

Validation data for HEK-Blue™ IL-2 cells

<https://www.invivogen.com/hek-blue-il2>

For research use only

Version 21124-NJ

HEK-Blue™ IL-2 cells allow the detection of bioactive human (h) / murine (m) IL-2, and hIL-15 by monitoring the activation of the STAT-5 pathway. IL-2 and IL-15 are closely related cytokines which share IL-2R β (CD122) and IL-2R γ (CD132). HEK-Blue™ IL-2 cells were generated by the stable transfection of HEK293 cells with the human IL-2R α (CD25), IL-2R β (CD122), and IL-2R γ (CD132) genes and a STAT5-inducible SEAP (secreted embryonic alkaline phosphatase) reporter gene. These cells can detect both hIL-2 and mIL-2, and also respond to hIL-15 (**Figures 1 and 2**). They can thus be used to screen for molecules that inhibit IL-2 or IL-15 signaling such as antibodies targeting hIL-2 or hIL-15. The response of HEK-Blue™ IL-2 cells to other cytokines has been determined. Of note, a weak response to other STAT5-activating cytokines such as IFN- γ is observed (**Figure 2**).

Interestingly, a HEK-Blue-derived cell line overexpressing only CD122 and CD132 (and not CD25) responds similarly to hIL-2 and hIL-15 as the HEK-Blue™ IL-2 cell line (**Figure 3A and C**). Addition of CD25 in HEK-Blue IL-2 cells does not increase the sensitivity of these cells to hIL-2, however it does increase their sensitivity to mIL-2 (**Figure 3B**).

Cellular response to IL-2 and IL-15

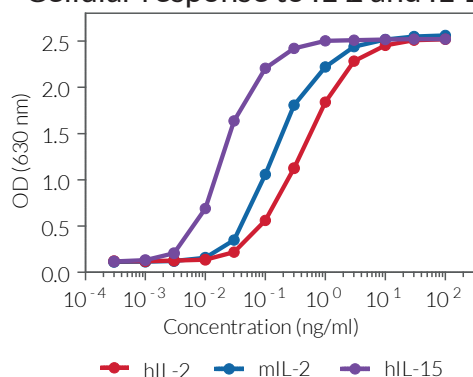


Figure 1. Dose-response of HEK-Blue™ IL-2 cells to recombinant IL-2 and IL-15. Cells were stimulated with increasing concentrations of recombinant human and murine IL-2, or human IL-15. After overnight incubation, the STAT5 response was determined using QUANTI-Blue™ Solution, a SEAP detection reagent, and reading the optical density (OD) at 630 nm.

Cell line specificity

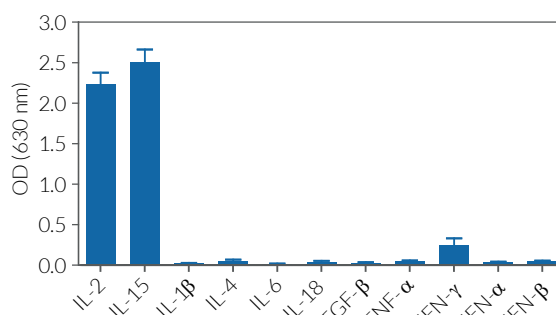


Figure 2. Cytokine response profile of HEK-Blue™ IL-2 cells. Cells were stimulated with various human recombinant cytokines; 1 ng/ml IL-2 or IL-15 and 10 ng/ml IL-1 β , IL-4, IL-6, IL-18, TGF- β , TNF- α , IFN- γ , and 1x10⁴ IU/ml IFN- α , IFN- β . After overnight incubation, SEAP activity was assessed using QUANTI-Blue™ Solution, a SEAP detection reagent, and reading the optical density (OD) at 630 nm.

HEK-Blue™ IL-2 and HEK-Blue™ CD122/CD132 response to IL-2 and IL-15

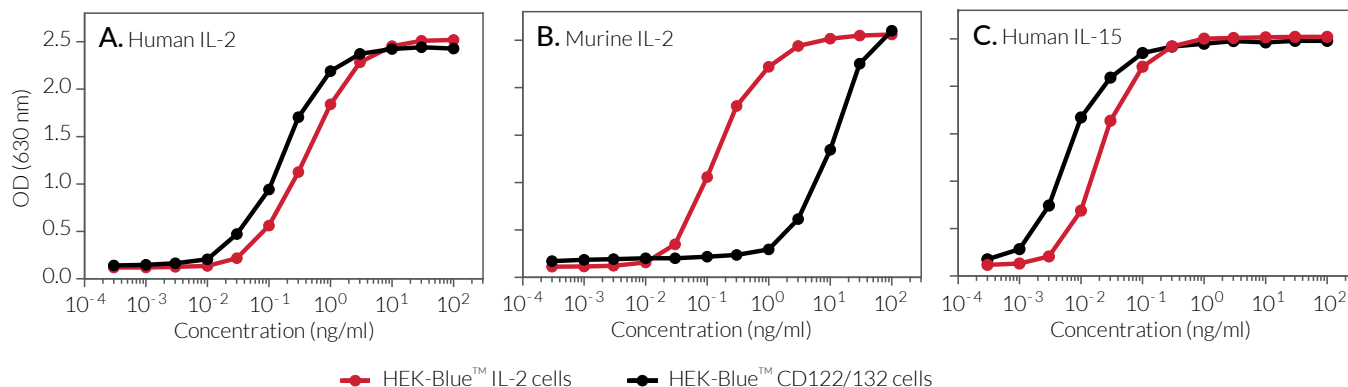


Figure 3. Dose-response of HEK-Blue™ IL-2 and HEK-Blue™ CD122/CD132 cells to recombinant IL-2 and IL-15. Cells were stimulated with increasing concentrations of recombinant human and murine IL-2, and human IL-15. After overnight incubation, the STAT5 response was determined using QUANTI-Blue™ Solution, a SEAP detection reagent, and reading the optical density (OD) at 630 nm.

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