

Validation data for HEK-Dual™ hTLR4 cells

<https://www.invivogen.com/hek-dual-htlr4>

For research use only

Version 24C19-AK

HEK-Dual™ hTLR4 cells were generated from the HEK-Dual™ cell line through the stable expression of the human Toll-like receptor 4 (hTLR4), as well as the adapter proteins myeloid differentiation factor 2 (MD-2), and cluster of differentiation 14 (CD14). These cells feature two reporter genes allowing the simultaneous study of NF- κ B- and IRF-induced responses, by monitoring the SEAP (secreted embryonic alkaline phosphatase) and Lucia luciferase activities, respectively. These cells show strong NF- κ B and IRF responses upon incubation with TLR4 ligands, such as Lipopolysaccharide (LPS) and monophosphoryl Lipid A (MPLA), when compared to their parental cells HEK-Dual™ (Figures 1 & 2). Of note, as HEK293 cells express endogenous levels of TLR3 and TLR5, HEK-Dual™ - derived cells respond to their cognate ligands. Using CLI-095, a small-molecule inhibitor of TLR4 signaling, completely abrogates NF- κ B and IRF responses in HEK-Dual™ hTLR4 cells (Figure 3).

Functional validation of HEK-Dual™-derived cells (NF- κ B responses)

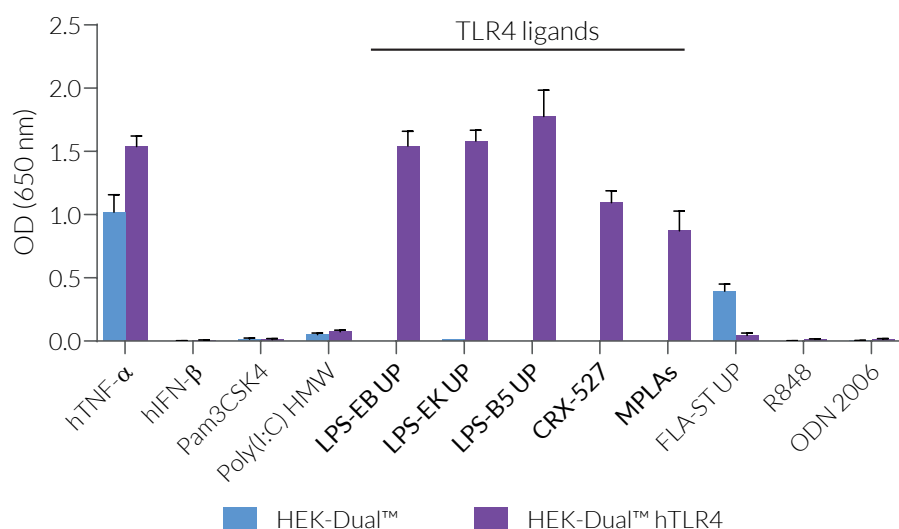


Figure 1. NF- κ B responses in HEK-Dual™-derived cells. HEK-Dual™ and HEK-Dual™ hTLR4 cells were incubated for 24 hours with cytokines and various TLR agonists: Human TNF- α (NF- κ B-positive control, 1 ng/ml), hIFN- β (IRF-positive control, 1000 U/ml), Pam3CSK4 (TLR2 ligand, 100 ng/ml), Poly(I:C) HMW (TLR3 ligand, 1 μ g/ml), LPS-EB Ultrapure (UP) (TLR4 ligand, 1 ng/ml), LPS-EK UP (TLR4 ligand, 1 ng/ml), LPS-B5 UP (TLR4 ligand, 100 pg/ml), CRX-527 (TLR4 ligand, 1 ng/ml), MPLAs (TLR4 ligand, 100 ng/ml), FLA-ST UP (TLR5 ligand, 100 ng/ml), R848 (TLR7/8 ligand, 10 μ g/ml), or ODN 2006 (TLR9 ligand, 10 μ g/ml). After 24h incubation, the NF- κ B-induced SEAP activity was assessed using QUANTI-Blue™. Data are shown as optical density (OD) at 650 nm (mean \pm SEM).

