

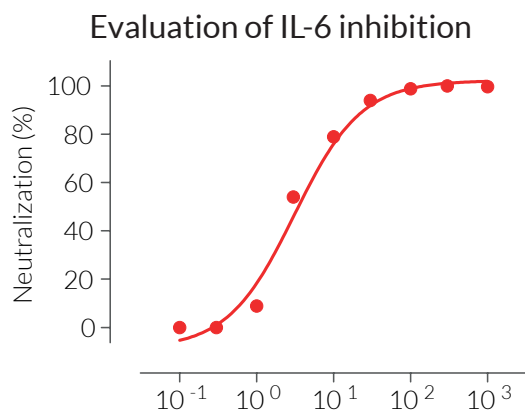
# Validation data for Anti-hIL-6-IgG

<https://www.invivogen.com/anti-hil6-igg>

For research use only

Version 23B13-MM

Anti-hIL-6-IgG is a neutralizing monoclonal antibody against human interleukin 6 (hIL-6). Neutralizing activity of this antibody was determined using recombinant hIL-6 and HEK-Blue™ IL-6 cells. These HEK293-derived cells allow the detection of bioactive human IL-6 by monitoring the expression of a STAT3-inducible SEAP (secreted embryonic alkaline phosphatase) reporter gene.



**Dose-dependent inhibition of IL-6 response using Anti-hIL-6-IgG neutralizing antibody in HEK-Blue™ IL-6 cells.** Anti-hIL-6-IgG was incubated with 1 ng/ml of hIL-6 for 30 minutes prior to the addition of HEK-Blue™ IL-6 cells. After overnight incubation, SEAP activity 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)