

# Validation data for HEK-Blue™ mTLR4 cells

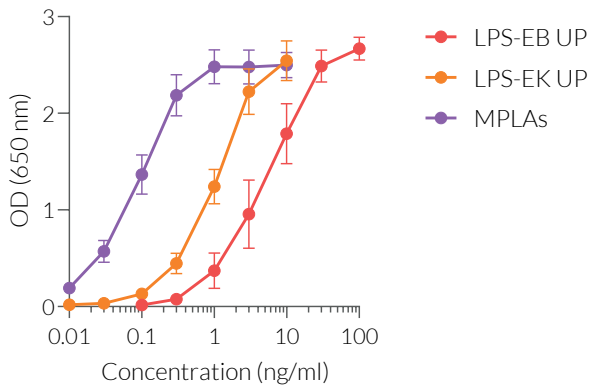
<https://www.invivogen.com/hek-blue-mtlr4>

For research use only

Version 24B29-AK

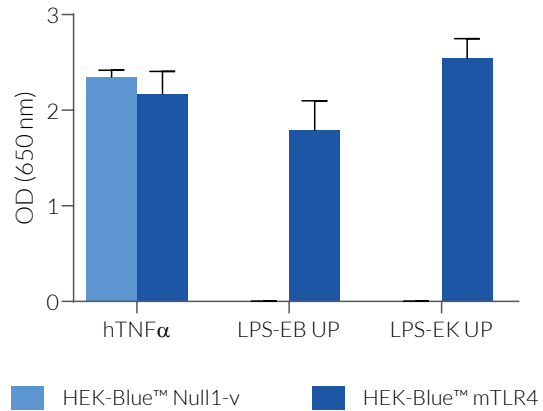
HEK-Blue™ mTLR4 cells are engineered HEK293 cells that feature the stable expression of the mouse Toll-like receptor 4 (mTLR4), as well as the adapter proteins mouse (m)MD-2 (myeloid differentiation factor 2) and mCD14 (cluster of differentiation 14). They also express an NF-κB-inducible SEAP (secreted embryonic alkaline phosphatase) reporter gene. These cells are highly responsive to TLR4 ligands, such as Lipopolysaccharide (LPS) and monophosphoryl Lipid A (MPLA) (Figure 1), when compared to their parental cell line HEK-Blue™ Null1-v (Figure 2). As HEK293 cells express endogenous levels of various pattern recognition receptors (PRRs), HEK-Blue™ mTLR4 cells respond to their cognate ligands (Figure 3).

## Dose-response of HEK-Blue™ mTLR4 cells to TLR4 agonists



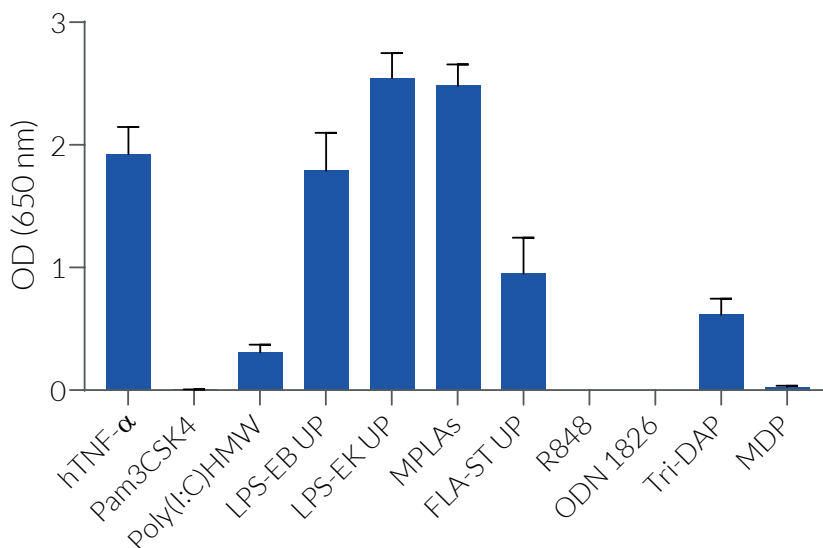
**Figure 1. Dose-response of HEK-Blue™ mTLR4 cells to TLR4 agonists.** Cells were cultured in HEK-Blue™ Detection reagent and stimulated with increasing concentrations of LPS-EB Ultrapure (UP), LPS-EK UP, and MPLAs (synthetic). After 24 hour incubation, the NF-κB-induced SEAP activity was determined by reading the OD at 650 nm (mean ± SEM).

## Response of HEK-Blue™-derived cells to TLR4 agonists



**Figure 2. Response of HEK-Blue™-derived cells to TLR4 agonists.** HEK-Blue™ Null1-v and HEK-Blue™ mTLR4 cells were cultured in HEK-Blue™ Detection reagent and incubated with 10 ng/ml of the TLR4 agonists LPS-EB Ultrapure (UP) and LPS-EK UP. Human TNF-α, (10 ng/ml) serves as an NF-κB-positive control. After 24h incubation, the NF-κB-induced SEAP activity was assessed by reading the OD at 650 nm (mean ± SEM).

## Response of HEK-Blue™ mTLR4 cells to various PRR agonists and cytokines



**Figure 3. Response of HEK-Blue™ mTLR4 cells to various PRR agonists and cytokines.** Cells were cultured in HEK-Blue™ Detection reagent and stimulated for 24 hours with cytokines and various PRR agonists: Human TNF-α (NF-κB-positive control, 10 ng/ml), Pam3CSK4 (TLR2 ligand, 100 ng/ml), Poly(I:C) HMW (TLR3 ligand, 1 μg/ml), LPS-EB UP, LPS-EK UP (TLR4 ligands, 10 ng/ml), MPLAs (synthetic TLR4 ligand, 1 ng/ml) FLA-ST UP (TLR5 ligand, 30 ng/ml), R848 (TLR7/8 ligand, 10 μg/ml), ODN 1826 (TLR9 ligand, 10 μg/ml), Tri-DAP (NOD1 ligand, 10 μg/ml), and MDP (NOD2 ligand, 10 μg/ml). After 24h incubation, the NF-κB-induced SEAP activity was assessed 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 3622-34-80  
 E-mail: [info@invivogen.com](mailto:info@invivogen.com)