Functional validation of HEK-Dual™-derived cells (IRF responses)

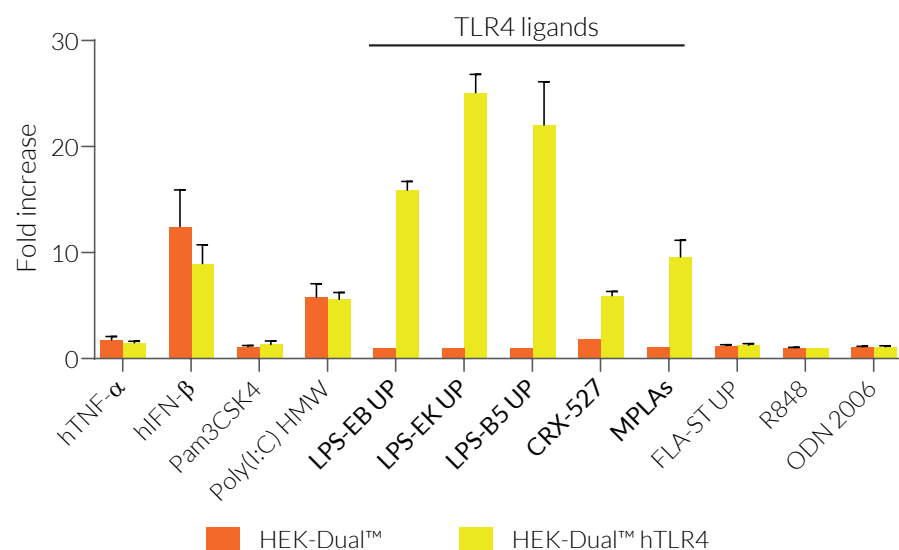


Figure 2. IRF responses in HEK-Dual™-derived cells. HEK-Dual™ and HEK-Dual™ hTLR4 cells were incubated for 24 hours with cytokines and various TLR agonists: Human TNF- α (NF- κ B-positive control, 10 ng/ml), hIFN- β (IRF-positive control, 10 U/ml), Pam3CSK4 (TLR2 ligand, 100 ng/ml), Poly(I:C) HMW (TLR3 ligand, 100 ng/ml), LPS-EB Ultrapure (UP) (TLR4 ligand, 10 ng/ml), LPS-EK UP (TLR4 ligand, 10 ng/ml), LPS-B5 UP (TLR4 ligand, 1 ng/ml), CRX-527 (TLR4 ligand, 10 ng/ml), MPLAs (TLR4 ligand, 100 ng/ml), FLA-ST UP (TLR5 ligand, 100 ng/ml), R848 (TLR7/8 ligand, 10 μ g/ml), or ODN 2006 (TLR9 ligand, 10 μ g/ml). After 24h incubation, the IRF response was assessed by measuring the activity of Lucia luciferase in the supernatant using QUANTI-Luc™. Data are shown in fold response over non-induced cells (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

Inhibition of TLR4-dependent NF- κ B and IRF responses

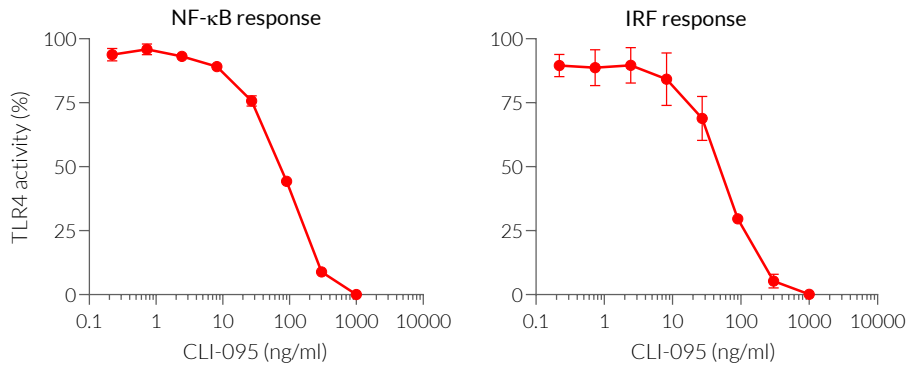


Figure 3. Inhibition of hTLR4 signaling pathway using CLI-095. HEK-Dual™ hTLR4 cells were incubated in the presence of increasing concentrations of CLI-095 for 3 hours before adding 10 ng/ml LPS-EK Ultrapure (TLR4 agonist). After overnight incubation, the inhibitory effect of CLI-095 on the NF- κ B and IRF pathways was determined by measuring the reduction of SEAP and Lucia production in the supernatant using the QUANTI-Blue™ and QUANTI-Lucia™ detection reagents, respectively. Data are shown as a percentage (%) of maximal TLR4 activation (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