

Validation data for recombinant human IFN- α 2b

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

For research use only

Version 22B28-NJ

Recombinant human interferon alpha 2b (hIFN- α 2b) 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 2b to trigger the IRF-dependent transcription pathway was validated using two reporter cells expressing secreted embryonic alkaline phosphatase (SEAP) (Figure 1A) or Lucia luciferase (Figure 1B), under the control of an ISG promoter.

Cellular response to recombinant human IFN- α 2b

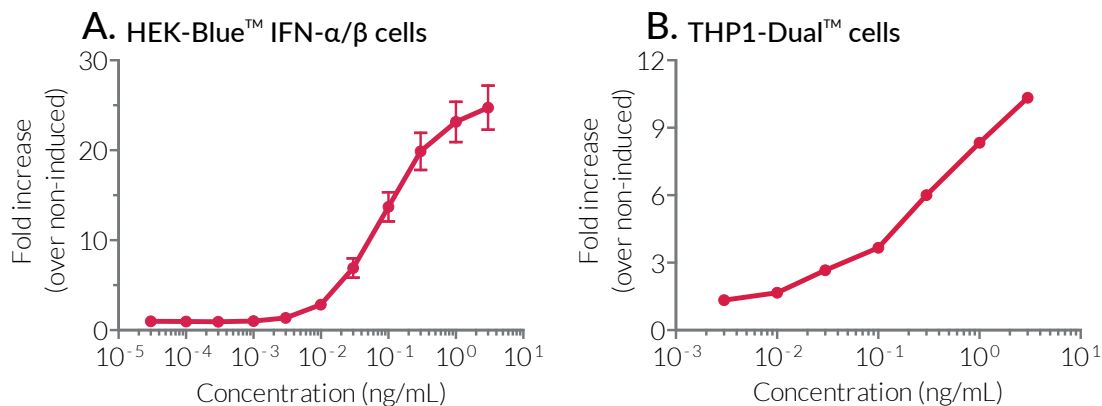


Figure 1. Dose-response of IRF-reporter cells to recombinant IFN- α 2 β .

HEK-Blue™ IFN- α / β cells (A) or THP1-Dual™ cells (B) were stimulated with increasing concentrations of recombinant human IFN- α 2b. After overnight incubation, ISG activation was assessed by measuring SEAP activity (A) or Lucia luciferase activity (B) in the supernatant using QUANTI-Blue™ Solution and QUANTI-Luc™ detection reagents, respectively.

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