

Validation data for Anti-hIFN- γ -IgA2

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

For research use only

Version 24H06-NJ

Anti-hIFN- γ -IgA2 is a recombinant monoclonal antibody (mAb) designed to efficiently neutralize human (h)IFN- γ *in vitro*. IFN- γ plays a role in activating lymphocytes to enhance anti-microbial and anti-tumor effects. The ability of Anti-hIFN- γ -IgA2 to neutralize the biological activity of hIFN- γ has been confirmed using InvivoGen's HEK-Blue™ IFN- γ cells (Figure 1).

Dose-dependent neutralization of hIFN- γ signaling

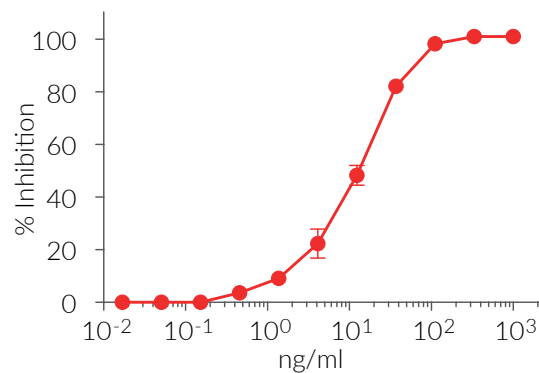


Figure 1. Anti-hIFN- γ -IgA2 is a potent neutralization antibody. Increasing concentrations of Anti-hIFN- γ -IgA2 were incubated for 30 minutes with the recombinant cytokine hIFN- γ (1 ng/ml) prior to the addition of HEK-Blue™ IFN- γ cells. After overnight incubation, SEAP activity in the cell culture supernatant was assessed using QUANTI-Blue™ Solution. Data are shown in percentage (%) of neutralization (mean \pm 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