

Validation data for Anti-hIFN- β -IgG

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

For research use only

Version 23C01-MM

Anti-hIFN- β -IgG is a recombinant monoclonal antibody (mAb) designed to efficiently neutralize human interferon beta (hIFN- β) *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- β -IgG to neutralize the biological activity of hIFN- β has been confirmed using InvivoGen's HEK-Blue™ IFN- α/β cells (Figure 1).

Dose-dependent neutralization of hIFN- β signaling

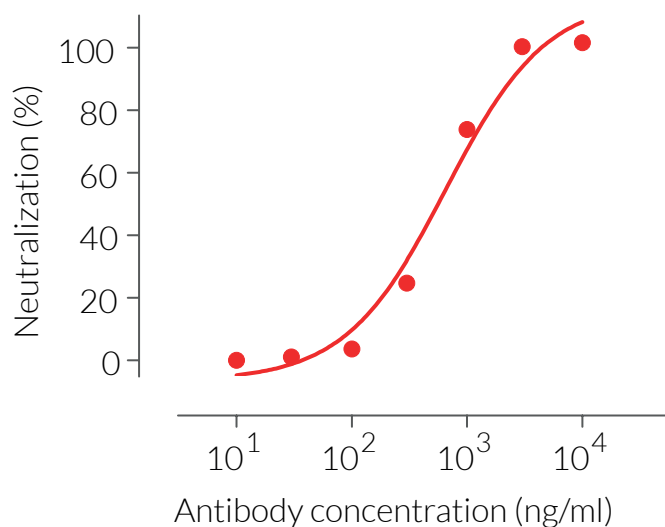


Figure 1. Anti-hIFN- β -IgG is a potent neutralization antibody. Increasing concentrations of Anti-hIFN- β -IgG were incubated for 30 minutes with the recombinant cytokine hIFN- β (0.3 U/ml) prior to the addition of HEK-Blue™ IFN- α/β cells. After overnight incubation, SEAP activity in the cell culture supernatant was assessed using QUANTI-Blue™ Solution. Data are 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