

Validation data for HEK-Blue™ TL1A Cells

<https://www.invivogen.com/hek-blue-tl1a>

For research use only

Version 24L18-NJ

HEK-Blue™ TL1A cells allow the detection of bioactive tumor necrosis factor (TNF)-like 1A (TL1A) by monitoring AP-1/NF- κ B activation. These human embryonic kidney 293 (HEK293)-derived cells express the human DR3 (Death Receptor 3) homotrimeric receptor for TL1A, as well as an AP-1/NF- κ B-inducible secreted embryonic alkaline phosphatase (SEAP) reporter. HEK-Blue™ TL1A cells respond to human and murine TL1A (Figure 1 and Figure 3). These cells can be used for the screening of antibodies targeting the TL1A pathway (Figure 2). Of note, HEK-Blue™ TL1A cells also respond to two other human AP-1/ NF- κ B-signaling cytokines, TNF- α and IL-1 β . However, they do not respond to other cytokines of the TNF superfamily: APRIL, BAFF, RANKL, and CD40L (Figure 3).

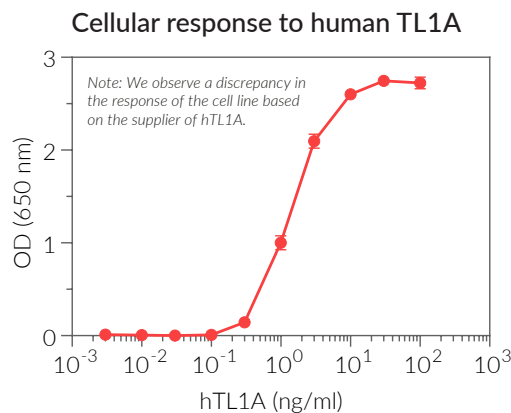


Figure 1. Dose-response of HEK-Blue™ TL1A cells to recombinant human TL1A. Cells were stimulated with increasing concentrations of recombinant human TL1A. After overnight incubation, the NF- κ B/AP1-induced SEAP activity was determined using QUANTI-Blue™, a SEAP detection reagent. Data are shown as optical density (OD) at 650 nm (mean \pm SEM).

Neutralization of hTL1A signaling using Tulsokibart

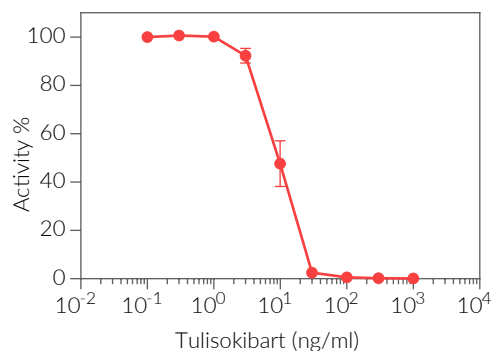


Figure 2. Dose-dependent inhibition of HEK-Blue™ TL1A cell response using Tulsokibart biosimilar. Increasing concentrations of Anti-hTL1A Tulsokibart biosimilar (0.1 ng/ml - 1 μ g/ml) were incubated with recombinant human TL1A (3 ng/ml) for 1 h before the addition of HEK-Blue™ TL1A cells. After overnight incubation, SEAP activity in the cell culture supernatant was assessed using QUANTI-Blue™ Solution. Data are shown in percentage of activity (mean \pm SEM).

Cell line specificity

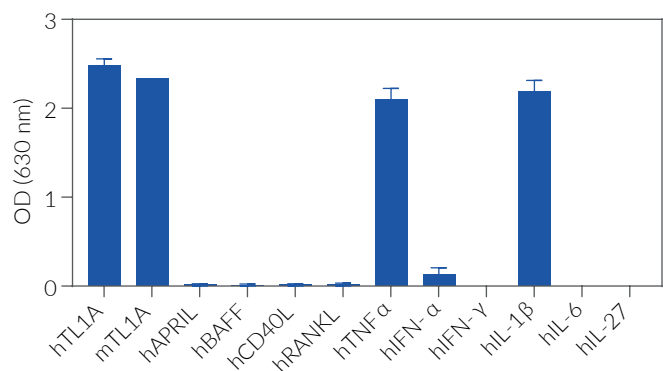


Figure 3. Response of HEK-Blue™ TL1A cells to a panel of cytokines. Cells were stimulated with various human and murine recombinant cytokines: 10 ng/ml of hTL1A, mTL1A, hAPRIL, hBAFF, hCD40L, hRANKL, hTNF- α , hIFN- γ , hIL-1 β , hIL-6, hIL-27 and 1000 U/ml hIFN- α . After overnight incubation, SEAP activity was assessed using QUANTI-Blue™. The OD at 630 nm is shown as 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



Any questions about our cell lines?
Visit our FAQ page.

 **InvivoGen**
www.invivogen.com