

Validation data for M5049

<https://www.invivogen.com/m5049>

For research use only

Version 23D27-NJ

M5049 (also known as Enpatoran) is a small molecule that functions as a dual and selective inhibitor of TLR7 and TLR8. The ability of M5049 to inhibit TLR7 and TLR8 signaling was validated using a panel of InvivoGen's reporter cell lines. M5049 efficiently inhibits human (h)TLR7, mouse (m)TLR7, hTLR8, but not mTLR8, 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 inhibition potency of M5049 for both NF- κ B and IRF signaling pathways downstream of human TLR7 and TLR8 was also confirmed using THP1-Dual™-derived cell lines expressing inducible SEAP and Lucia luciferase reporter genes (Figure 2). The specific inhibition of TLR7 and TLR8 signaling by M5049 has been verified (Figure 3).

Inhibition of TLR7 and TLR8 signaling by M5049 in HEK-Blue™ cells

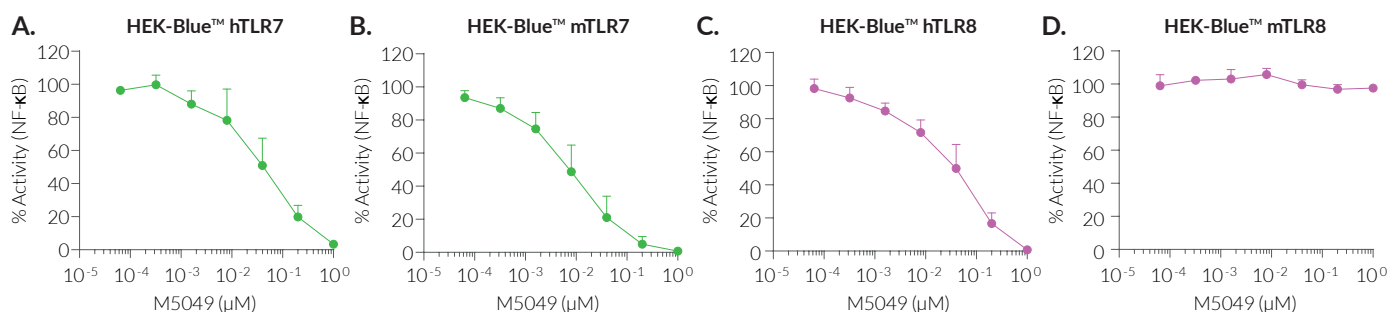


Figure 1: M5049 is a potent inhibitor of human TLR7, mouse TLR7, human TLR8, but not mouse TLR8 signaling pathways. HEK-Blue™ cells expressing hTLR7 (A), mTLR7 (B), hTLR8 (C), or mTLR8 (D), were cultured with increasing concentrations of M5049. After 3 hours of incubation, the following ligands were added: 30 ng/ml R848, a TLR7/8 agonist (A and B), or 30 ng/ml (C) or 300 ng/ml (D) TL8-506, a TLR8 agonist. After overnight incubation, the neutralizing activity of M5049 was determined by measuring the reduction of SEAP production in the supernatant using the HEK-Blue™ detection reagent. Data are shown as a percentage (%) of maximal TLR activation with each agonist (without inhibitor; mean + SEM).

