

Validation data for recombinant human IL-1 β

For research use only

Version # 16A05-MM

Recombinant human interleukin-1 beta (IL-1 β) is intended for use in cell culture applications. Recombinant human IL-1 β is produced in *Escherichia coli*. The biological activity of this cytokine has been confirmed using HEK-Blue™ IL-1 β cells. They express an NF- κ B/AP-1-inducible SEAP reporter gene, which enables the detection of bioactive IL-1 β by monitoring the activation of the NF- κ B and AP-1 pathways. Binding of IL-1 β to its receptor IL-1R on the surface of HEK-Blue™ IL-1 β cells triggers a signaling cascade leading to the activation NF- κ B and the subsequent production of SEAP.

Response of HEK-Blue™ IL-1 β cells to recombinant human IL-1 β

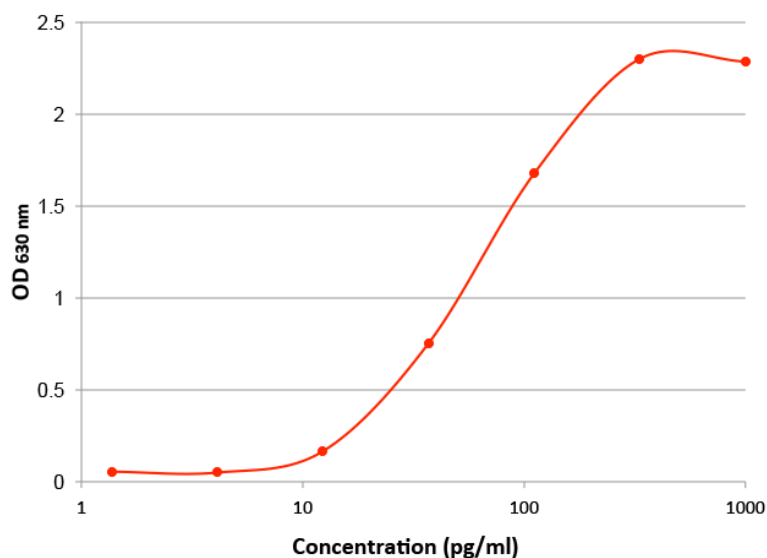


Figure 1: HEK-Blue™ IL-1 β cells were stimulated with increasing concentrations of recombinant human IL-1 β . After a 24h incubation, the levels of SEAP were determined using QUANTI-Blue™, a SEAP detection reagent, and by reading the optical density (OD) at 630 nm.

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 3-622-34-80

E-mail: info@invivogen.com