

Validation data for TL8-506

<https://www.invivogen.com/tl8-506>

For research use only

Version 23K30-AK

TL8-506 is a benzodiazepine compound, analog of the Toll-like receptor 8 (TLR8) agonist VTX-2337. The ability of TL8-506 to activate TLR8 signaling was validated using a panel of InvivoGen's reporter cell lines. TL8-506 efficiently activates human (h)TLR8, murine (m)TLR7 and mTLR8, but not hTLR7, as assessed by the expression of an NF- κ B-inducible secreted embryonic alkaline phosphatase (SEAP) reporter in HEK-Blue™-derived cell lines (Figure 1). The induction of the NF- κ B and IRF pathways by TL8-506 has been tested using InvivoGen's THP1-Dual™ cells featuring two reporter genes, the NF- κ B-inducible SEAP and IRF-inducible Lucia luciferase, as well as the overexpression and/or a knockout (KO) of a TLR gene (Figure 2). Due to the low level of endogenous TLR8 in THP-1-derived cells, high concentrations of TL8-506 can stimulate THP1-Dual™ and THP1-Dual™ hTLR7 cells.

Dose-dependent NF- κ B response in HEK-Blue™-derived cells

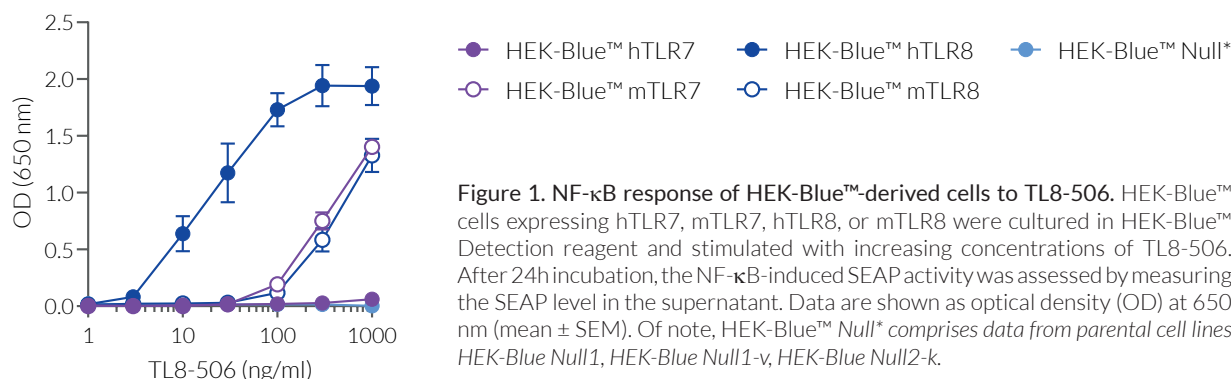


Figure 1. NF- κ B response of HEK-Blue™-derived cells to TL8-506. HEK-Blue™ cells expressing hTLR7, mTLR7, hTLR8, or mTLR8 were cultured in HEK-Blue™ Detection reagent and stimulated with increasing concentrations of TL8-506. After 24h incubation, the NF- κ B-induced SEAP activity was assessed by measuring the SEAP level in the supernatant. Data are shown as optical density (OD) at 650 nm (mean \pm SEM). Of note, HEK-Blue™ Null* comprises data from parental cell lines HEK-Blue Null1, HEK-Blue Null1-v, HEK-Blue Null2-k.

Dose-dependent NF- κ B and IRF responses in THP1-Dual™-derived cells

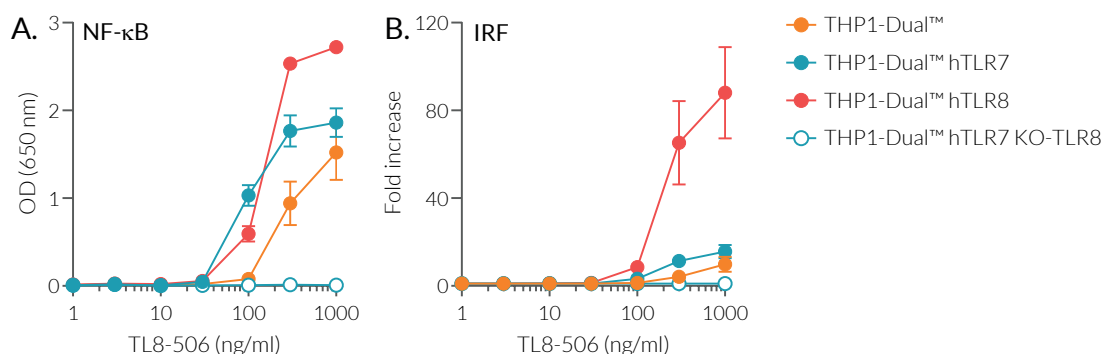


Figure 2. NF- κ B and IRF responses of THP1-Dual™-derived cells to TL8-506. THP1-Dual™, THP1-Dual™ hTLR7 cells, THP1-Dual™ hTLR8 cells, and THP1-Dual™ hTLR7 KO-TLR8 cells were incubated for 24 hours with increasing concentrations of TL8-506. After 24h incubation, the (A) NF- κ B-induced SEAP activity was assessed using QUANTI-Blue™. Data are shown as optical density (OD) at 650 nm (mean \pm SEM). (B) The IRF response was assessed by measuring the activity of Lucia luciferase in the supernatant using QUANTI-Luc™. Data are shown in fold response over non-induced cells (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