

Validation data for THP1-Dual™ MD2-CD14 KO-TLR4 Cells

<https://www.invivogen.com/thp1-dual-mc-ko-tnfr4>

For research use only

Version 22L01-AK

THP1-Dual™ MD2-CD14 KO-TLR4 cells were generated from THP1-Dual™ cells as a control cell line for the THP1-Dual™ MD2-CD14-TLR4 cells, through a biallelic knockout of the human (h)TLR4 gene (Figure S1). Additionally, they express two co-adaptors important for lipopolysaccharide (LPS) signaling, MD-2 and CD14. They also feature two reporter genes, allowing the simultaneous study of NF-κB- and IRF-induced responses by monitoring the SEAP (secreted embryonic alkaline phosphatase) and Lucia luciferase activities, respectively. Upon stimulation with smooth and rough LPS, we observe the complete loss of NF-κB-mediated responses when compared to the parental THP1-Dual™ cells (Figure 1). There is no notable difference in the IRF-mediated responses between THP1-Dual™ MD2-CD14 KO-TLR4 cells and their parental cell line (Figure 2). However, there are strong NFκB and IRF responses to both smooth and rough LPS in the cell line overexpressing TLR4, THP1-Dual™ MD2-CD14-TLR4 (Figures 1&2).

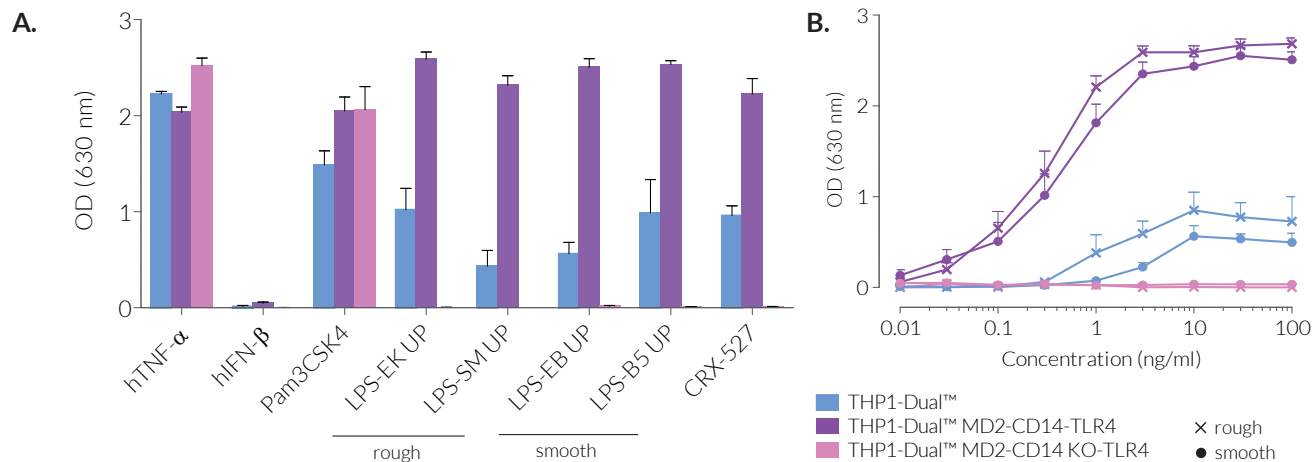


Figure 1: NF-κB responses in THP1-Dual™-derived cells. **A.** The cells were incubated with a variety of rough and smooth LPS: LPS-EK Ultrapure (UP) (rough; 10 ng/ml), LPS-SM UP (rough; 10 ng/ml), LPS-EB UP (smooth; 10 ng/ml), LPS-B5 UP (smooth; 1 ng/ml) and CRX-527 (synthetic; 10 ng/ml). Human (h)TNF-α (10 ng/ml) and hIFN-β (1x10³ U/ml) were used as a NF-κB-SEAP positive control and as a IRF positive control, respectively. The TLR2-TLR1 agonist Pam3CSK4 (10 ng/ml) served as a functional control. **B.** The cells were incubated with increasing concentration of rough (LPS-EK UP) and smooth (LPS-EB UP) LPS. After 24 hours, the activation of NF-κB was assessed by measuring the activity of SEAP in the supernatant using QUANTI-Blue™ Solution. Data are shown as optical density (OD) at 630 nm (mean ± SEM).

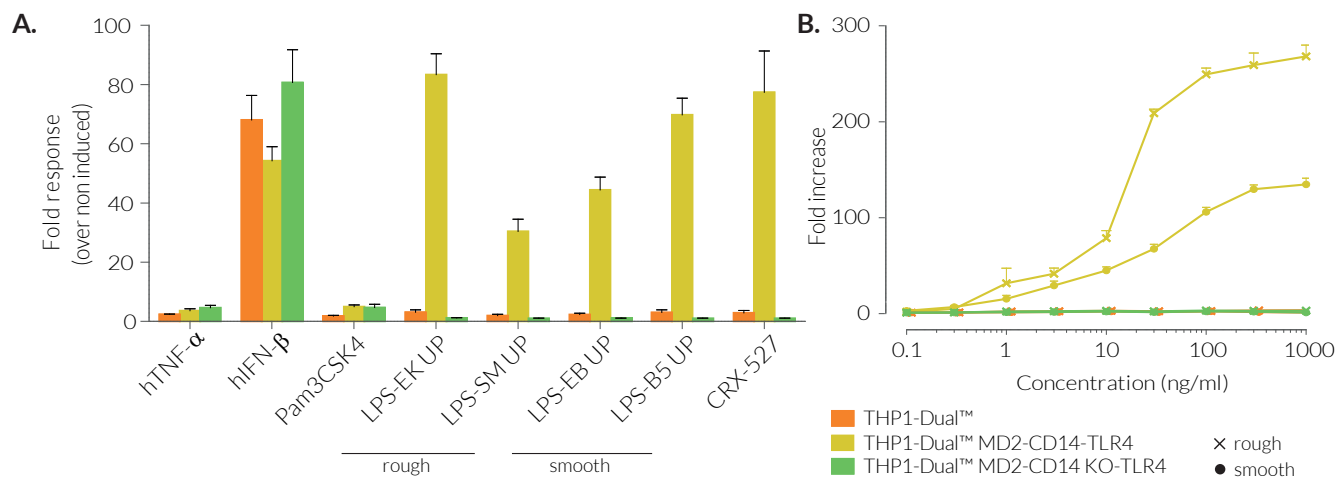


Figure 2: IRF responses in THP1-Dual™-derived cells. The cells were treated as described above (see figure 1). After 24 hours, the IRF response was assessed by measuring the activity of Lucia luciferase in the supernatant using QUANTI-Luc™. Data are shown in fold response over non-induced cells (mean ± 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

Validation of TLR4 knockout

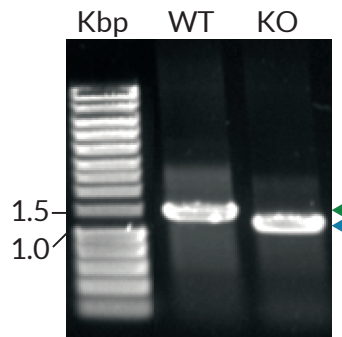


Figure S1: Validation of TLR4 KO. The targeted TLR4 region in parental THP1-Dual™ cells (WT; green arrow) and THP1-Dual™ MD2-CD14 KO-TLR4 (KO; blue arrow) cells was amplified by PCR. THP1-Dual™ MD2-CD14 KO-TLR4 cells were generated by a biallelic deletion causing the inactivation of hTLR4.

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