

# Validation data for Anti-hTNF- $\alpha$ -hIgA2

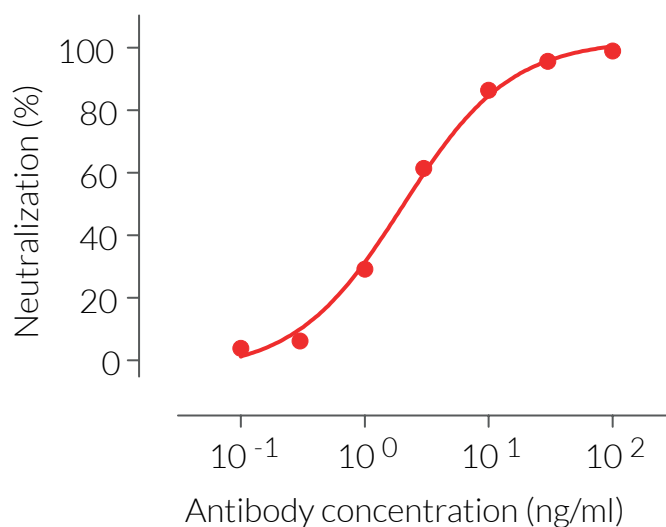
<https://www.invivogen.com/anti-htnfa-higa2>

For research use only

Version 23C01-MM

Anti-hTNF- $\alpha$ -IgA2 is a recombinant monoclonal antibody (mAb) designed to efficiently neutralize human necrosis factor-alpha (hTNF- $\alpha$ ) *in vitro*. TNF- $\alpha$  is a pleiotropic cytokine that notably induces NF- $\kappa$ B-mediated production of pro-inflammatory cytokines. The ability of Anti-hTNF- $\alpha$ -IgA2 to neutralize the biological activity of hTNF- $\alpha$  has been confirmed using InvivoGen's HEK-Blue™ TNF- $\alpha$  cells (Figure 1).

## Dose-dependent neutralization of hTNF- $\alpha$ signaling



**Figure 1. Anti-hTNF- $\alpha$ -IgA2 is a potent neutralization antibody.** Increasing concentrations of Anti-hTNF- $\alpha$ -IgA2 were incubated for 30 minutes with the recombinant cytokine hTNF- $\alpha$  (30 pg/ml) prior to the addition of HEK-Blue™ TNF- $\alpha$  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](mailto:info@invivogen.com)