

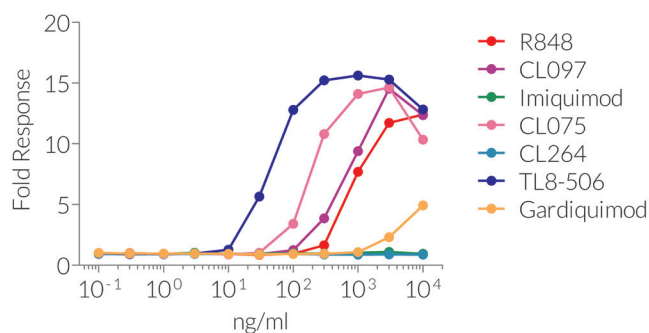
# Validation data for HEK-Blue™ hTLR8 cells

<https://www.invivogen.com/hek-blue-htlr8>

For research use only  
Version 23K27-AK

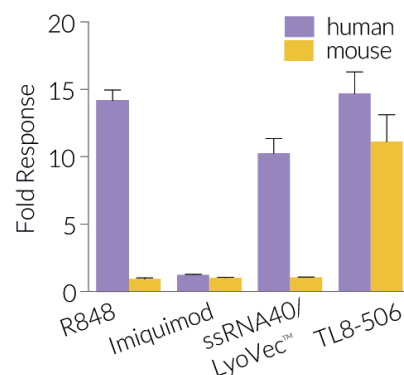
HEK-Blue™ hTLR8 cells are designed for studying the human Toll-like receptor 8 (hTLR8) signaling by monitoring the activation of NF-κB/AP1. They express the hTLR8 gene, and an NF-κB/AP1-inducible secreted embryonic alkaline phosphatase (SEAP) reporter gene. Levels of SEAP produced upon TLR8 activation can be easily determined in real-time with HEK-Blue™ Detection cell culture medium. HEK-Blue™ hTLR8 cells respond in a dose-dependent manner to synthetic base analogs, while they not respond to TLR7-specific base analogs (Figure 1). Importantly, HEK-Blue™ hTLR8 cells respond to the TLR8-specific agonist ssRNA40 (a single-stranded RNA sequence from HIV-1) and this response is potentiated by the addition of Poly(dT) (Figure 2 and data not shown). Of note, there are discrepancies in the functional activities between human and mouse TLR8 (Figure 2). As HEK293 cells express endogenous levels of various pattern recognition receptors, HEK-Blue™ hTLR8 cells might respond to the cognate ligands (Figure 3).

## Cellular response to synthetic base analogs



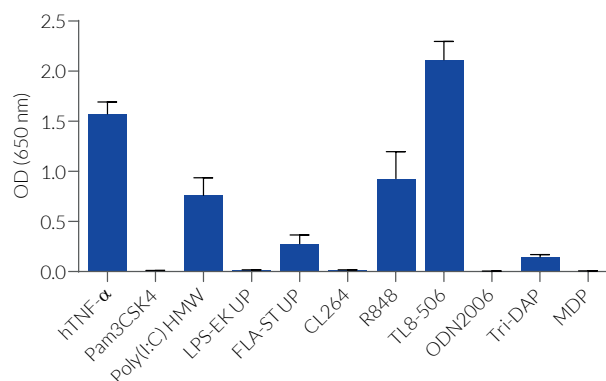
**Figure 1. Dose-response of HEK-Blue™ hTLR8 cells to synthetic base analogs.** Cells were cultured in HEK-Blue™ Detection medium with increasing concentrations of a TLR8 agonist (TL8-506), various TLR7/8 agonists (R848, CL097, CL075), or TLR7 agonists (CL264, Imiquimod, Gardiquimod). After 24h incubation, TLR8-induced NF-κB/AP1-induced SEAP activity was determined by reading the optical density (OD) at 650 nm. OD fold increase over non-induced cells is shown.

## Human and Mouse TLR8-induced responses



**Figure 2. Species-driven TLR8 differential responses.** HEK-Blue™ hTLR8 or mTLR8 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 as described before. OD fold increase over non-induced cells is shown (mean ± SEM).

## Response to various PRR agonists and cytokines



**Figure 3. Response of HEK-Blue™ hTLR8 cells to various PRR agonists and cytokines.** Cells were cultured in HEK-Blue™ Detection medium and stimulated for 24 hours with cytokines and various TLR agonists: Human TNF-α (NF-κB-positive control, 1 ng/ml), Pam3CSK4 (TLR2 ligand, 100 ng/ml), Poly(I:C) HMW (TLR3 ligand, 100 ng/ml), LPS-EK Ultrapure (UP) (TLR4 ligand, 100 ng/ml), FLA-ST UP (TLR5 ligand, 10 ng/ml), C264 (TLR7 ligand, 1 μg/ml), R848 (TLR7/8 ligand, 1 μg/ml), ODN 2006 (TLR9 ligand, 10 μg/ml), Tri-DAP (NOD1 ligand, 100 ng/ml), or MDP (NOD2 ligand, 100 ng/ml). After 24h incubation, TLR8-induced NF-κB/AP1 responses were assessed as described before. 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)