

# Validation data for Anti-mIL-13-mIgG1 InvivoFit™

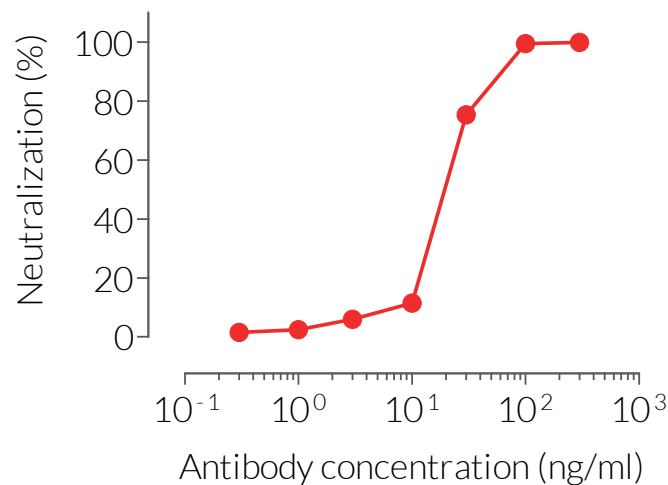
<https://www.invivogen.com/anti-mil13-igg>

For research use only

Version 22J12-AK

Anti-mIL-13-mIgG1 InvivoFit™ is a recombinant monoclonal antibody (mAb) designed to efficiently neutralize mIL-13. 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- $\alpha$ . The ability of Anti-mIL-13-mIgG1 InvivoFit™ 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 InvivoFit™ is a potent neutralization antibody.** Increasing concentrations of Anti-mIL-13-mIgG1 InvivoFit™ were incubated for 30 minutes with the recombinant cytokine mIL-13 (5 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](mailto:info@invivogen.com)