

# Validation data for Anti-hIFN- $\gamma$ -IgA2 (H7WM120)

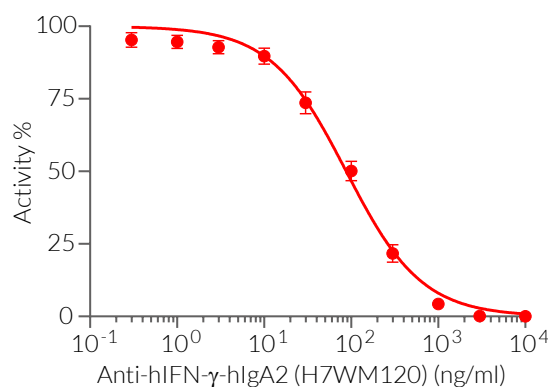
<https://www.invivogen.com/anti-hifng-iga>

For research use only

Version 25C20-NJ

Anti-hIFN- $\gamma$ -IgA2 (H7WM120) is a recombinant monoclonal antibody (mAb) designed to efficiently neutralize human (h)IFN- $\gamma$  *in vitro*. IFN- $\gamma$  plays a role in activating lymphocytes to enhance anti-microbial and anti-tumor effects. The ability of Anti-hIFN- $\gamma$ -IgA2 (H7WM120) to neutralize the biological activity of hIFN- $\gamma$  is confirmed using InvivoGen's HEK-Blue™ IFN- $\gamma$  cells (**Figure 1**).

## Dose-dependent neutralization of hIFN- $\gamma$ signaling



**Figure 1. Anti-hIFN- $\gamma$ -IgA2 (H7WM120) is a potent neutralization antibody.** Increasing concentrations of Anti-hIFN- $\gamma$ -IgA2 (H7WM120) were incubated for 30 minutes with recombinant human IFN- $\gamma$  (1 ng/ml) prior to the addition of HEK-Blue™ IFN- $\gamma$  cells. After overnight incubation, SEAP activity in the cell culture supernatant was assessed using QUANTI-Blue™ Solution. Data are shown in percentage of activity (%) (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](mailto:info@invivogen.com)