

Validation data for CL307

<https://www.invivogen.com/cl307>

For research use only

Version 23L18-AK

CL307 was generated by covalently linking spermine to the adenine analog CL264. CL307 is a potent and specific agonist of Toll-like receptor 7 (TLR7). The ability of CL307 to activate TLR7, but not TLR8 signaling was validated using a panel of InvivoGen's reporter cell lines. CL307 efficiently activates human (h) and mouse (m)TLR7, but not hTLR8, 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 CL307 has been tested using InvivoGen's THP1-Dual™-derived monocytic cell lines, that features two reporter genes, the NF- κ B-inducible SEAP and IRF-inducible Lucia luciferase, as well as the overexpression of TLR7 or TLR8 (Figure 2).

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

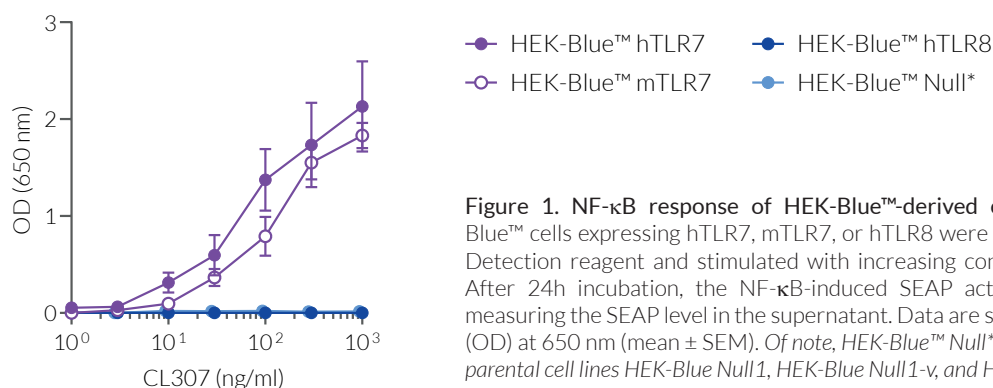


Figure 1. NF- κ B response of HEK-Blue™-derived cells to CL307. HEK-Blue™ cells expressing hTLR7, mTLR7, or hTLR8 were cultured in HEK-Blue™ Detection reagent and stimulated with increasing concentrations of CL307. After 24h incubation, the NF- κ B-induced SEAP activity was assessed by measuring the 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 the parental cell lines HEK-Blue Null1, HEK-Blue Null1-v, and HEK-Blue Null2-k.

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

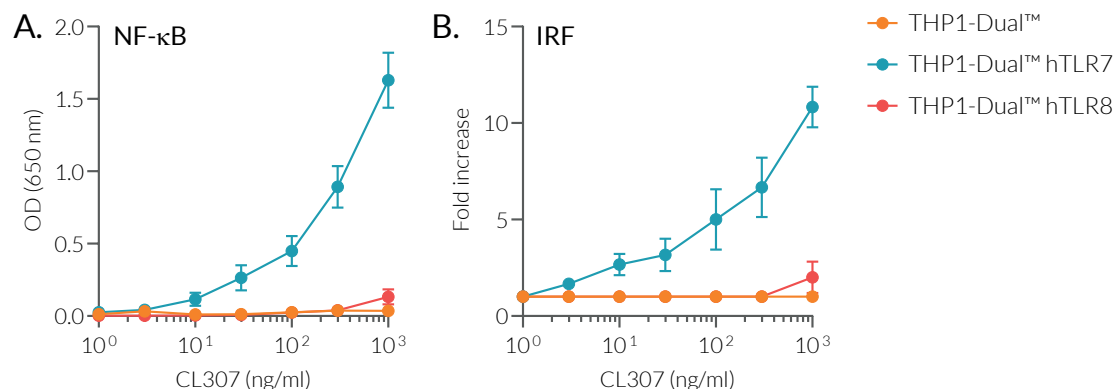


Figure 2. NF- κ B and IRF responses of THP1-Dual™-derived cells to CL307. THP1-Dual™, THP1-Dual™ hTLR7, and THP1-Dual™ hTLR8 cells were incubated for 24 hours with increasing concentrations of CL307. After 24h incubation, the (A) NF- κ B-induced SEAP activity was assessed using QUANTI-Blue™. Data are 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