

# Validation data for THP1-Dual™ KO-SAMHD1 Cells

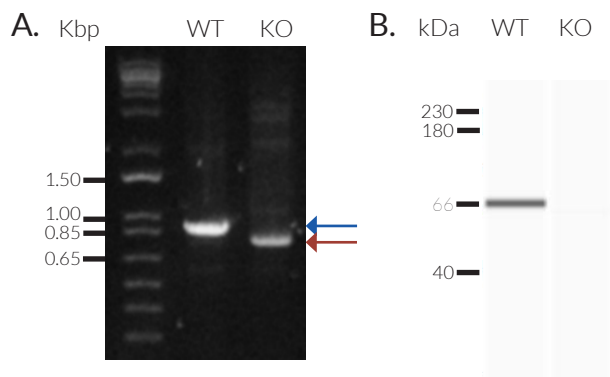
<https://www.invivogen.com/thpd-kosamhd1>

For research use only

Version 22F16-AK

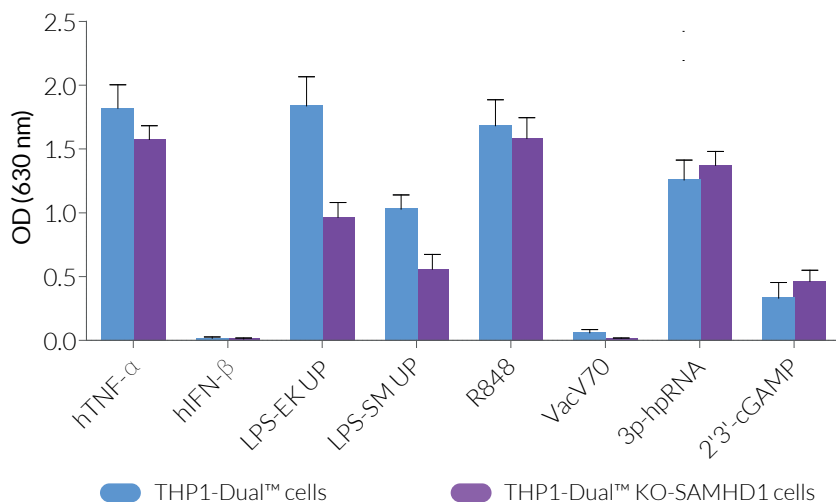
THP1-Dual™ KO-SAMHD1 cells were generated from the THP1-Dual™ cell line through the stable biallelic knockout of the SAMHD1 gene, as verified by PCR and western blot (Figure 1, A and B). Additionally, these cells feature two reporter genes allowing the simultaneous study of NF- $\kappa$ B- and IRF-induced responses by monitoring the SEAP (secreted embryonic alkaline phosphatase) and Lucia luciferase activities, respectively (Figures 2 and 3). We observe a reduction of NF- $\kappa$ B-mediated responses induced by LPS-EK, and to a lesser extent by LPS-SM, two TLR4 agonists (Figure 2). There is no notable difference in the IRF-mediated responses between THP1-Dual™ KO-SAMHD1 and their parental cell line, upon incubation with TLR4, TLR7/8, CDS, RIG-I, or STING agonists (Figure 3).

## Validation of SAMHD1 knockout



**Figure 1: Validation of SAMHD1 KO.** (A) The targeted SAMHD1 region in THP1-Dual™ (WT; blue arrow) parental cells and THP1-Dual™ KO-SAMHD1 (KO; red arrow) cells was amplified by PCR. THP1-Dual™ KO-SAMHD1 cells were generated by a biallelic deletion of 132 bp in exon 7, causing the inactivation of SAMHD1. The WT PCR product is 868 bp, whereas the truncated KO band measures only 736 bp. (B) Lysates from THP1-Dual™ (WT) and THP1-Dual™ KO-SAMHD1 (KO) cells were analyzed using an anti-human SAMHD1 antibody, followed by an HRP-conjugated anti-rabbit secondary antibody (JESS™ system). As expected a band was detected at ~70 kDa in the WT cells only (green arrow).

## Functional validation of SAMHD1 knockout (NF- $\kappa$ B response)

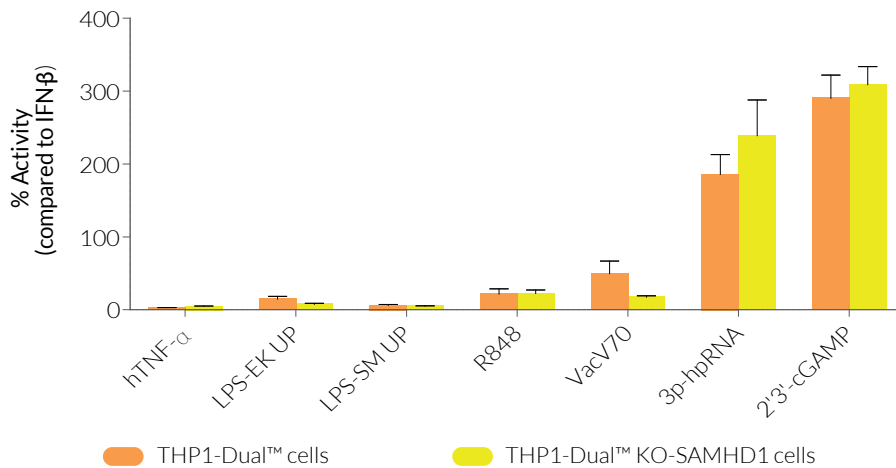


**Figure 2: NF- $\kappa$ B responses in THP1-Dual™-derived cells.** THP1-Dual™ and THP1 Dual™ KO-SAMHD1 cells were incubated with 1 ng/ml human (h)TNF- $\alpha$  (NF- $\kappa$ B-SEAP positive control), 1000 U/ml hIFN- $\beta$  (IRF-Lucia positive control), 100 ng/ml LPS-EK Ultrapure and 100  $\mu$ g/ml LPS-SM Ultrapure (both TLR4 agonists), 10  $\mu$ g/ml R848 (TLR7/8 agonist), 1  $\mu$ g/ml VacV70 complexed with LyoVec™ (CDS agonist), 1  $\mu$ g/ml 3p-hpRNA complexed with LyoVec™ (RIG-I agonist) and 30  $\mu$ g/ml, 2'3'-cGAMP (STING agonist). After overnight incubation, the activation of NF- $\kappa$ 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 (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](mailto:info@invivogen.com)

### Functional validation of SAMHD1 knockout (IRF response)



**Figure 3: IRF responses in THP1-Dual™-derived cells.** THP1-Dual™ and THP1-Dual™ KO-SAMHD1 cells were incubated with 1 ng/ml human (h)TNF- $\alpha$  (NF- $\kappa$ B-SEAP positive control), 100 ng/ml LPS-EK Ultrapure and 1  $\mu$ g/ml LPS-SM Ultrapure (both TLR4 agonists), 10  $\mu$ g/ml R848 (TLR7/8 agonist), 1  $\mu$ g/ml VacV70 complexed with LyoVec™ (CDS agonist), 1  $\mu$ g/ml 3p-hpRNA complexed with LyoVec™ (RIG-I agonist) and 30  $\mu$ 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™. The IRF induction of each ligand is expressed relative to that of hIFN- $\beta$  at  $1 \times 10^3$  U/ml (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](mailto:info@invivogen.com)