

Validation data for recombinant human IL-2

Mammalian cell-expressed human interleukin 2 (IL-2)

<https://www.invivogen.com/human-il2>

For research use only

Version 24I03-AK

Interleukin 2 (IL-2) is a pleiotropic cytokine showing immunostimulatory or immunoinhibitory activity. InvivoGen offers a recombinant human IL-2, produced in mammalian Chinese hamster ovary (CHO) cells. The size and purity of hIL-2 was assessed using SDS PAGE (Figure 1). Its predicted size is ~15 kDa. The biological activity of IL-2 has been confirmed using InvivoGen's HEK-Blue™ IL-2 cells (Figure 2). Binding of IL-2 to its receptor on the surface of HEK-Blue™ IL-2 cells triggers a signaling cascade leading to the activation of STAT5 and the subsequent production of SEAP. Levels of SEAP can be easily measured using QUANTI-Blue™, a detection reagent. IL-2 was also successfully tested for T cell expansion using InvivoGen's bispecific antibody (bsAb) CD3-CD28 [data not shown].

Detection by SDS-PAGE

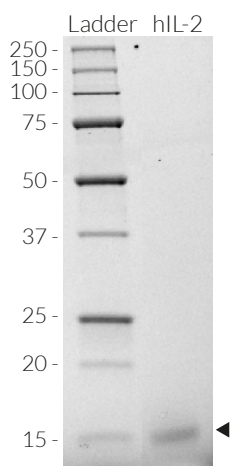


Figure 1. SDS PAGE of the recombinant human (h)IL-2 protein. 1.5 µg of hIL-2 was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gel (Bio-Rad). Detection was performed as per the manufacturer's instructions.

Dose-response in HEK cells to human IL-2

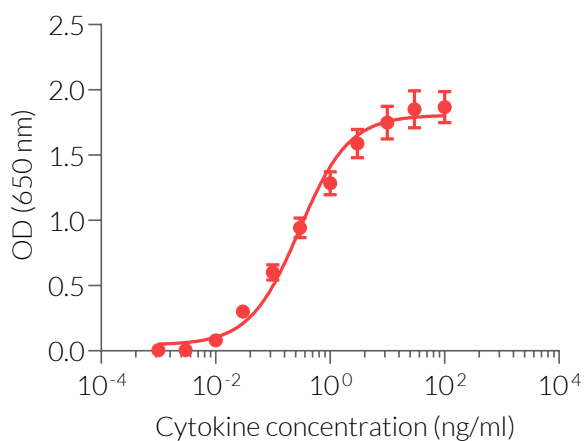


Figure 2. Dose-response in HEK-Blue™ IL-2 cells to recombinant IL-2 cytokine. Cells were stimulated with increasing concentrations of recombinant human IL-2. After overnight incubation, the STAT5-induced SEAP activity was determined using QUANTI-Blue™, a SEAP detection reagent. Data are shown as optical density (OD) at 650 nm (mean ± 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