

# Validation data for THP1-Dual™ KO-MDA5 cells

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

For research use only

Version 21C29-NJ

THP1-Dual™ KO-MDA5 cells were generated from the THP1-Dual™ cell line through the stable knockout (KO) of the *MDA5* 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-κB pathway by monitoring the activity of an inducible SEAP (secreted embryonic alkaline phosphatase). As expected, the IRF and NF-κB responses are severely impaired in THP1-Dual™ KO-MDA5 cells upon incubation with Poly(I:C) HMW, a RLR inducer. Of note, the responses to 3p-hpRNA, a RIG-I-specific ligand, are also decreased (Figures 2-3). These cells retain the ability to respond to other NF-κB activating ligands such as recombinant human TNF-α, as well as IRF-activating ligands such as type I interferons (IFNs) (Figures 2-3).

## Validation of MDA5 knockout

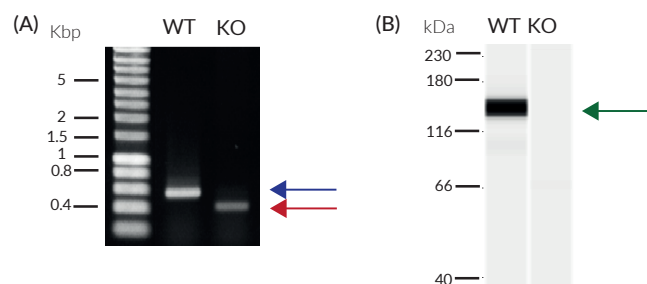


Figure 1: Validation of MDA5 KO.

(A) The targeted MDA5 region in THP1-Dual™ (WT; blue arrow) parental cells and THP1-Dual™ KO-MDA5 (KO; red arrow) cells was amplified by PCR. THP1-Dual™ KO-MDA5 cells feature a frameshift deletion, causing an early stop codon and inactivation of MDA5. (B) Lysates from THP1-Dual™ (WT) and THP1-Dual™ KO-MMDA5(KO) cells were analyzed using an anti-human MDA5 antibody (green arrow), followed by a HRP-conjugated anti-rabbit secondary antibody (WES assay). As expected, a band was detected at ~135 KDa in the WT cells only.

## Functional validation of MDA5 knockout (IRF response)

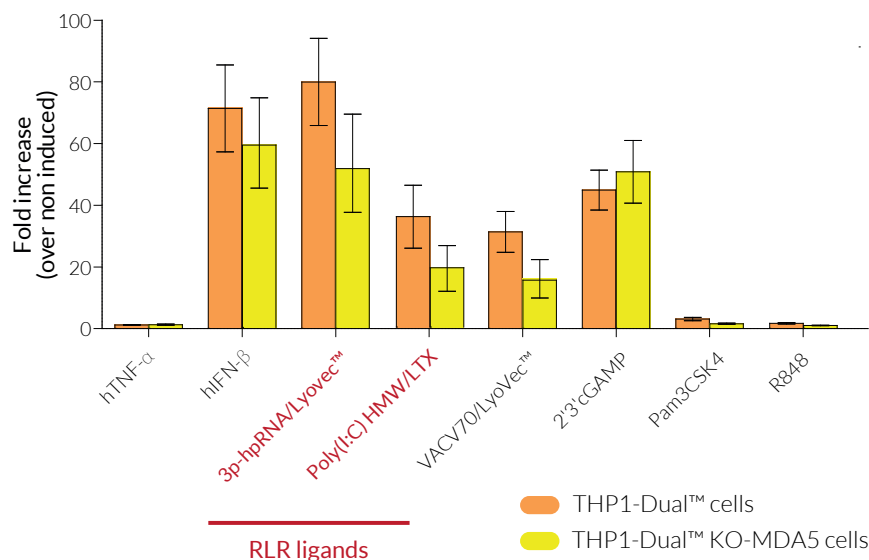


Figure 2: IRF responses in THP1-Dual™-derived cells.

THP1-Dual™ (WT) and THP1-Dual™ KO-MDA5 cells were incubated with 1 ng/ml human TNF-α (hTNF-α), or 10<sup>4</sup> U/ml human IFN-β (hIFN-β), 300 ng/ml 3p-hpRNA/Lyovec, 1 μg/ml Poly(I:C) HMW/LTX, 1 μg/ml VACV-70/Lyovec, 3 μg/ml 2'3'-cGAMP, 3 ng/ml Pam3CSK4, or 1 μg/ml R848. After overnight incubation, the IRF response was assessed by measuring Lucia luciferase activity in the supernatant using QUANTI-Luc™. Data are shown as a fold change (mean ± SEM) over non-induced cells.

### TECHNICAL SUPPORT

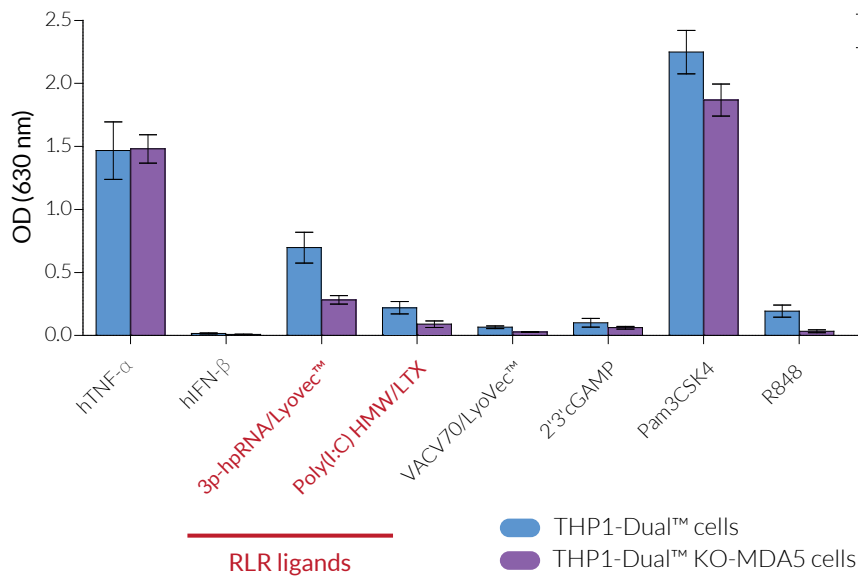
InvivoGen USA (Toll-Free): 888-457-5873  
InvivoGen USA (International): +1 (858) 457-5873  
InvivoGen Europe: +33 (0) 5-62-71-69-39  
InvivoGen Hong Kong: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)



Any questions about our cell lines?  
Visit our FAQ page.

**InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)

## Functional validation of MDA5 knockout (NF-κB response)



**Figure 3: NF-κB responses in THP1-Dual™-derived cells.**

THP1-Dual™ (WT) and THP1-Dual™ KO-MDA5 cells were incubated with 1 ng/ml human TNF-α (hTNF-α), or 10<sup>4</sup> U/ml human IFN-β (hIFN-β), 300 ng/ml 3p-hpRNA/Lyovec, 1 μg/ml Poly(I:C) HMW/LTX, 1 μg/ml VACV-70/Lyovec, 3 μg/ml 2'3'-cGAMP, 3 ng/ml Pam3CSK4, or 1 μg/ml R848. After overnight incubation, the NF-κB 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 ± 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 Hong Kong: +852 3622-3480  
 E-mail: [info@invivogen.com](mailto:info@invivogen.com)



Any questions about our cell lines?  
 Visit our FAQ page.