

Validation data for Anti-hTLR6-IgG

<https://www.invivogen.com/anti-htlr6-igg>

For research use only

Version 23C30-MM

Anti-hTLR6-IgG is a monoclonal antibody (mAb) designed to efficiently neutralize human Toll-like receptor 6 (hTLR6) *in vitro*. TLR6 co-operates with TLR2 to recognize diacylated lipoproteins (LPs) found in mycoplasma such as FSL-1. Recognition of diacylated LP by the TLR2/TLR6 heterodimer induces a signaling cascade leading to AP-1 and NF- κ B activation and cytokine production. The ability of Anti-hTLR6-IgG to neutralize the biological activity of hTLR6 has been confirmed using InvivoGen's HEK-Blue™ hTLR2 cells (Figure 1).

Dose-dependent neutralization of hTLR6 signaling

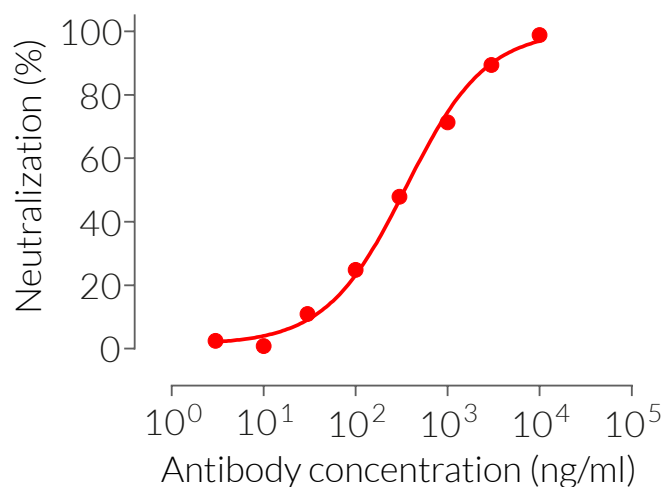


Figure 1. Anti-hTLR6-IgG is a potent neutralization antibody. Increasing concentrations of Anti-hTLR6-IgG were incubated for 30 minutes with FSL-1 (0.1 ng/ml) prior to the addition of HEK-Blue™ hTLR2 cells. After overnight incubation, secreted embryonic alkaline phosphatase (SEAP) activity in the cell 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