

Validation data for Loxoribine

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

For research use only

Version 23L21-AK

Loxoribine is a guanosine analog derivatized at position N⁷ and C⁸. It is a potent agonist of the Toll-like receptors 7 (TLR7). The ability of loxoribine to activate TLR7 signaling was validated using a panel of InvivoGen's reporter cell lines. Loxoribine efficiently activates human (h) and mouse (m)TLR7, but not TLR8, 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).

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

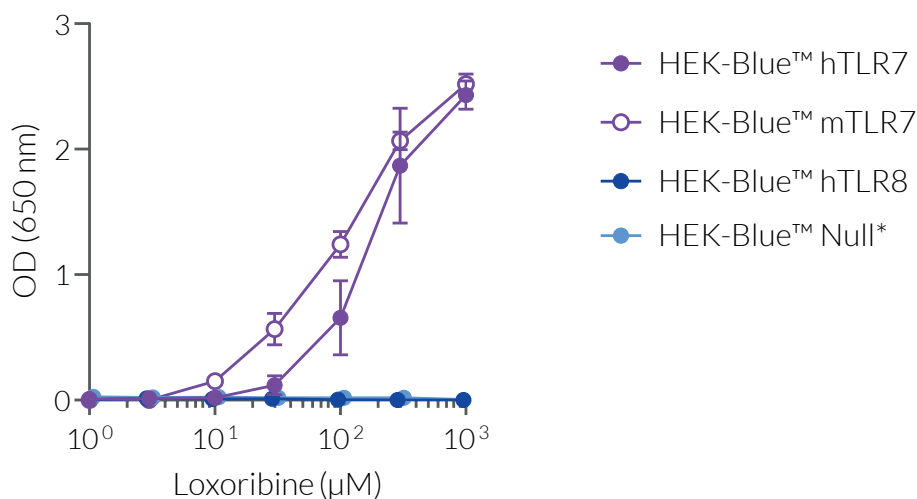


Figure 1. NF- κ B response of HEK-Blue™-derived cells to loxoribine. HEK-Blue™ cells expressing hTLR7, mTLR7, or hTLR8 were cultured in HEK-Blue™ Detection reagent and stimulated with increasing concentrations of loxoribine. After 24h incubation, the NF- κ B-induced SEAP activity was assessed by measuring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

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