

# Validation data for Anti-hTLR2-IgA2

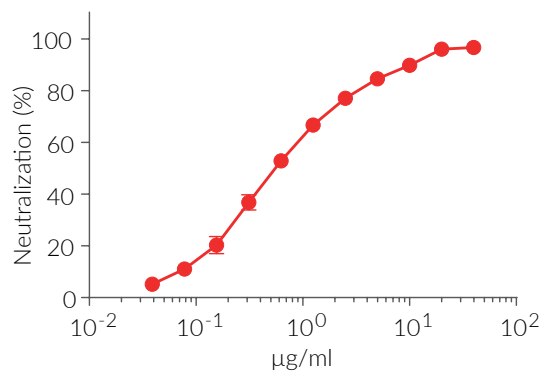
<https://www.invivogen.com/anti-htr2-iga>

For research use only

Version 24H06-NJ

Anti-hTLR2-IgA2 is a recombinant monoclonal antibody (mAb) designed to efficiently neutralize human Toll-like receptor 2 (hTLR2) *in vitro*. TLR2 is involved in the recognition of a wide array of microbial molecules. Activation of the receptor stimulates the production of proinflammatory cytokines. The ability of Anti-hTLR2-IgA2 to neutralize the biological activity of hTLR2 is validated using InvivoGen's HEK-Blue™ hTLR2 cells (Figure 1).

## Dose-dependent neutralization of hTLR2 signaling



**Figure 1. Anti-hTLR2-IgA2 is a potent neutralization antibody.** Increasing concentrations of Anti-hTLR2-IgA2 were incubated for 1 hour with HEK-Blue™ hTLR2 cells prior to the addition of 1 ng/ml FSL-1, a synthetic diacylated lipoprotein and TLR2 agonist. After overnight incubation, SEAP activity in the cell culture supernatant was assessed using QUANTI-Blue™ Solution. Data are shown in percentage (%) of neutralization (mean ± SEM).

### 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)