

# Validation data for recombinant human IFN- $\gamma$

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

For research use only

Version 24L17-AK

Interferon-gamma (IFN- $\gamma$ ) plays a role in activating lymphocytes to enhance anti-microbial and anti-tumor effects. Its size and purity was assessed using SDS-PAGE (Figure 1). The biological activity of IFN- $\gamma$  is confirmed using HEK-Blue™ IFN- $\gamma$  cells featuring a STAT1-inducible SEAP reporter (Figure 2).

Human IFN- $\gamma$  SDS-PAGE analysis

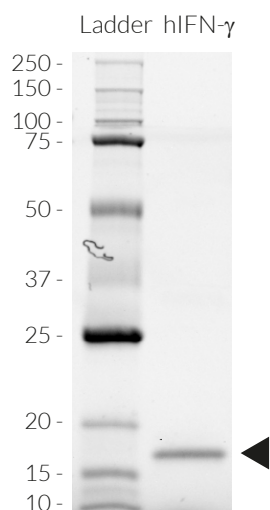


Figure 1. SDS-PAGE analysis of the recombinant human (h)IFN- $\gamma$ . 1.5  $\mu$ g of IFN- $\gamma$  was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gel (Bio-Rad). Detection was performed as per the manufacturer's instructions. A band was detected at ~17 kDa.

Cellular response to human IFN- $\gamma$

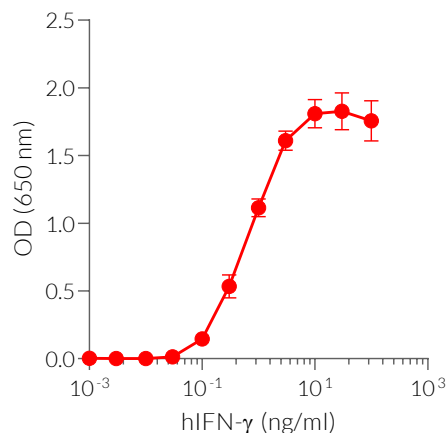


Figure 2. Dose-response of HEK-Blue™ IFN- $\gamma$  cells to human recombinant IFN- $\gamma$ . Cells were stimulated with increasing concentrations of recombinant human (h)IFN- $\gamma$ . After overnight incubation, the STAT1-induced SEAP activity was determined using QUANTI-Blue™, a SEAP detection reagent. 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](mailto:info@invivogen.com)