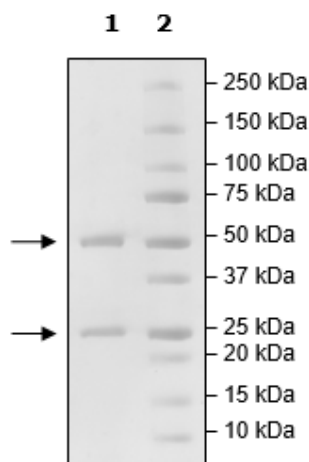


Product Information

Description:	Neutralizing recombinant human chimeric antibody recognizing human IL-17A. This antibody has not been tested for cross reactivity with other species.
Concentration:	1.21 mg/ml
Species:	Human
Isotype:	IgG1κ
Formulated In:	8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	Heavy Chain: 53 kDa; Light Chain: 26 kDa
Purity:	≥90%
Assay Conditions:	Antibody was added at various concentrations using the IL17RA[Biotin]:IL-17A Inhibitor Screening Assay Kit (BPS Bioscience #72060). Assay was performed according to the recommended protocol.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



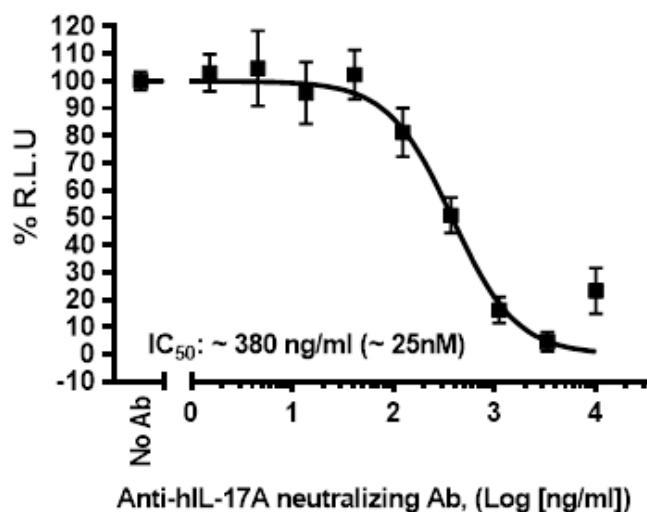
Inhibition of NF- κ B luciferase reporter activity by anti-IL-17 antibody

Figure 1. NF- κ B luciferase reporter HEK293 cells (BPS Bioscience #60650) were plated at a density of 40,000 cells/well in a 96-well plate. The following day, cells were incubated with increasing concentrations of the anti-IL-17A Neutralizing Antibody for 1 hour before stimulation with 100 ng/ml IL-17 for 5-6 hours. Luciferase activity was measured using the ONE-Step™ Luciferase Assay system (BPS Bioscience #60690). Background (cell-free wells) was subtracted from all other values. Results are expressed as percent of positive control (IL-17-stimulated in the absence of antibody).

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

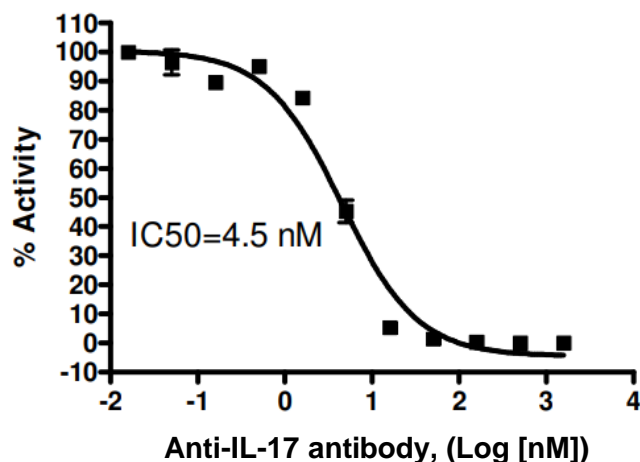


Figure 2. Wells were coated with 100 ng of purified IL-17A (BPS Bioscience #91014). After washing and blocking, 10 ng of Biotin-labeled IL17RA, (BPS Bioscience #91013) and varying concentrations of IL-17A Neutralizing Antibody (BPS Bioscience #91015) were added and the plate was incubated at room temperature for 2 hours. Streptavidin-HRP was added, the plate was washed, followed by addition of the HRP substrate, and chemiluminescence detection.