

Validation data for recombinant human IFN- α

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

For research use only

Version 23K02-AK

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 1, IFN- α 2b, and IFN- α 10 were assessed using SDS PAGE (Figure 1). Their biological activities have been confirmed using InvivoGen's HEK-Blue™ IFN- α / β cells (Figure 2). Binding of each IFN- α subtype to the surface receptor of HEK-Blue™ IFN- α / β cells triggers a signaling cascade leading to the activation NF- κ B and the subsequent production of SEAP. Levels of SEAP can be easily measured using QUANTI-Blue™, a detection reagent.

Detection by SDS-PAGE

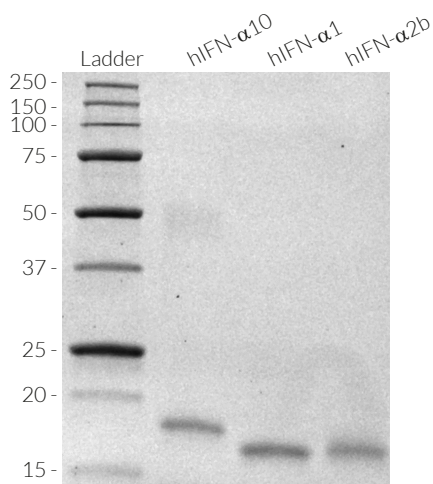


Figure 1. SDS PAGE of the recombinant human (h)IFN- α cytokines. 1 μ g of recombinant human (h)IFN- α 10, hIFN- α 1, and hIFN- α 2b were loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gel (Bio-Rad). Detection was performed as per the manufacturer's instructions. The bands were detected at ~17 - 19 kDa.

Dose-response in HEK cells to human IFN- α subtypes

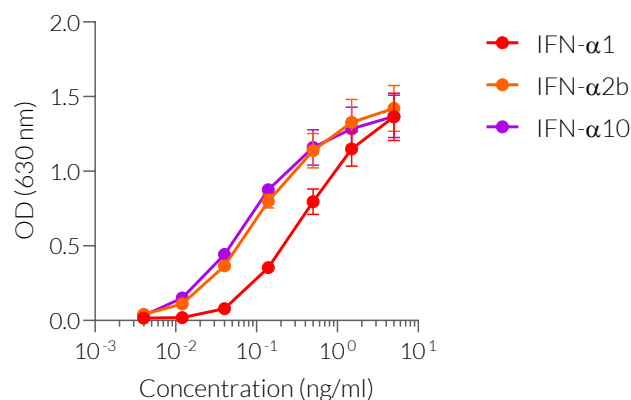


Figure 2. Dose-response in HEK-Blue™ IFN- α / β cells to recombinant hIFN- α cytokines. Cells were stimulated with increasing concentrations of recombinant human (h)IFN- α 1, hIFN- α 2b, and hIFN- α 10. After overnight incubation, the NF- κ B-induced SEAP activity was determined using QUANTI-Blue™, a SEAP detection reagent. Data are shown as optical density (OD) at 650 nm (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