

Validation data for THP1-Dual™ KO-IRF7 cells

<https://www.invivogen.com/thp1-dual-ko-irf7>

For research use only

Version 21L02-NJ

THP1-Dual™ KO-IRF7 cells were generated from the THP1-Dual™ cell line through the stable knockout (KO) of the *IRF7* gene, as verified by Western blot (Figure 1). These cells 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. We observe an impact of the IRF-7 knock-out on NF-κB-mediated responses (Figure 2), and no effect on IRF-mediated responses (Figure 3), in THP1-Dual™ KO-IRF7 cells upon incubation with TLR agonists (LPS-EK, R848) when compared to the THP1-Dual™ cells.

Validation of *IRF7* knockout

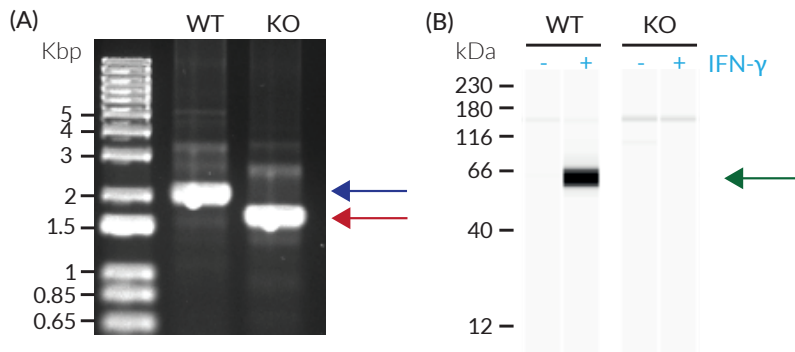


Figure 1: Validation of IRF7 KO. (A) The targeted *IRF7* region in THP1-Dual™ (WT; blue arrow) parental cells and THP1-Dual™ KO-IRF7 (KO; red arrow) cells was amplified by PCR. THP1-Dual™ KO-IRF7 cells feature a frameshift deletion, causing an early stop codon and inactivation of IRF7. (B) THP1-Dual™ (WT) and THP1-Dual™ KO-IRF7 (KO) cells were treated with 20 ng/ml interferon (IFN)- γ to induce the expression of IRF7. Lysates from non-treated (-) and treated (+) WT and KO cells were analyzed using an anti-human IRF7 antibody, followed by an HRP-conjugated anti-rabbit secondary antibody (WEST™ assay). As expected, a band was detected in the IFN- γ treated WT cells only (green arrow). The predicted molecular weight of IRF7 is ~66 kDa.

Functional validation of *IRF7* knockout (NF-κB response)

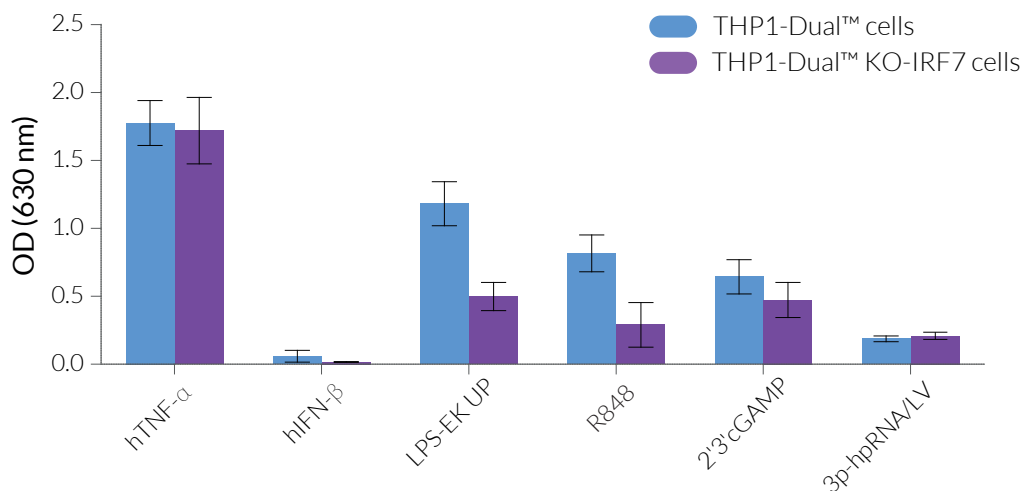


Figure 2: NF-κB responses in THP1-Dual™-derived cells. THP1-Dual™ and THP1-Dual™ KO-IRF7 cells were incubated with 1 ng/ml human (h)TNF- α (NF-κB-SEAP positive control), 1000 U/ml hIFN- β (IRF-Lucia positive control), 10 ng/ml LPS-EK Ultrapure (UP; TLR4 agonist), 3 μ g/ml R848 (TLR7/8 agonist), 30 μ g/ml 2'3'-cGAMP (STING agonist), and 1 μ g/ml 3p-hpRNA complexed with LyoVec™ (LV; RIG-I agonist). After overnight incubation, 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 \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

Functional validation of *IRF7* knockout (IRF response)

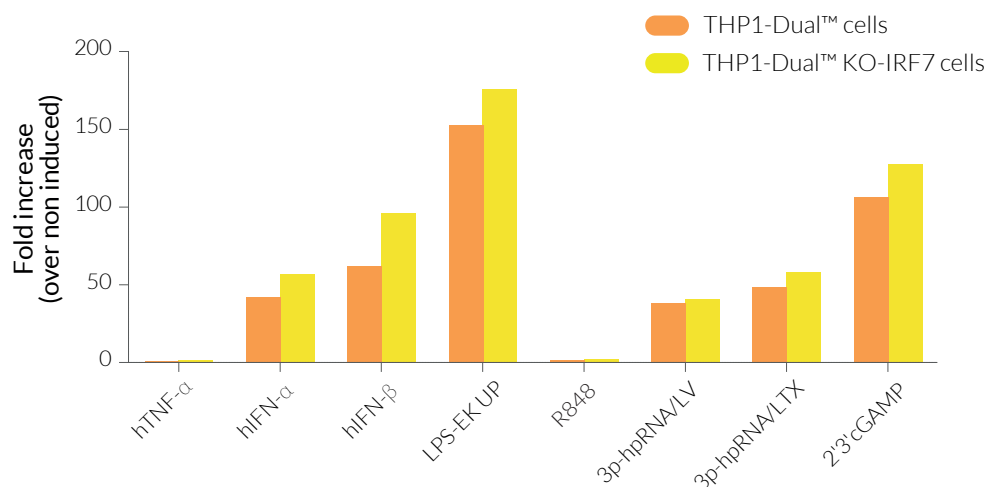


Figure 3: IRF responses in PMA-differentiated THP1-Dual™-derived cells. THP1-Dual™ and THP1-Dual™ KO-IRF7 cells were pretreated with PMA (Phorbol 12-myristate 13-acetate; 10 ng/ml for 3 hours). After 3 days recovery, the cells were incubated with 1 ng/ml human (h) TNF- α (NF- κ B-SEAP positive control), 1000 U/ml hIFN- α or hIFN- β (IRF-Lucia positive control), 10 ng/ml LPS-EK Ultrapure (UP; TLR4 agonist), 1 μ g/ml R848 (TLR7/8 agonist), 100 ng/ml 3p-hpRNA complexed with either LyoVec™ (LV) or LTX (RIG-I agonist), and 10 μ g/ml 2'3'-cGAMP (STING 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