

Validation data for HEK-Blue-Lucia™ TNF-α Cells

<https://www.invivogen.com/hek-blue-lucia-tnf-a>

For research use only

Version 24G05-NJ

HEK-Blue-Lucia™ TNF-α cells are designed to detect the bioactive human tumor necrosis factor-alpha (hTNF-α) by monitoring the activation of the NF-κB pathway. They express two different reporter proteins: secreted embryonic alkaline phosphatase (SEAP) and secreted Lucia luciferase. Stimulation of HEK-Blue-Lucia™ TNF-α cells with recombinant hTNF-α triggers the NF-κB-mediated production of SEAP and Lucia luciferase (figures 1-3). Of note, these cells do not respond to human IL-1β (figures 2 & 3).

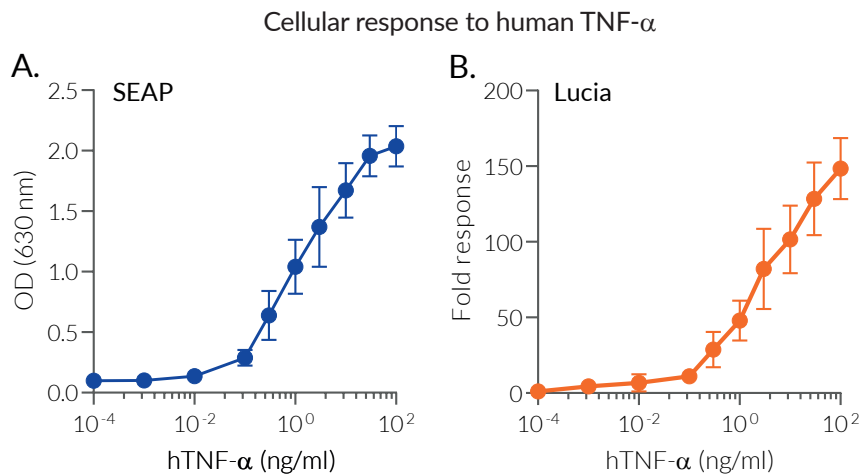


Figure 1. Dose-response of HEK-Blue-Lucia™ TNF-α cells to recombinant human TNF-α. Cells were stimulated overnight with increasing concentrations of recombinant hTNF-α. The NF-κB-induced SEAP (A) and Lucia (B) activities were assessed using QUANTI-Blue™ and QUANTI-Luc™ 4 Lucia/Gaussia detection reagents, respectively. Data are shown as optical density (OD) at 630 nm, or fold response (mean ± SEM).

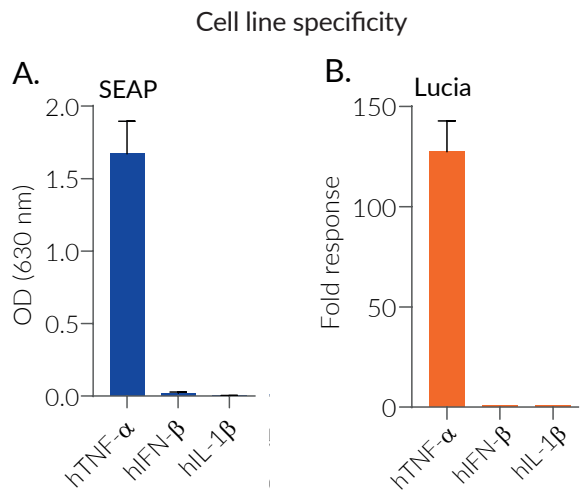


Figure 2. Response profile of HEK-Blue-Lucia™ TNF-α cells. Cells were incubated overnight with recombinant hTNF-α (10 ng/ml), hIFN-β (1000 U/ml), or hIL-1β (1 μg/ml). The NF-κB-induced SEAP (A) and Lucia (B) activities were assessed using QUANTI-Blue™ and QUANTI-Luc™ 4 Lucia/Gaussia detection reagents, respectively. Data are shown as optical density (OD) at 630 nm, or fold response (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