

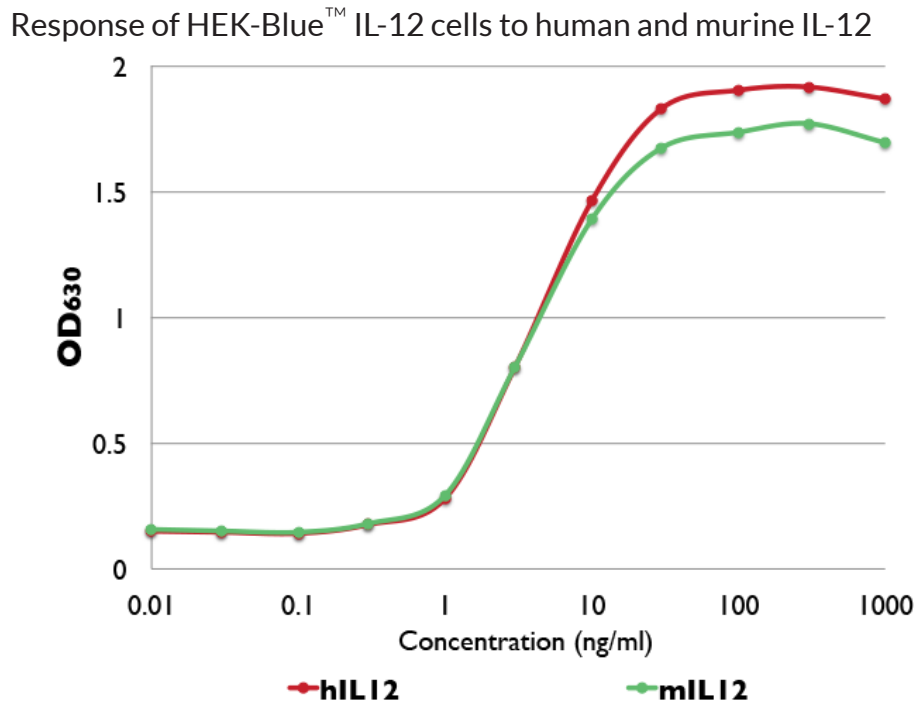
# Validation data for HEK-Blue™ IL-12 cells

<https://www.invivogen.com/hek-blue-il12>

For research use only

Version 15L03-MM

HEK-Blue™ IL-12 cells are designed to detect bioactive human and murine IL-12 by monitoring the activation of the STAT-4 pathway. They were generated by stably introducing the human genes for the IL-12 receptor and the genes of the IL-12 signaling pathway into HEK293 cells. Furthermore, these cells express a STAT4-inducible SEAP reporter gene. Stimulation of HEK-Blue™ IL-12 cells with human or murine IL-12 triggers the activation of the STAT4-inducible promoter and the production of SEAP (see figure 1).



**Figure 1:** Response of HEK-Blue™ IL-12 cells to increasing concentrations of human and murine recombinant IL-12. After a 24h incubation, the levels of SEAP were determined using QUANTI-Blue™ Solution, a SEAP detection reagent, and by reading the optical density (OD) at 630 nm.

Ligand	EC50	Response ratio
Human IL-12	4.2 ng/ml	13
Murine IL-12	4.1 ng/ml	11

The response ratio was calculated by dividing the OD at 630 nm for the treated cells by the OD at 630 nm for the untreated cells.

## 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)