

Validation data for recombinant human IFN- α 2b

<https://www.invivogen.com/human-ifna2b>

For research use only

Version 24L13-NJ

The interferon-alpha (IFN- α) family comprises 12 members playing a non-redundant role in anti-viral and anti-tumor immune responses. The size and purity of IFN- α 2b was assessed using SDS-PAGE (Figure 1). The biological activity of IFN- α 2b is confirmed using HEK-Blue™ IFN- α / β cells featuring an ISGF3-inducible SEAP reporter (Figure 2). Recombinant human IFN- α 2b can be used together with HEK-Blue™ IFN- α / β cells for the screening of inhibitory molecules, such as Anifrolumab, a therapeutic monoclonal antibody targeting the IFNAR1 subunit of the IFN- α receptor (Figure 3).

Human IFN- α 2b SDS-PAGE analysis

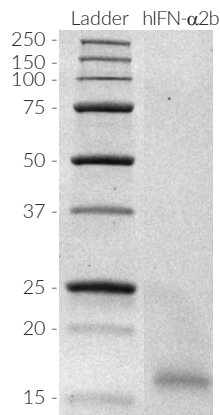


Figure 1. SDS-PAGE analysis of the recombinant human (h)IFN- α . 1 μ g of recombinant human (h)IFN- α 2b was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gel (Bio-Rad). Detection was performed as per the manufacturer's instructions. The band is detected at ~17 kDa.

Cellular response to human IFN- α 2b

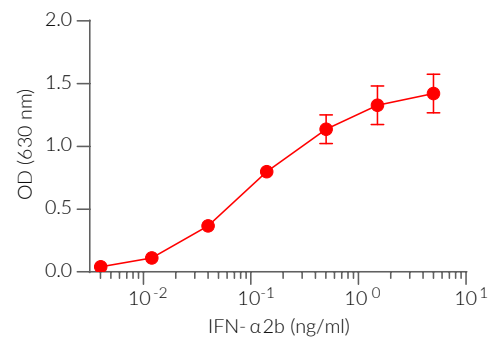


Figure 2. Dose-response of HEK-Blue™ IFN- α / β cells to recombinant human IFN- α 2b. Cells were stimulated with increasing concentrations of recombinant human (h)IFN- α 2b. After overnight incubation, the ISGF3-induced SEAP activity was determined using QUANTI-Blue™, a SEAP detection reagent. Data is shown as optical density (OD) at 630 nm (mean \pm SEM).

Dose-dependent neutralization of hIFN- α signaling

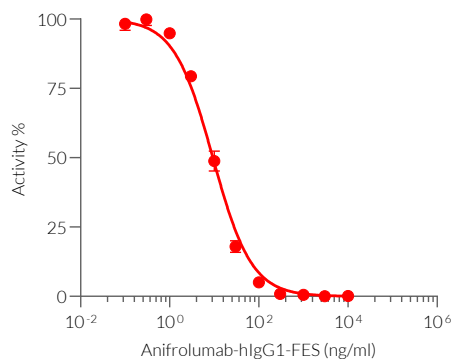


Figure 3. Anti-hIFNAR-hIgG1-FES is a potent neutralization antibody. HEK-Blue™ IFN- α / β cells were incubated with increasing concentrations of Anti-hIFNAR-hIgG1-FES (0.1 ng/ml - 10 μ g/ml) for 1 h prior to the addition of recombinant human IFN- α 2b (100 pg/ml). After overnight incubation, SEAP activity 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

 **InvivoGen**
www.invivogen.com