

# Validation data for recombinant human IL-7

Mammalian cell-expressed human interleukin 7 (IL-7)

<https://www.invivogen.com/cytokines-human-il-7-recombinant-protein>

For research use only

Version 24L17-AK

Interleukin 7 (IL-7) is a hematopoietic growth factor that is involved in lymphocyte differentiation, proliferation, and survival. InvivoGen offers a recombinant human IL-7, produced in mammalian Chinese hamster ovary (CHO) cells. The size and purity of hIL-7 was assessed using SDS PAGE (Figure 1). The gel shows a characteristic smear due to the heterogeneous glycosylation of IL-7. The biological activity of IL-7 has been confirmed using InvivoGen's HEK-Blue™ IL-7 cells featuring a STAT5-inducible SEAP reporter (Figure 2).

Human IL-7 SDS-PAGE analysis

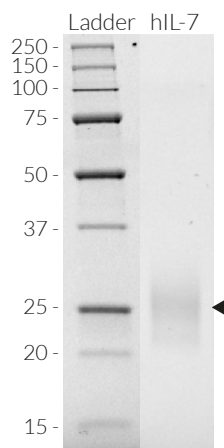


Figure 1. SDS-PAGE of the recombinant human (h)IL-7 protein. 3 µg of hIL-7 was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gel (Bio-Rad). Detection was performed as per the manufacturer's instructions.

Cellular response to human IL-7

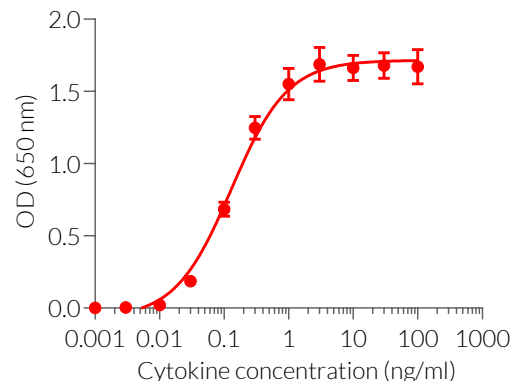


Figure 2. Dose-response of HEK-Blue™ IL-7 cells to recombinant human IL-7 cytokine. Cells were stimulated with increasing concentrations of recombinant human IL-7. After overnight incubation, the STAT5-induced SEAP activity was determined using QUANTI-Blue™, a SEAP detection reagent. Data are shown as 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](mailto:info@invivogen.com)

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)