

Validation data for Anti-mTLR5-IgG

<https://www.invivogen.com/anti-mtlr5-igg>

For research use only

Version 23D04-MM

Anti-mTLR5-IgG is a monoclonal antibody (mAb) designed to efficiently neutralize murine Toll-like receptor 5 (mTLR5) *in vitro*. TLR5 recognizes flagellin from Gram-positive and Gram-negative bacteria. Activation of the receptor stimulates the production of proinflammatory cytokines. The ability of Anti-mTLR5-IgG to neutralize the biological activity of mTLR5 has been confirmed using InvivoGen's HEK-Blue™ mTLR5 cells (Figure 1).

Dose-dependent neutralization of mTLR5 signaling

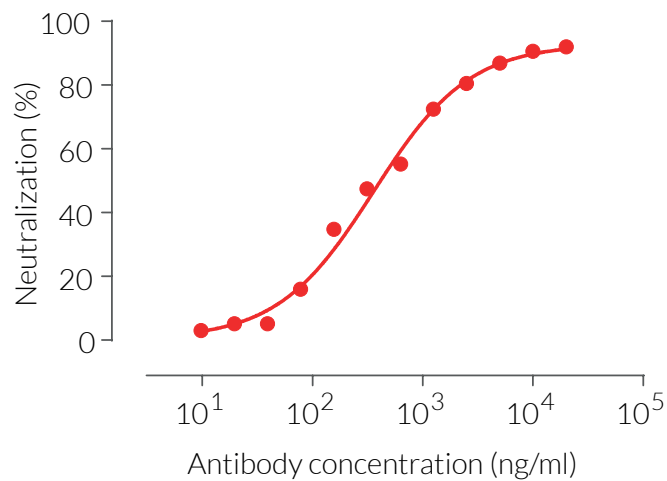


Figure 1. Anti-mTLR5-IgG is a potent neutralization antibody. Increasing concentrations of Anti-mTLR5-IgG were incubated for 30 minutes with RecFLA-ST (30 pg/ml) prior to the addition of HEK-Blue™ mTLR5 cells. After overnight incubation, secreted embryonic alkaline phosphatase (SEAP) activity in the 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