

# Validation data for recombinant human IFN- $\alpha$ 10

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

For research use only

Version 23A16-NJ

Recombinant human interferon alpha 10 (hIFN- $\alpha$  C) is one subset of the IFN- $\alpha$  family. It binds to the heterodimeric receptor IFNAR1/R2 and triggers the JAK1/TYK2 and ISGF3 (STAT1/STAT2/IRF9) pathway, inducing the expression of various interferon stimulated genes (ISGs). The ability of InvivoGen's recombinant hIFN- $\alpha$ 10 to trigger the IRF-dependent transcription pathway was validated using THP1-Dual™ reporter cells expressing the Lucia luciferase under the control of an ISG promoter (Figure 1).

## Cellular response to recombinant human IFN- $\alpha$ 10

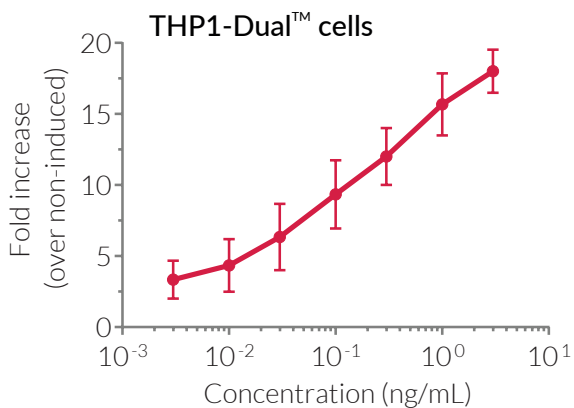


Figure 1. Dose-response of IRF-reporter cells to recombinant IFN- $\alpha$ 10.

THP1-Dual™ cells were stimulated with increasing concentration of recombinant human IFN- $\alpha$ 10. After overnight incubation, ISG activation was assessed by measuring Lucia luciferase activity in the supernatant using QUANTI-Luc™ detection reagent. Data is shown as 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)