

Validation data for HEK-Blue™ IL-19/IL-20 cells

<https://www.invivogen.com/hek-blue-il20>

For research use only

Version 23F29-AK

HEK-Blue™ IL-19/IL-20 reporter cells have been specifically designed to detect bioactive human (h) and murine (m) IL-19/IL-20, IL-22, and IL-24 by monitoring the activation of the JAK-STAT pathway. HEK-Blue™ IL-19/IL-20 cells were generated by stable transfection of the human embryonic kidney HEK293 cell line with the genes encoding human *IL-20R α* , *IL-20R β* , and *STAT3* genes to obtain a fully active IL-19/IL-20 signaling pathway, as verified by functional assays (Figure 1). Since IL-19 and IL-20 belong to the IL-10 cytokine family, the response of HEK-Blue™ IL-19/IL-20 cells to members of this cytokine family as well as other cytokines has been determined (Figure 2). The ability to block the signaling of hIL-24 in HEK-Blue™ IL-19/IL-20 cells was assessed using an anti-hIL-24 neutralization monoclonal antibody (mAb) (Figure 3).

Dose response to human and murine recombinant cytokines

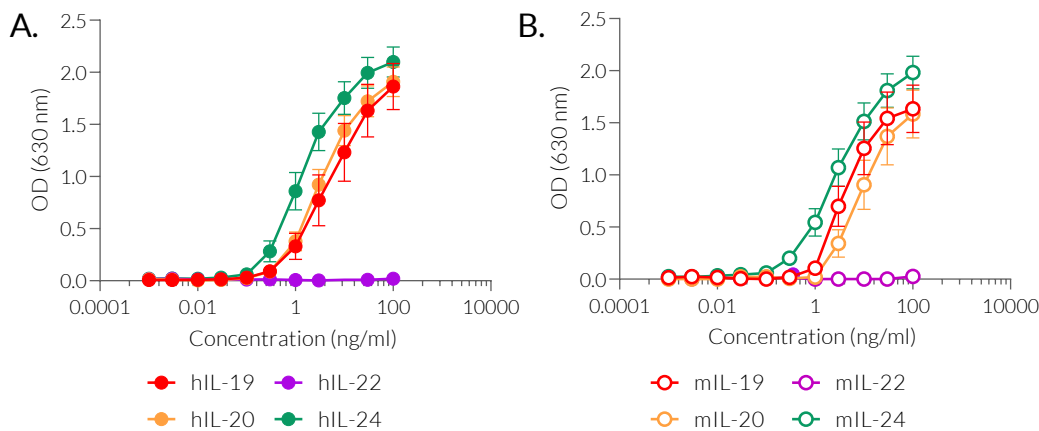


Figure 1. Dose-response of HEK-Blue™ IL-19/IL-20 cells to recombinant cytokines. Cells were stimulated with increasing concentrations of recombinant (A) human or (B) murine IL-19, IL-19/IL-20, IL-22, and IL-24. After overnight incubation, the NF- κ B-induced SEAP activity was determined using QUANTI-Blue™ Solution, a SEAP detection reagent. Data are shown as optical density (OD) at 630 nm (mean \pm SEM).

Cytokine response profile of HEK-Blue™ IL-19/IL-20 cells

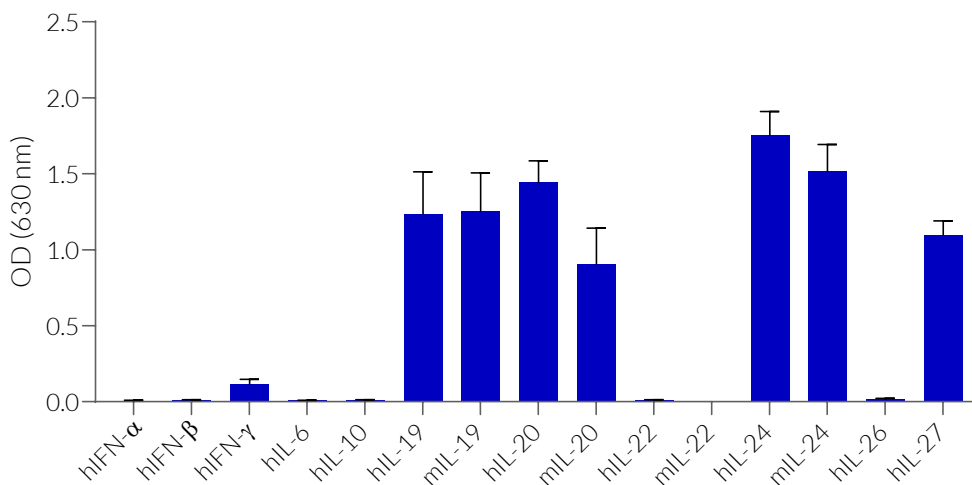


Figure 2. Cytokine response profile of HEK-Blue™ IL-19/IL-20 cells. Cells were stimulated with various human and murine recombinant cytokines: hIFN- α (1000 U/ml), hIFN- β (1000 U/ml), hIFN- γ (10 ng/ml), hIL-6 (10 ng/ml), hIL-10 (10 ng/ml), hIL-19 (10 ng/ml), mL-19 (10 ng/ml), hIL-19/IL-20 (10 ng/ml), mL-19/IL-20 (10 ng/ml), hIL-22 (10 ng/ml), mL-22 (10 ng/ml), hIL-24 (10 ng/ml), mL-24 (10 ng/ml), hIL-26 (10 ng/ml), and hIL-27 (10 ng/ml). After overnight incubation, the NF- κ B-induced SEAP activity was determined using QUANTI-Blue™ Solution. Data are shown as optical density (OD) at 630 nm (mean \pm SEM).

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
 InvivoGen USA (International): +1 (858) 457-5873
 InvivoGen Europe: +33 (0) 5-62-71-69-39
 InvivoGen Asia: +852 3622-3480
 E-mail: info@invivogen.com

Inhibition of hIL-24 signaling in HEK-Blue™ IL-19/IL-20 cells

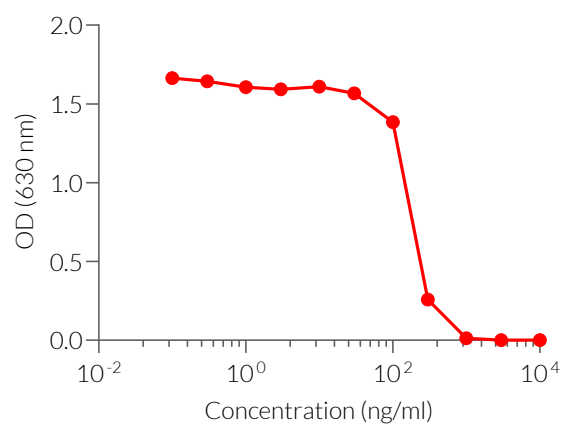


Figure 3. Inhibition of hIL-24 signaling in HEK-Blue™ IL-19/IL-20 cells. Increasing concentrations of an anti-hIL-24 neutralization mAb were incubated for 30 minutes with the recombinant cytokine hIL-24 (3 ng/ml) prior to the addition of HEK-Blue™ IL-19/IL-20 cells. After overnight incubation, SEAP activity in the cell culture supernatant was assessed using QUANTI-Blue™ Solution. Data are shown as optical density (OD) at 630 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