

# Validation data for Anti-hIL-1 $\alpha$ -hIgG

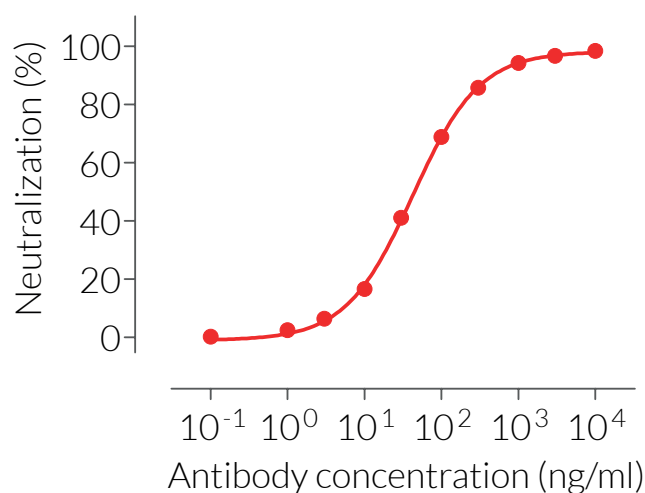
<https://www.invivogen.com/anti-hil1a-igg>

For research use only

Version 23B21-MM

Anti-hIL-1 $\alpha$ -IgG is a monoclonal antibody (mAb) designed to efficiently neutralize human interleukin 1 alpha (hIL-1 $\alpha$ ) *in vitro*. IL-1 $\alpha$  is a pro-inflammatory cytokine that plays a central role in the regulation of immune responses and inflammation. The ability of Anti-hIL-1 $\alpha$ -IgG to neutralize the biological activity of hIL-1 $\alpha$  has been confirmed using InvivoGen's HEK-Blue™ IL-1 $\beta$  cells which respond to both human IL-1 $\alpha$  and IL-1 $\beta$  (Figure 1).

## Dose-dependent neutralization of hIL-1 $\alpha$ signaling



**Figure 1. Anti-hIL-1 $\alpha$ -IgG is a potent neutralization antibody.** Increasing concentrations of Anti-hIL-1 $\alpha$ -IgG were incubated for 30 minutes with the recombinant cytokine hIL-1 $\alpha$  (1 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)

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)