

# Validation data for BMS-986256

<https://www.invivogen.com/tlr-7-8-inhibitor-bms986256-afimetroan>

For research use only

Version 24J07-NJ

BMS-986256 (also known as Afimetroan) is a small molecule that functions as a dual and selective inhibitor of TLR7 and TLR8. The ability of BMS-986256 to inhibit TLR7 and TLR8 signaling was validated using a panel of InvivoGen's cell lines featuring an NF- $\kappa$ B-inducible secreted embryonic alkaline phosphatase (SEAP) reporter. BMS-986256 efficiently inhibits human (h)TLR7, mouse (m)TLR7, hTLR8, but not mTLR8, in HEK-Blue™-derived cell lines overexpressing TLR7 or TLR8 (Figure 1). The inhibition potency of BMS-986256 for both NF- $\kappa$ B and IRF signaling pathways downstream of hTLR7 and hTLR8 was also confirmed using THP1-Dual™-derived cell lines expressing NF- $\kappa$ B-inducible SEAP and IRF-inducible Lucia luciferase reporter genes (Figure 2). The specific inhibition of TLR7 and TLR8 signaling by BMS-986256 has been verified (Figure 3).

## Inhibition of TLR7 and TLR8 signaling by BMS-986256 in HEK-Blue™ cells

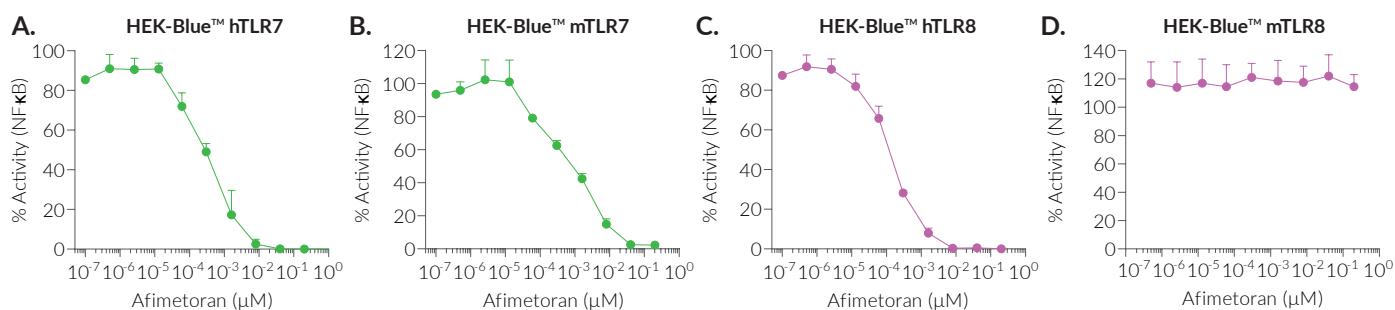


Figure 1: BMS-986256 is a potent inhibitor of human TLR7, mouse TLR7, and human TLR8, but not mouse TLR8 signaling pathways in HEK cells overexpressing TLR7 or TLR8. HEK-Blue™ cells overexpressing hTLR7 (A), mTLR7 (B), hTLR8 (C), or mTLR8 (D), were cultured with increasing concentrations of BMS-986256. After 3 hours of incubation, the following ligands were added: 100 ng/ml (A) or 300 ng/ml (B) R848, a TLR7/8 agonist, or 300 ng/ml (C) or 1 μg/ml (D) TL8-506, a TLR8 agonist. After overnight incubation, the neutralizing activity of BMS-986256 was determined by measuring the reduction of SEAP production in the supernatant using the QUANTI-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 BMS-986256 in THP1-Dual™ cells

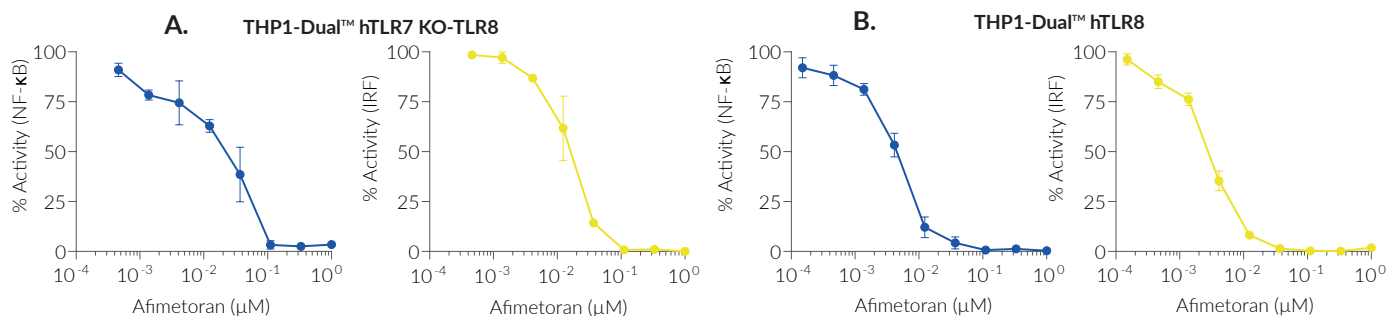


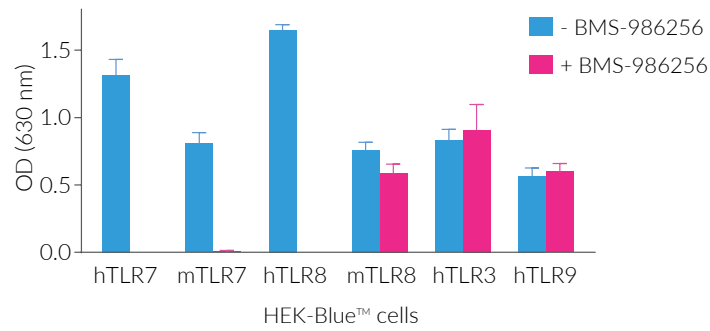
Figure 2: BMS-986256 is a potent inhibitor of NF- $\kappa$ B and IRF signaling pathways downstream of human TLR7 and TLR8. THP1-Dual™ hTLR7 KO-TLR8 (A) and THP1-Dual™ hTLR8 (B) cells were cultured in the presence of increasing concentrations of BMS-986256. After 3 hours of incubation, the following ligands were added: 1 μg/ml R848 (TLR7/8 agonist) (A), or 1 μg/ml TL8-506 (TLR8 agonist) (B). After overnight incubation, the neutralizing activity of BMS-986256 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; 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](mailto:info@invivogen.com)

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)

## Specific inhibition of TLR7 and TLR8 signaling by BMS-986256



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: BMS-986256 is a specific dual inhibitor of TLR7 and TLR8.** HEK-Blue™ cells overexpressing hTLR7, mTLR7, hTLR8, mTLR8, hTLR3, or hTLR9 were incubated with BMS-986256 (1  $\mu$ M). After 3 hours of incubation, the following ligands were added: R848 30 ng/ml (hTLR7) or 300 ng/ml (mTLR7), TL8-506 100 ng/ml (hTLR8) or 3  $\mu$ g/ml (mTLR8), Poly(I:C) HMW 50 ng/ml (hTLR3), and ODN 2006 500 ng/ml (hTLR9). After overnight incubation, the neutralizing activity of BMS-986256 was determined by measuring the reduction of SEAP production in the supernatant using QUANTI-Blue™ detection reagent. Data are shown as optical density (OD) at 630 nm (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](mailto:info@invivogen.com)