

# Validation data for HEK-Blue™ IFN-λ cells

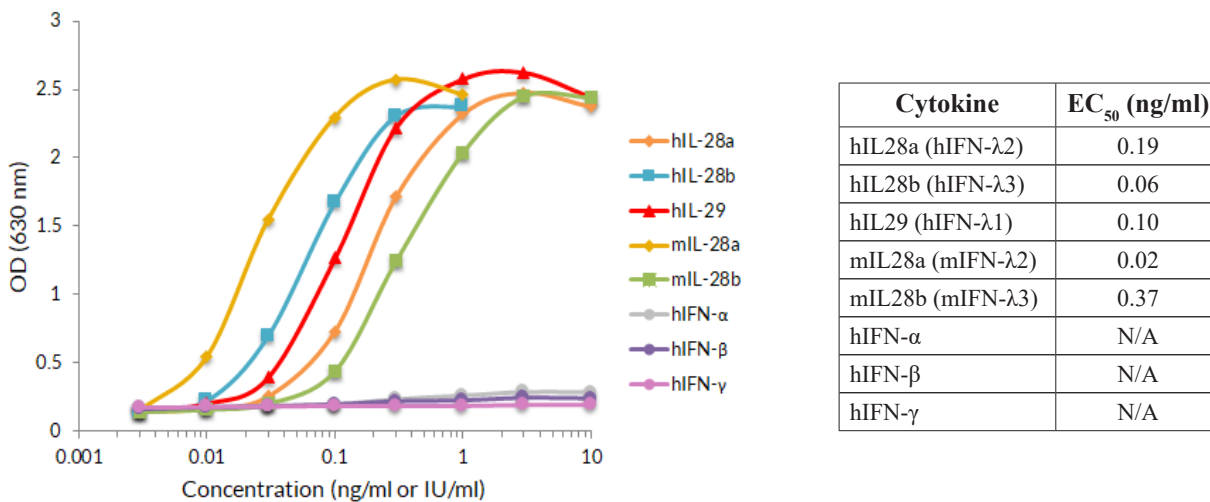
<http://www.invivogen.com/hek-blue-ifnl>

For research use only

Version 18H24-MM

HEK-Blue™ IFN-λ cells allow the detection of bioactive interferon lambda (IFN-λ; also known as type III IFNs) by monitoring the activation of the JAK/STAT/ISGF3 pathway. These cells were generated by the stable transfection of HEK293 cells with the human IFNLR and IL10R receptor genes, along with the human STAT2 and IRF9 genes to obtain a fully active IFN-λ signaling pathway. They also express a SEAP reporter gene under the control of the IFN-inducible ISG54 promoter. The response to type I, II, and III IFNs was tested. As expected, human and murine interleukin-28a (IL-28a; IFN-λ2) and IL-28b (IFN-λ3) and human IL-29 (IFN-λ1) induced the production of SEAP. These cells do not respond to type I and type II IFNs due to IFNAR2 and IFNGR1 receptor gene knockouts.

## Response of HEK-Blue™ IFN-λ cells to IFNs



**Figure 1. Response of HEK-Blue™ IFN-λ cells to type I, II and III IFNs:** HEK-Blue™ IFN-λ cells were incubated with increasing concentrations of recombinant human or mouse IFN-λ (ng/ml) or human IFN-α, IFN-β or IFN-γ (IU/ml). After 24h incubation, ISG activation was assessed by measuring SEAP levels in the supernatant using Quanti-Blue™. EC<sub>50</sub> is indicated for each cytokine (N/A: non applicable).

### 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 Hong Kong: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)