

# Validation data for Anti-mIL-1 $\beta$ -mIgG1 InvivoFit™

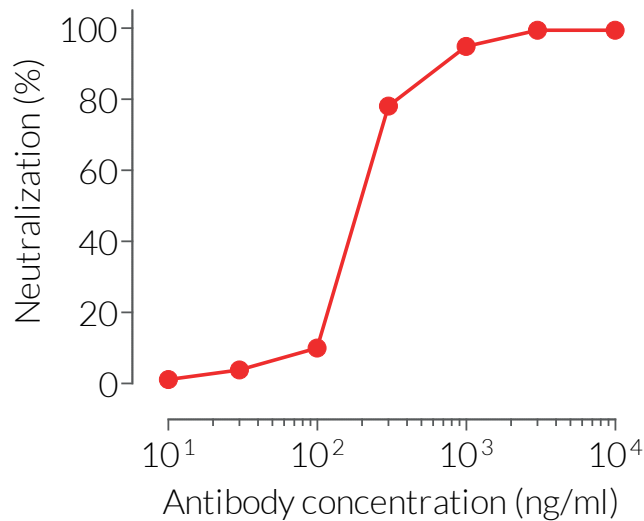
<https://www.invivogen.com/anti-mil1b-igg>

For research use only

Version 23A06-AK

Anti-mIL-1 $\beta$ -mIgG1 InvivoFit™ is a recombinant mouse monoclonal antibody (mAb) designed to efficiently neutralize mIL-1 $\beta$  *in vitro*. IL-1 $\beta$  is a soluble pro-inflammatory cytokine that plays a critical role in the host's response to infection and injury. This antibody features a 100% murine constant and variable region. Its ability to neutralize the biological activity of mIL-1 $\beta$  has been confirmed using InvivoGen's HEK-Blue™ IL-1 $\beta$  cells (Figure 1).

## Dose-dependent neutralization of mIL-1 $\beta$ signaling



**Figure 1. Anti-mIL-1 $\beta$ -mIgG1 InvivoFit™ is a potent neutralization antibody.** Increasing concentrations of Anti-mIL-1 $\beta$ -mIgG1 InvivoFit™ were incubated for 30 minutes with the recombinant cytokine mIL-1 $\beta$  (50 ng/ml) prior to the addition of HEK-Blue™ IL-1 $\beta$  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)