

Validation data for recombinant human TGF- β 1

<https://www.invivogen.com/human-tgfb1>

For research use only

Version 23J27-AK

Transforming growth factor-beta (TGF- β 1) is a secreted cytokine that regulates numerous cellular functions, such as cell proliferation, apoptosis, differentiation and migration. Its size and purity was assessed using SDS PAGE (**Figure 1**). The biological activity of TGF- β 1 has been confirmed using InvivoGen's HEK-Blue™ TGF- β 1 cells (**Figure 2**). Binding of TGF- β 1 to its receptor on the surface of HEK-Blue™ TGF- β 1 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

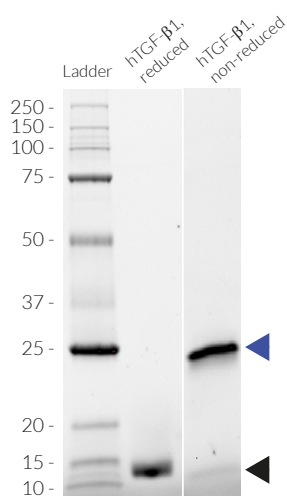


Figure 1. SDS PAGE of the recombinant human (h)TGF β 1 protein. 1.5 μ g of hTGF- β 1, reduced or non-reduced, was loaded on a 12% Mini-PROTEAN® TGX Stain-Free™ Precast Gel (Bio-Rad). Detection was performed as per the manufacturer's instructions. The bands were detected at ~15 kDa (reduced condition) and ~25 kDa (non-reduced condition).

Dose-response in HEK cells to human TGF- β 1

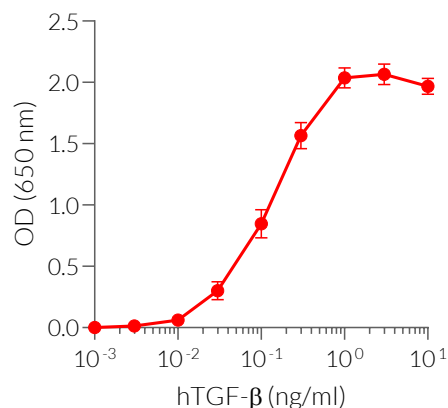


Figure 2. Dose-response in HEK-Blue™ TGF- β 1 cells to recombinant TGF- β 1 cytokine. Cells were stimulated with increasing concentrations of recombinant human TGF- β 1. After overnight incubation, the NF- κ B-induced SEAP activity was determined using QUANTI-Blue™, a SEAP detection reagent. Data are 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) 5-182-71-189-39
InvivoGen Asia: +852 3622-3480
E-mail: info@invivogen.com