

# Validation data for THP1-Dual™ KO-*IKKε* cells

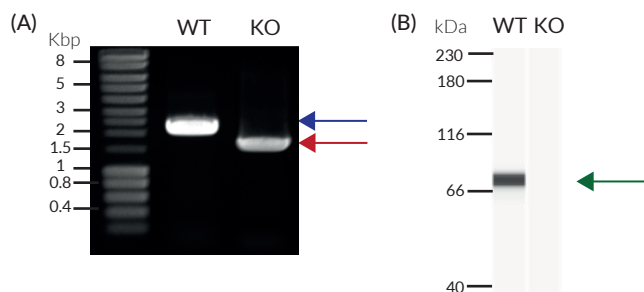
<https://www.invivogen.com/thp1-dual-ko-ikke>

For research use only

Version 21B04-ED

THP1-Dual™ KO-*IKKε* cells were generated from the THP1-Dual™ cell line through the stable knockout (KO) of the *IKKε* gene, as verified by PCR and Western blot (Figure 1). These cells feature two reporter genes allowing the simultaneous study of the IRF pathway by monitoring the activity of an inducible secreted Lucia luciferase, and the NF- $\kappa$ B pathway by monitoring the activity of an inducible SEAP (secreted embryonic alkaline phosphatase). Importantly, comparison to our THP1-Dual™ KO-TBK1 cells, which feature a KO of the highly similar IKK-related kinase TBK1, allows overlap and differences in signaling functions between *IKKε* and TBK1 to be fully elucidated (Figure 2-3).

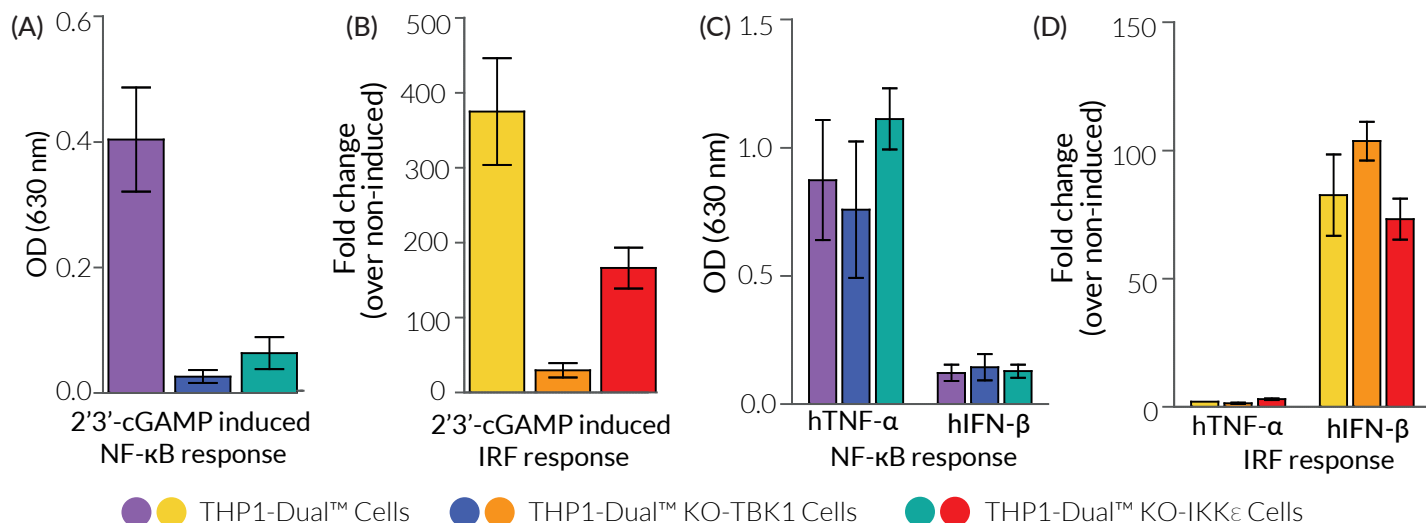
## Validation of *IKKε* knockout



**Figure 1: Validation of *IKKε* KO.** (A) The targeted *IKKε* region in THP1-Dual™ (WT; blue arrow) parental cells and THP1-Dual™ KO-*IKKε* (KO; red arrow) cells was amplified by PCR. THP1-Dual™ KO-*IKKε* cells feature a frameshift deletion, causing an early stop codon and inactivation of *IKKε*. (B) Lysates from THP1-Dual™ (WT) and THP1-Dual™ KO-*IKKε* (KO) cells were analyzed using an anti-human *IKKε* antibody (green arrow), followed by a HRP-conjugated anti-rabbit secondary antibody (WES assay). As expected a band was detected at ~80 Da in the WT cells only.

