

Validation data for Anti-hIFN- α -hIgG

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

For research use only

Version 23112-AK

Anti-hIFN- α -IgG is a recombinant monoclonal antibody (mAb) designed to efficiently neutralize hIFN- α *in vitro*. Type I interferons, in particular IFN- α and IFN- β , play a vital role in host resistance to viral infections. The ability of Anti-hIFN- α -IgG to neutralize the biological activity of hIFN- α 2b has been confirmed using InvivoGen's HEK-Blue™ IFN- α / β cells (Figure 1).

Dose-dependent neutralization of hIFN- α 2b signaling

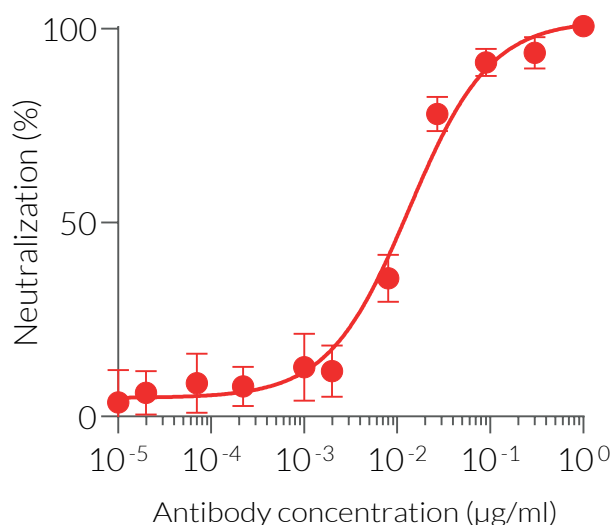


Figure 1. Anti-hIFN- α -IgG is a potent neutralization antibody. Increasing concentrations of Anti-hIFN- α -IgG were incubated for 30 minutes with the recombinant cytokine hIFN- α 2b (500 U/ml) prior to the addition of HEK-Blue™ IFN- α / β cells. After overnight incubation, SEAP activity in the cell culture supernatant was assessed using QUANTI-Blue™ Solution. Data are presented as percentage of neutralization (mean).

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