

# Validation data for Anti-hIL-33-IgG

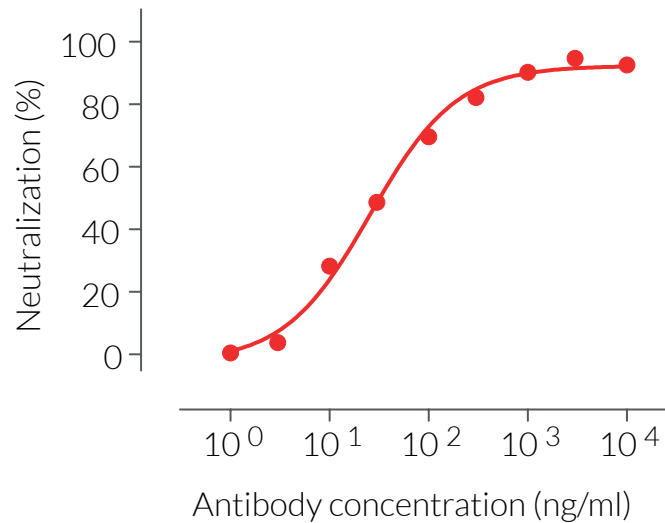
<https://www.invivogen.com/anti-hil33-igg>

For research use only

Version 23B21-MM

Anti-hIL-33-IgG is a monoclonal antibody (mAb) designed to efficiently neutralize human interleukin 33 (hIL-33) *in vitro*. IL-33 is a pro-inflammatory cytokine that shares structural and functional characteristics with the IL-1 cytokine family. The ability of Anti-hIL-33-IgG to neutralize the biological activity of hIL-33 has been confirmed using InvivoGen's HEK-Blue™ IL-33 cells (Figure 1).

## Dose-dependent neutralization of hIL-33 signaling



**Figure 1. Anti-hIL-33-IgG is a potent neutralization antibody.** Increasing concentrations of Anti-hIL-33-IgG were incubated for 30 minutes with the recombinant cytokine hIL-33 (1 ng/ml) prior to the addition of HEK-Blue™ IL-33 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)