

Validation data for Anti-mIL-1 β -mIgG1

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

For research use only

Version 22H24-AK

Anti-mIL-1 β -mIgG1 is a recombinant mouse monoclonal antibody (mAb) designed to efficiently neutralize mIL-1 β *in vitro*. IL-1 β 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 β has been confirmed using InvivoGen's HEK-Blue™ IL-1 β cells (Figure 1).

Dose-dependent neutralization of mIL-1 β signaling

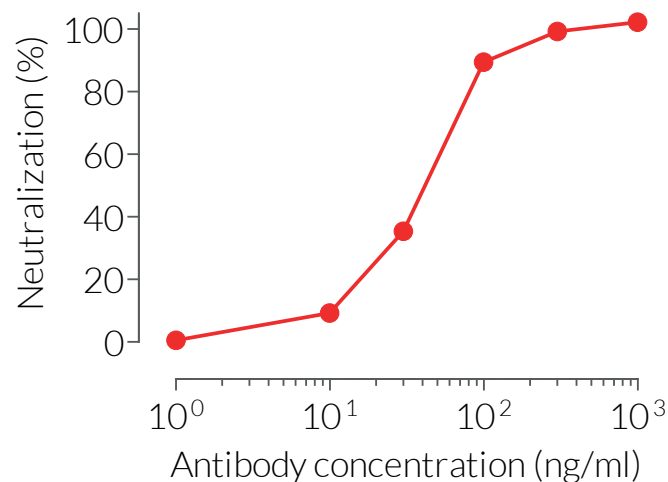


Figure 1. Anti-mIL-1 β -mIgG1 is a potent neutralization antibody. Increasing concentrations of anti-mIL-1 β -mIgG1 were incubated for 30 minutes with the recombinant cytokine mIL-1 β (10 ng/ml) prior to the addition of HEK-Blue™ IL-1 β cells. After overnight incubation, SEAP activity in the cell culture 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