

Validation data for Anti-mIL-6-mIgG1e3 InvivoFit™

<https://www.invivogen.com/anti-mil6-igg1e3>

For research use only

Version 22J11-AK

Anti-mIL-6-mIgG1e3 InvivoFit™ is a recombinant mouse monoclonal antibody (mAb) designed to efficiently neutralize mIL-6. IL-6 is a multifunctional cytokine that plays an important role in the regulation of the immune response, hematopoiesis, and inflammation. This recombinant mouse mAb consists of the variable region of rat anti-mIL-6 mAb and the engineered murine IgG1e3 constant region. Its ability to neutralize the biological activity of mIL-6 has been confirmed (Figure 1).

Dose-dependent neutralization of mIL-6 signaling

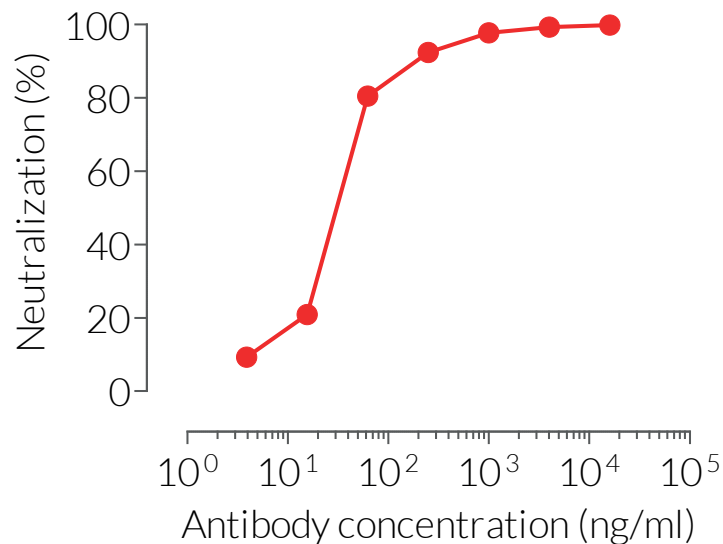


Figure 1. Anti-mIL-6-mIgG1e3 InvivoFit™ is a potent neutralization antibody. Increasing concentrations of Anti-mIL-6-mIgG1e3 InvivoFit™ were incubated for 30 minutes with the recombinant cytokine mIL-6 (30 ng/ml) prior to the addition of HEK-Lucia™ cells, which stably express the mIL-6R. After overnight incubation, Lucia activity in the cell culture supernatant was assessed using QUANTI-Luc™ detection reagent. 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