

Validation data for Anti-mIL-28b-mIgG1 InvivoFit™

<https://www.invivogen.com/anti-mil28b-igg>

For research use only

Version 22J18-AK

Anti-mIL-28b-mIgG1 InvivoFit™ is a recombinant monoclonal antibody (mAb) designed to efficiently neutralize mIL-28B. Interleukin 28B (IL-28B) is a member of the type III IFN family, playing an important role in antiviral and antitumor activity. The ability of Anti-mIL-28b-mIgG1 InvivoFit™ to neutralize the biological activity of mIL-28B has been confirmed using InvivoGen's HEK-Blue™ IFN- α/β cells (Figure 1).

Dose-dependent neutralization of mIL-28B signaling

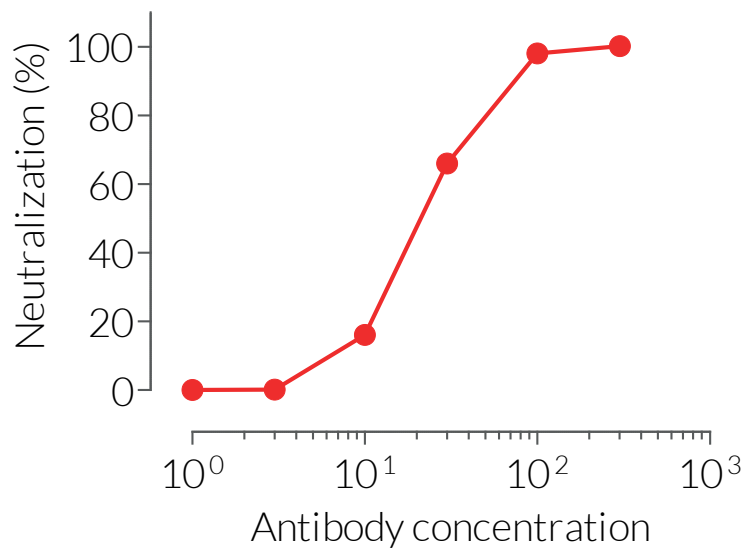


Figure 1. Anti-mIL-28b-mIgG1 InvivoFit™ is a potent neutralization antibody. Increasing concentrations of Anti-mIL-28b-mIgG1 InvivoFit™ were incubated for 30 minutes with the recombinant cytokine mIL-28B (30 ng/ml) prior to the addition of HEK-Blue™ IFN- α/β cells. After overnight incubation, SEAP activity 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