

Validation data for Anti-mIL-13-mIgG1

<https://www.invivogen.com/anti-mil13-igg>

For research use only

Version 22I27-AK

Anti-mIL-13-mIgG1 is a recombinant monoclonal antibody (mAb) designed to efficiently neutralize mIL-13 *in vitro*. Interleukin 13 (IL-13) is a cytokine that plays an important role in regulating inflammation and immune responses by inhibiting the production of pro-inflammatory cytokines, such as TNF- α . The ability of Anti-mIL-13-mIgG1 to neutralize the biological activity of mIL-13 has been confirmed using InvivoGen's HEK-Blue™ IL-4/IL-13 cells (**Figure 1**).

Dose-dependent neutralization of mIL-13 signaling

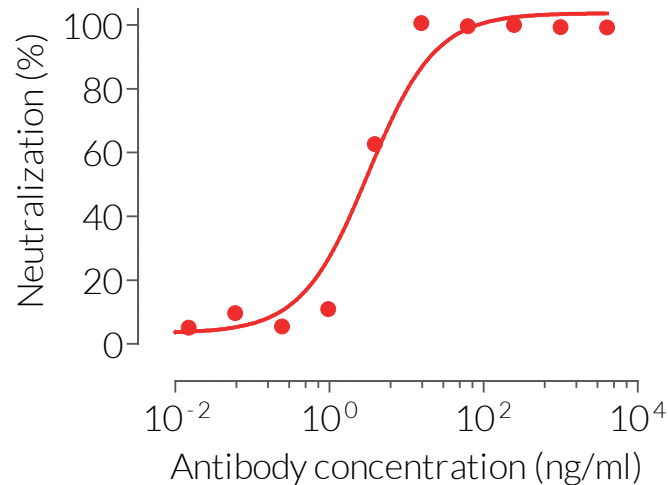


Figure 1. Anti-mIL-13-mIgG1 is a potent neutralization antibody. Increasing concentrations of Anti-mIL-13-mIgG1 were incubated for 30 minutes with the recombinant cytokine mIL-13 (1 ng/ml) prior to the addition of HEK-Blue™ IL-4/IL-13 cells. After overnight incubation, SEAP activity in the cell culture supernatant was assessed using QUANTI-Blue™ Solution. Data are shown in percentage (%) of neutralization.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Asia: +852 3622-3480
E-mail: info@invivogen.com