## Functional validation of *IKKε* knockout (NF- $\kappa$ B and IRF response)

*IKKε* and TBK1 have both been shown to coordinate the activation of IRF3 and NF- $\kappa$ B in the innate immune response. Thus, as expected the NF- $\kappa$ B-dependent response in THP1-Dual™ KO-*IKKε* upon stimulation with the STING agonist, 2'3'-cGAMP, is greatly reduced and comparable to THP1-Dual™ KO-TBK1 (A). However, *IKKε* seems to play a lesser role in IRF-dependent STING signaling (B). Notably, THP1-Dual™ KO-*IKKε* cells retain a full ability to respond to (C) NF- $\kappa$ B- and (D) IRF-dependent cytokines such as human hTNF- $\alpha$  and type I interferons (hIFN- $\beta$ ), respectively, for which signaling is *IKKε*-independent.



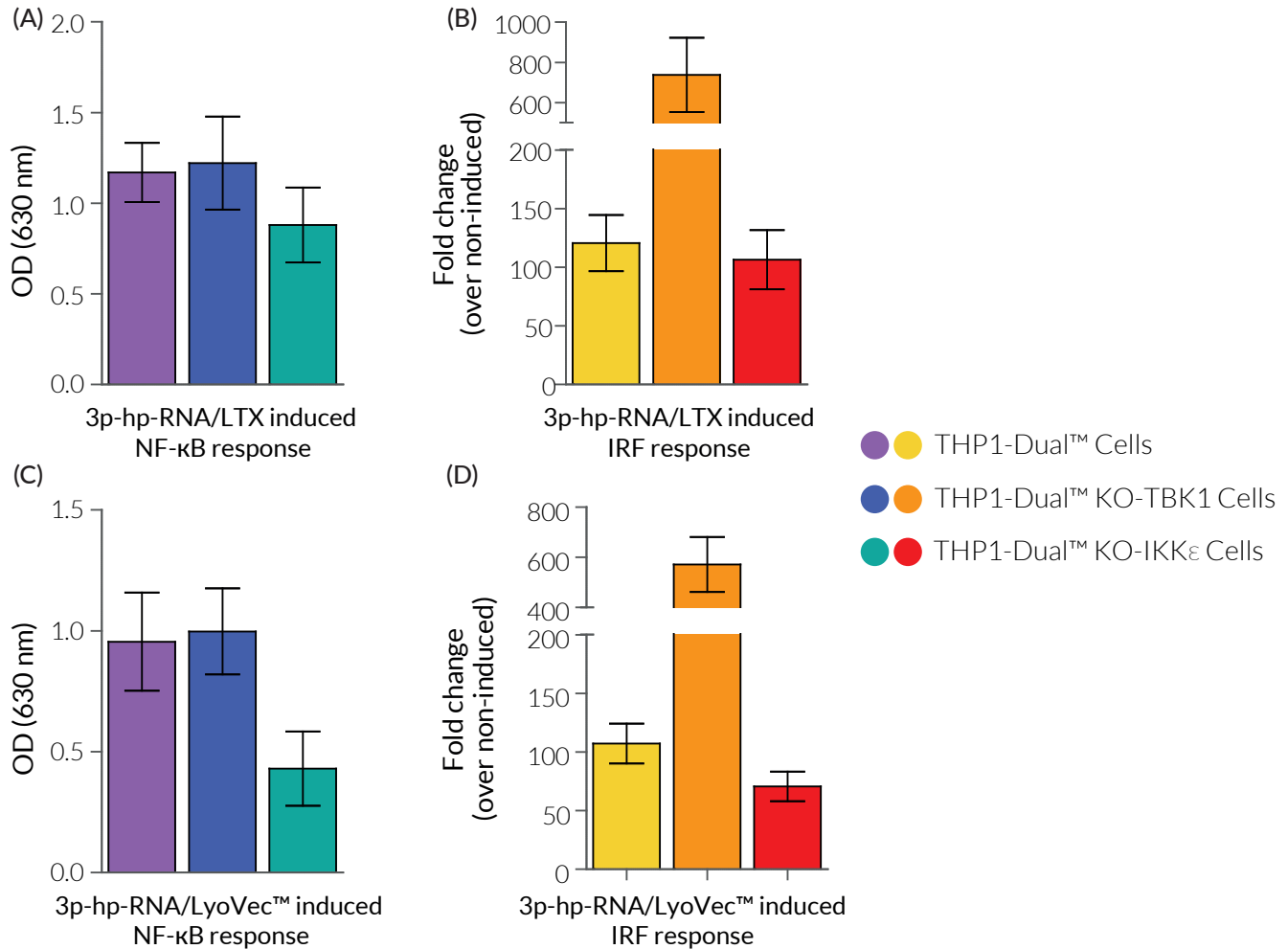
**Figure 2: NF- $\kappa$ B and IRF responses in THP1-Dual™-derived cells.** THP1-Dual™ (WT) and THP1-Dual™ KO-TBK1, and THP1-Dual™ KO-*IKKε* cells were incubated with (A-B) 30  $\mu$ g/ml 2'3'-cGAMP (STING agonist), (C-D) 0.3 ng/ml human TNF- $\alpha$  (hTNF- $\alpha$ ), or 10<sup>4</sup> U/ml human IFN- $\beta$  (hIFN- $\beta$ ). After overnight incubation, the NF- $\kappa$ B activity was assessed by measuring the SEAP activity in the supernatant using QUANTI-Blue™ Solution. Data are shown as optical density (OD) at 630 nm (mean  $\pm$  SEM) (A, C). The IRF response was assessed by measuring Lucia luciferase activity in the supernatant using QUANTI-Luc™. Data are shown as a fold change (mean  $\pm$  SEM) over non-induced cells (B, D).

