

Validation data for THP1-Dual™ KO-TRIF Cells

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

For research use only

Version 22L12-AK

THP1-Dual™ KO-TRIF cells were generated from the THP1-Dual™ cell line through the verified stable knockout of the human (h)TRIF gene (Figure 1). These cells 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. Lucia luciferase and SEAP activities are readily assessable in the cell culture supernatant using QUANTI-Luc™ and QUANTI-Blue™ Solution detection reagents, respectively. As expected, the prior differentiation treatment using PMA (Phorbol 12-myristate 13-acetate) does not rescue the IRF-response in LPS-stimulated TRIF-KO cells in comparison to their parental cell line THP1-Dual™ (Figure 2). The NF-κB response is slightly diminished in PMA-differentiated THP1-Dual™ KO-TRIF cells upon incubation with TLR4-specific ligands such as LPS and CRX-527. Of note, the remaining NF-κB-dependent signal to TLR4 ligand stimulation is most likely due to the more dominant activation of the TLR4-MyD88 signaling pathway (Figure 3).

Validation of TRIF knockout

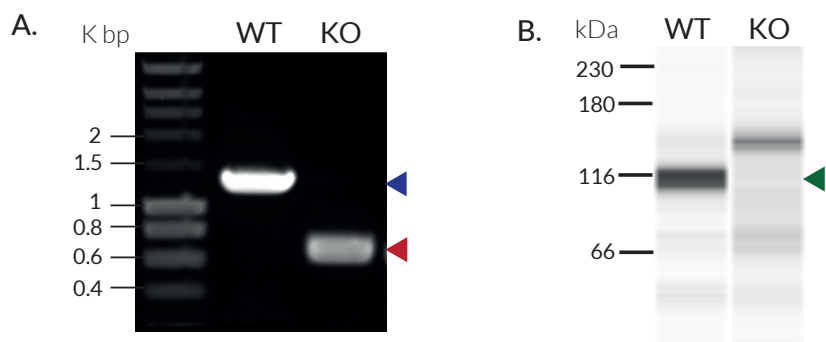


Figure 1: Validation of TRIF KO. (A) The targeted TRIF region in THP1-Dual™ (WT; blue arrow) parental cells and THP1-Dual™ KO-TRIF (KO; red arrow) cells was amplified by PCR. THP1-Dual™ KO-TRIF cells feature a frameshift deletion, causing an early stop codon and inactivation of TRIF. (B) Lysates from THP1-Dual™ (WT) and THP1-Dual™ KO-TRIF (KO) cells were analyzed using an anti-human TRIF antibody (green arrow), followed by an HRP-conjugated anti-rabbit secondary antibody. As expected a band was detected at ~100 kDa in the WT cells only.

Functional validation of TRIF knockout upon PMA differentiation (IRF response)

