL-17A

Catalog #: 91015

Lot #: 151110

Conc.: 1.21 mg/ml

Host Species | Isotype:

Human | IgG1k

Formulated in: 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. Store
diluted antibody with carrier protein (BSA 0.1
– 0.5%) at 4 °C, stable for 2 weeks.*

References:

1. Lubberts, E., *et al.*, *Arthritis Rheum.* 2004; **50(2)**: 650-659.
2. Cheng, X., *et al.*, *Atherosclerosis.* 2011; **215(2)**: 471-474.

Description: Neutralizing recombinant human chimeric antibody recognizing human IL-17A. This antibody has not been tested for cross-reactivity with other species.

IC50: 4.5 nM using the IL17-RA[Biotinylated]:IL-17 Inhibitor Screening Assay Kit (BPS Cat. #72060).

Assay Conditions: Antibody was added at various concentrations using the IL-17RA[Biotin]:IL-17A Inhibitor Screening binding assay (BPS Cat. #72060). Assay was performed according to the recommended protocol.

Purification: Purified using protein A affinity chromatography from HEK293 supernatants.

Application: Useful for binding studies and for the inhibition of IL-17RA:IL-17A interactions.

Image(s):

Inhibition of IL-17RA[B]-IL-17A by Anti-IL-17A

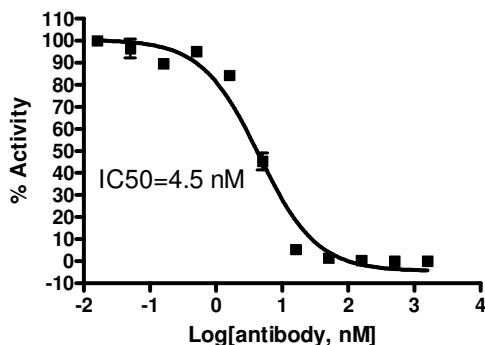


Figure 1. Wells were coated with 100 ng of IL-17A (BPS Cat. #91014). 10 ng of IL-17RA, Biotin-labeled (BPS Cat. #91013) and varying concentrations of IL-17A Neutralizing Antibody (BPS Cat. #91015) were added and plate was incubated at room temperature for 2 hours. Streptavidin-HRP was added, followed by HRP substrate, and chemiluminescence detection.

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