

Validation data for recombinant human IFN- α 1

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

For research use only

Version 23A16-NJ

Recombinant human interferon alpha 1 (hIFN-alpha D) is one subset of the IFN- α 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- α 1 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- α 1

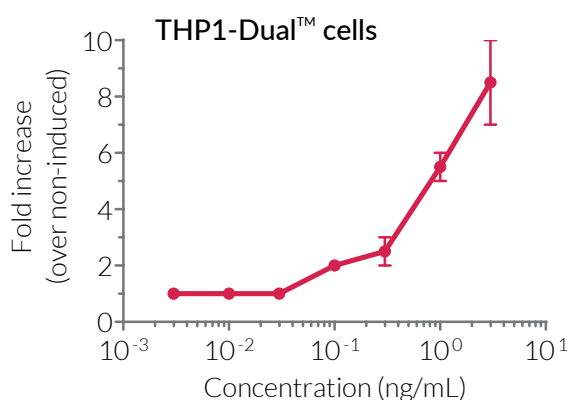


Figure 1. Dose-response of IRF-reporter cells to recombinant IFN- α 1.

THP1-Dual™ cells were stimulated with increasing concentration of recombinant human IFN- α 1. 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