

# Validation data for Anti-hTLR5-IgA

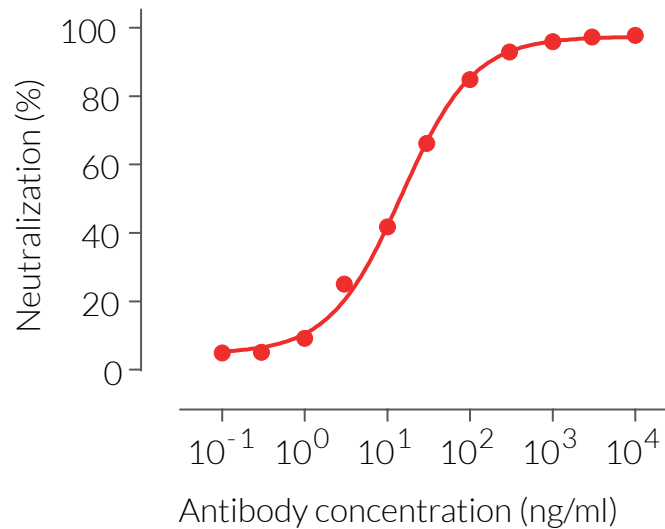
<https://www.invivogen.com/anti-htlr5-iga>

For research use only

Version 23C13-MM

Anti-hTLR5-IgA is a monoclonal antibody (mAb) designed to efficiently neutralize human Toll-like receptor 5 (hTLR5) *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-hTLR5-IgA to neutralize the biological activity of hTLR5 has been confirmed using InvivoGen's HEK-Blue™ hTLR5 cells (Figure 1).

## Dose-dependent neutralization of hTLR5 signaling



**Figure 1. Anti-hTLR5-IgA is a potent neutralization antibody.** Increasing concentrations of Anti-hTLR5-IgA were incubated for 30 minutes with FLA-ST Ultrapure (1 ng/ml) prior to the addition of HEK-Blue™ hTLR5 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](mailto:info@invivogen.com)