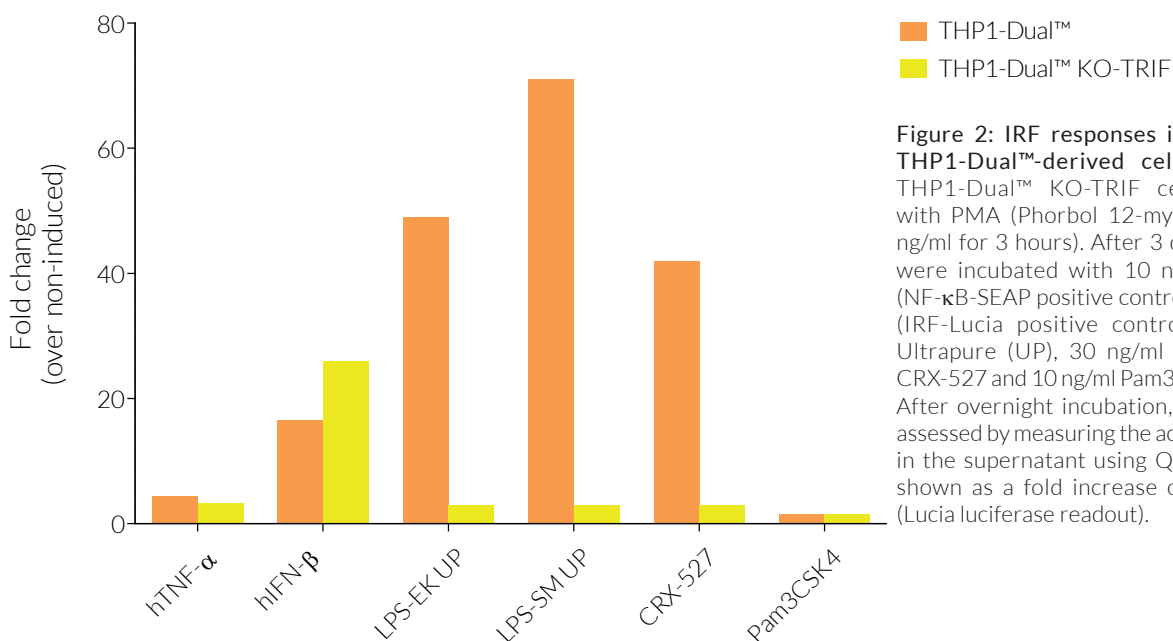


Figure 2: IRF responses in PMA-differentiated THP1-Dual™-derived cells. THP1-Dual™ and THP1-Dual™ KO-TRIF cells were pretreated with PMA (Phorbol 12-myristate 13-acetate; 10 ng/ml for 3 hours). After 3 days recovery, the cells were incubated with 10 ng/ml human (h)TNF-α (NF-κB-SEAP positive control), 1 x 10⁴ U/ml hIFN-β (IRF-Lucia positive control), 30 ng/ml LPS-EK Ultrapure (UP), 30 ng/ml LPS-SM-UP, 30 ng/ml CRX-527 and 10 ng/ml Pam3CSK4 (TLR2/1 agonist). After overnight incubation, the IRF response was assessed by measuring the activity of Lucia luciferase in the supernatant using QUANTI-Luc™. Data are shown as a fold increase over non-induced cells (Lucia luciferase readout).

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

Functional validation of TRIF knockout upon PMA differentiation (NF- κ B response)

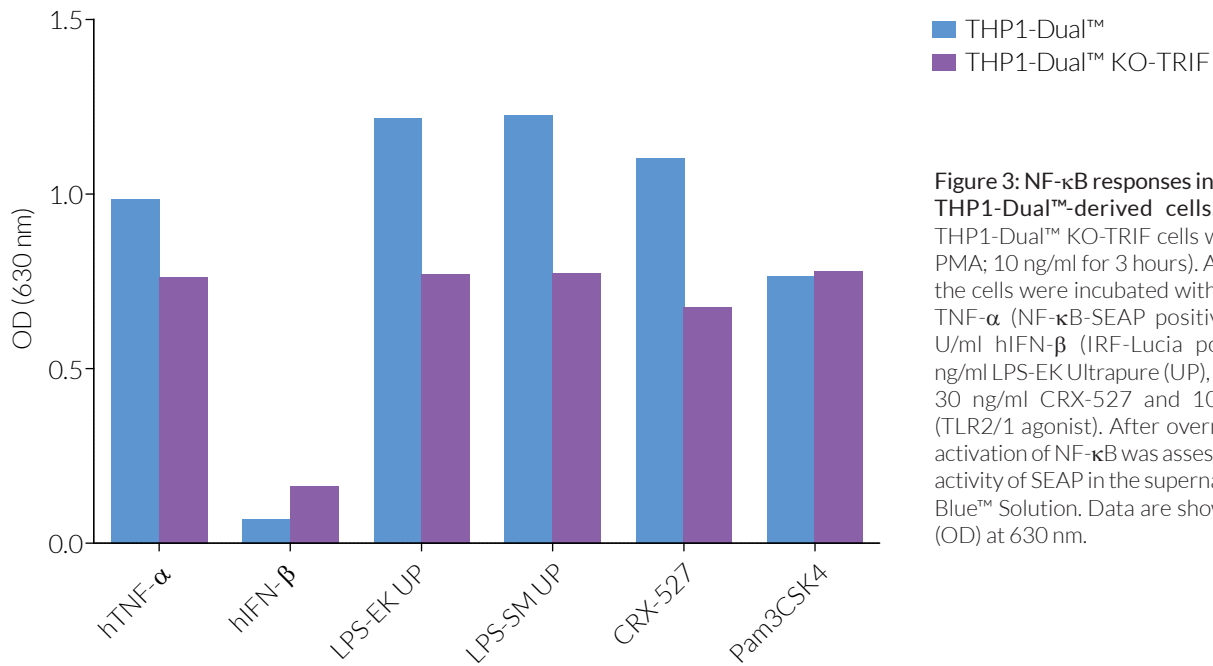


Figure 3: NF- κ B responses in PMA-differentiated THP1-Dual™-derived cells. THP1-Dual™ and THP1-Dual™ KO-TRIF cells were pretreated with PMA; 10 ng/ml for 3 hours). After 3 days recovery, the cells were incubated with 10 ng/ml human (h) TNF- α (NF- κ B-SEAP positive control), 1×10^4 U/ml hIFN- β (IRF-Lucia positive control), 30 ng/ml LPS-EK Ultrapure (UP), 30 ng/ml LPS-SM UP, 30 ng/ml CRX-527 and 10 ng/ml Pam3CSK4 (TLR2/1 agonist). After overnight incubation, the activation of NF- κ B 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.

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