

# Validation data for HEK-Blue™ mTLR7 cells

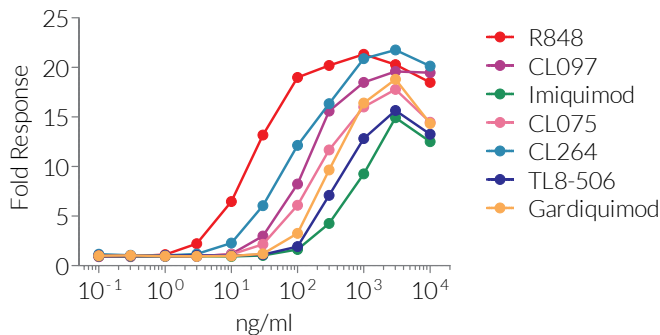
<https://www.invivogen.com/hek-blue-mltr7>

For research use only

Version 22K27-AK

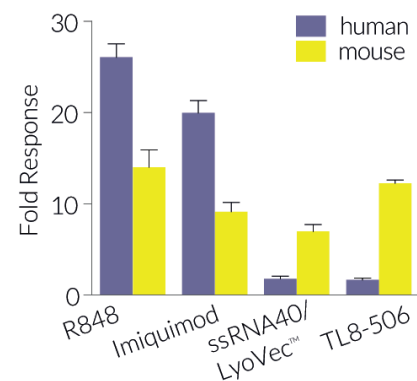
HEK-Blue™ mTLR7 cells are designed for studying the murine Toll-Like Receptor 7 (mTLR7) signaling by monitoring the activation of NF-κB/AP-1. They express the *mTLR7* gene as well as an NF-κB/AP1-inducible secreted embryonic alkaline phosphatase (SEAP) reporter gene. Levels of SEAP produced upon TLR7 activation can be easily determined in real-time with HEK-Blue™ Detection cell culture medium. HEK-Blue™ mTLR7 cells respond in a dose-dependent manner to synthetic base analogs such as TLR7/8-specific base analogs including R848 (Resiquimod), CL097, and CL075 (Figure 1). Importantly, HEK-Blue™ hTLR7 cells also respond to the TLR8-specific agonists TL8-506 and ssRNA40 (a single-stranded RNA sequence from HIV-1) (Figure 2). Of note, there are discrepancies in the functional activities between human and mouse TLR7 (Figure 2). Indeed, while HEK-Blue™ hTLR7 cells do not respond to the TLR8 agonists ssRNA40 and TL8-506, HEK-Blue™ mTLR7 cells do respond to these ligands. As HEK293 cells express endogenous levels of various pattern recognition receptors, HEK-Blue™ mTLR7 cells might respond to the cognate ligands (Figure 3).

## Cellular response to synthetic base analogs



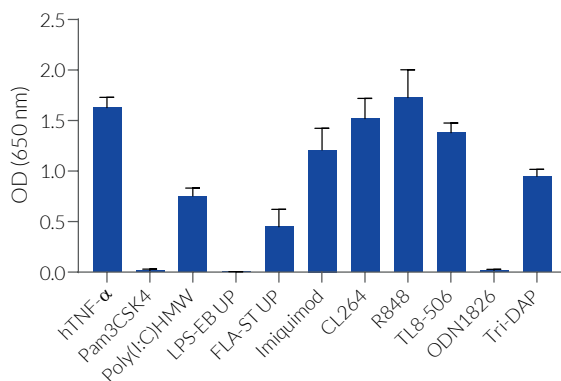
**Figure 1. Dose-response of HEK-Blue™ mTLR7 cells to synthetic base analogs.** HEK-Blue™ mTLR7 cells were cultured in HEK-Blue™ Detection medium with increasing concentrations of TLR7 agonists (CL264, Imiquimod, Gardiquimod), TLR7/8 agonists (R848, CL097, CL075), or a TLR8 agonist (TL8-506). After 24h incubation, TLR7-induced NF-κB/AP-1 responses were assessed by measuring SEAP levels in the supernatant by reading the optical density (OD) at 650 nm. OD fold increase over non-induced cells is shown.

## Human and Mouse TLR7-induced responses



**Figure 2. Species-driven TLR7 differential responses.** HEK-Blue™ hTLR7 or mTLR7 were cultured in HEK-Blue™ Detection medium with 1 μg/ml R848, 3 μg/ml Imiquimod, 5 μg/ml ssRNA40/LyoVec™ (referred as human TLR8 agonist), or 1 μg/ml TL8-506. After 24h incubation, TLR8-induced NF-κB/AP1 responses were assessed by measuring SEAP levels in the supernatant and reading the OD at 650 nm. OD fold increase over non-induced cells is shown (mean ± SEM). *Note: The HEK-Blue™ hTLR7 cell line used in this experiment is a different clone than the one on sale.*

## Response to various PRR agonists and cytokines



**Figure 3. Response of HEK-Blue™ mTLR7 cells to various PRR agonists and cytokines.** Cells were cultured in HEK-Blue™ Detection medium and incubated for 24 hours with cytokines and various TLR agonists: Human TNF-α (NF-κB-positive control, 1 ng/ml), Pam3CSK4 (TLR2 ligand, 1 μg/ml), Poly(I:C) HMW (TLR3 ligand, 1 μg/ml), LPS-EK Ultrapure (UP) (TLR4 ligand, 10 μg/ml), FLA-ST UP (TLR5 ligand, 10 μg/ml), C264 (TLR7 ligand, 1 μg/ml), R848 (TLR7/8 ligand, 1 μg/ml), ODN 2006 (TLR9 ligand, 10 μg/ml), or Tri-DAP (NOD1 ligand, 1 μg/ml). After 24h incubation, TLR8-induced NF-κB/AP1-induced SEAP activity was determined by measuring the SEAP level in the supernatant. Data are shown as OD at 650 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](mailto:info@invivogen.com)