Inhibition of TLR7 and TLR8 signaling by M5049 in THP1-Dual™ cells

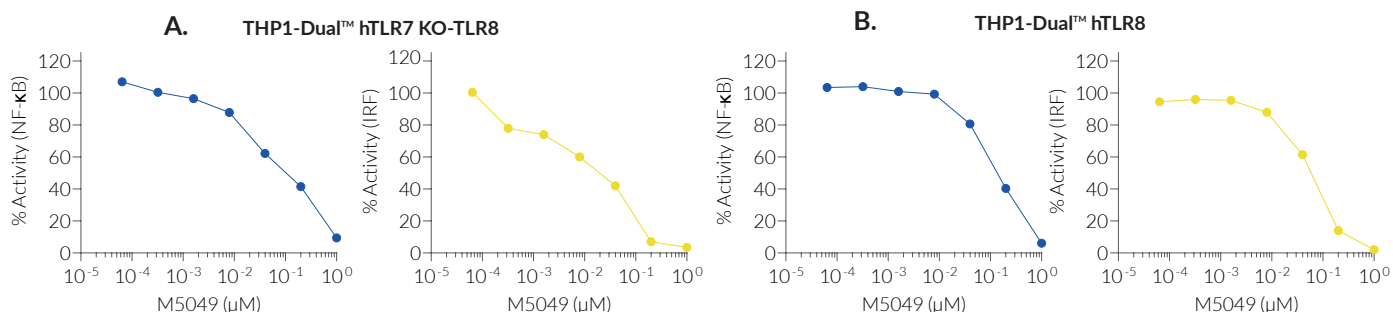


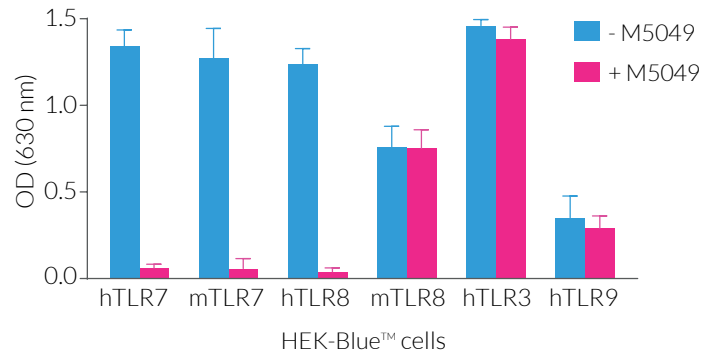
Figure 2: M5049 is a potent inhibitor of NF- κ B and IRF signaling pathways downstream of human TLR7 and TLR8. THP1-Dual™ hTLR7 KO-TLR8 (A) and THP1-Dual™ hTLR8 (B) cells were cultured in the presence of increasing concentrations of M5049. After 3 hours of incubation, the following ligands were added: 1 μ g/ml R848 (TLR7/8 agonist) (A), or 1 μ g/ml TL8-506 (TLR8 agonist) (B). After overnight incubation, the neutralizing activity of M5049 was determined by measuring the reduction of SEAP and Lucia luciferase production in the supernatant using the QUANTI-Blue™ and QUANTI-Luc™ 4 detection reagents, respectively. Data are shown as a percentage (%) of maximal TLR activation with each agonist (without inhibitor).

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

 **InvivoGen**
www.invivogen.com

Specific inhibition of TLR7 and TLR8 signaling by M5049



Abbreviations: hTLR7 = HEK-Blue™ hTLR7, mTLR7 = HEK-Blue™ mTLR7, hTLR8 = HEK-Blue™ hTLR8, mTLR8 = HEK-Blue™ mTLR8, hTLR3 = HEK-Blue™ hTLR3, hTLR9 = HEK-Blue™ hTLR9 cells.

Figure 3: M5049 is a specific dual inhibitor of TLR7 and TLR8. HEK-Blue™ cells expressing hTLR7, mTLR7, hTLR8, mTLR8, hTLR3, or hTLR9 were incubated with M5049 (1 μ M). After 3 hours of incubation, the following ligands were added: R848 30 ng/ml (hTLR7) or 100 ng/ml (mTLR7), TL8-506 30 ng/ml (hTLR8) or 300 ng/ml (mTLR8), Poly(I:C) HMW 50 ng/ml (hTLR3), LPS-EK 1 ng/ml (hTLR4), and ODN 2006 500 ng/ml (hTLR9). After overnight incubation, the neutralizing activity of M5049 was determined by measuring the reduction of SEAP production in the supernatant using the HEK-Blue™ detection reagent. 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