### 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 Hong Kong: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)

## Evaluation of IRF and NF- $\kappa$ B responses upon RNA transfection

3p-hpRNA is an agonist of RIG-I that must be complexed with a transfection reagent to enter the cytoplasm. Upon stimulation with 3p-hp-RNA complexed to LTX, the NF- $\kappa$ B-dependent response in THP1-Dual™ KO- $\text{IKK}\epsilon$  is barely affected, similar to THP1-Dual™ KO-TBK1 (A). However, when complexed to LyoVec™, a reduction in activity is noted in THP1-Dual™ KO- $\text{IKK}\epsilon$  cells only (C). Interestingly, with either transfection reagent (B: LTX or D: LyoVec™), the IRF-dependant response of THP1-Dual™ KO- $\text{IKK}\epsilon$  is comparable to the parental cell line THP1-Dual™. This is noticeably different to THP1-Dual™ KO-TBK1, where the IRF-response is unexpectedly increased. This opposing observation between  $\text{IKK}\epsilon$  and TBK1 in the IRF-dependent response to RNA ligands may suggest that in the absence of TBK1,  $\text{IKK}\epsilon$  homodimers play a greater role in the RNA-sensing signaling pathway.



**Figure 3: RNA-induced IRF and NF- $\kappa$ B responses in THP1-Dual™-derived cells.** THP1-Dual™, THP1-Dual™ KO-TBK1, or THP1-Dual™ KO- $\text{IKK}\epsilon$  cells were transfected with 1  $\mu\text{g}/\text{ml}$  3p-hpRNA complexed with LTX (A, B) or LyoVec™ (C, D). After overnight incubation, the NF- $\kappa$ B activity was assessed by measuring the SEAP activity in the supernatant using QUANTI-Blue™ Solution. Data are shown as optical density (OD) at 630 nm (mean  $\pm$  SEM) (A, C). The IRF response was assessed by measuring Lucia luciferase activity in the supernatant using QUANTI-Luc™. Data are shown as a fold change (mean  $\pm$  SEM) over non-transfected cells (B, D).

### 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 Hong Kong: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)