

Validation data for recombinant human IL-4

<https://www.invivogen.com/human-il4>

For research use only

Version 23J27-AK

Interleukin 4 (IL-4) is a cytokine that plays an important role in regulating inflammation and immune responses. Its size and purity was assessed using SDS PAGE (**Figure 1**). The biological activity of IL-4 has been confirmed using InvivoGen's HEK-Blue™ IL-4/IL-13 cells (**Figure 2**). Binding of IL-4 to its receptor on the surface of HEK-Blue™ IL-4/IL-13 cells triggers a signaling cascade leading to the activation NF- κ B and the subsequent production of SEAP. Levels of SEAP can be easily measured using QUANTI-Blue™, a detection reagent.

Detection by SDS-PAGE

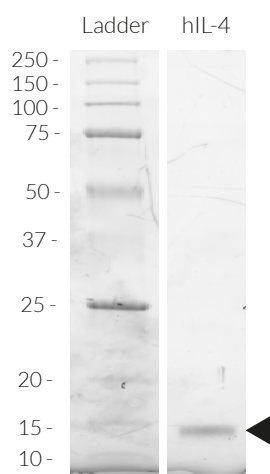


Figure 1. SDS PAGE of the recombinant human (h)IL-4 protein. 1.5 μ g of hIL-4 was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gel (Bio-Rad). Detection was performed as per the manufacturer's instructions. A band was detected at ~15 kDa.

Dose-response in HEK cells to human IL-4

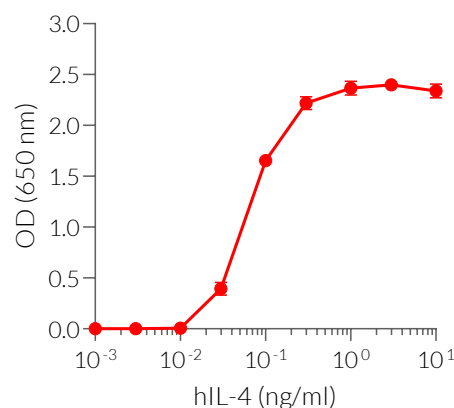


Figure 2. Dose-response in HEK-Blue™ IL-4/IL-13 cells to recombinant IL-4 cytokine. Cells were stimulated with increasing concentrations of recombinant human IL-4. After overnight incubation, the NF- κ B-induced SEAP activity was determined using QUANTI-Blue™, a SEAP detection reagent. Data are shown as optical density (OD) at 650 nm (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