

Validation data for THP1-Dual™ KO-DNase2 Cells

<https://www.invivogen.com/thpd-kodnase2>

For research use only

Version 22F15-AK

THP1-Dual™ KO-DNase2 cells were generated from the THP1-Dual™ cell line through the stable biallelic knockout of the *DNASE2* gene, as verified by PCR (Figure 1). Additionally, these cells feature two reporter genes allowing the simultaneous study of NF- κ B- and IRF-induced responses by monitoring the SEAP (secreted embryonic alkaline phosphatase) and Lucia luciferase activities, respectively. There are no notable differences in the NF- κ B- and IRF-mediated responses in THP1-Dual™ KO-DNase2 cells when compared to their parental cell line (Figures 2 and 3).

Validation of DNASE2 knockout

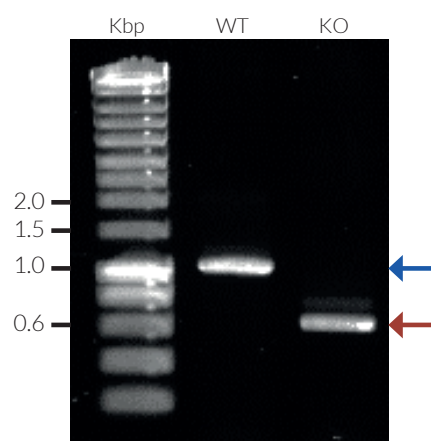


Figure 1: Validation of DNase2 KO. The targeted DNase2 region in THP1-Dual™ (WT; blue arrow) parental cells and THP1-Dual™ KO-DNase2 (KO; red arrow) cells was amplified by PCR. THP1-Dual™ KO-DNase2 cells were generated by a biallelic deletion of 443 bp in exon 2 and parts of exon 1 and 3, causing the inactivation of DNase2. The WT PCR product is 1133 bp, whereas the truncated KO band measures only 690 bp.

Functional validation of DNASE2 knockout (NF- κ B response)

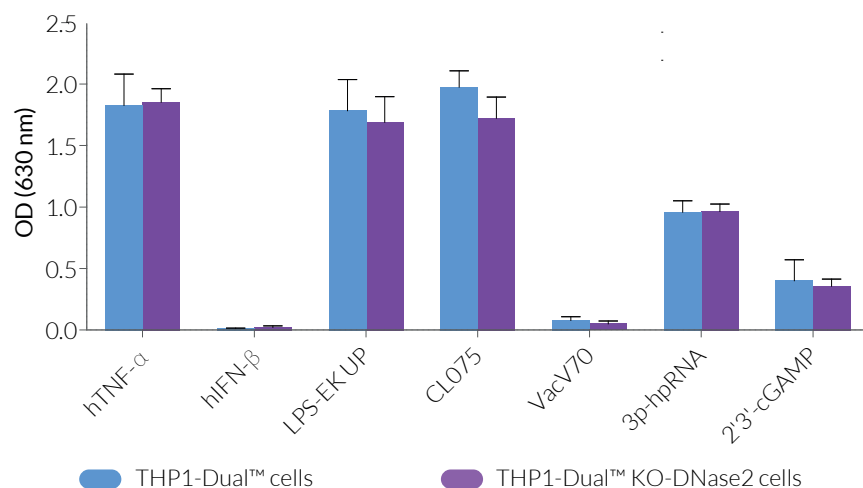


Figure 2: NF- κ B responses in THP1-Dual™-derived cells. THP1-Dual™ and THP1-Dual™ KO-DNase2 cells were incubated with 1 ng/ml human (h)TNF- α (NF- κ B-SEAP positive control), 1000 U/ml hIFN- β (IRF-Lucia positive control), 10 ng/ml LPS-EK Ultrapure (UP; TLR4 agonist), 10 μ g/ml CLO75 (TLR7/8 agonist), 1 μ g/ml VacV70 complexed with LyoVec™ (CDS agonist), 1 μ g/ml 3p-hpRNA complexed with LyoVec™ (RIG-I agonist) and 30 μ g/ml 2'3'-cGAMP (STING 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 (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 DNASE2 knockout (IRF response)

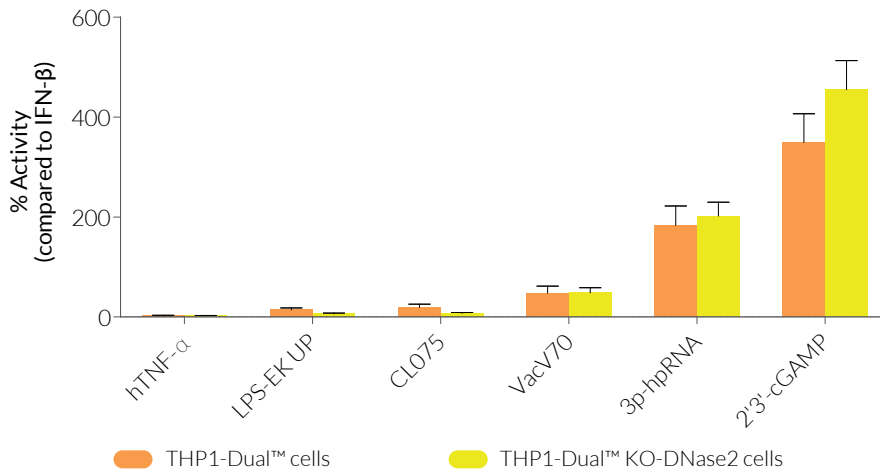


Figure 3: IRF responses in THP1-Dual™-derived cells. THP1-Dual™ and THP1-Dual™ KO-DNase2 cells were incubated with 1 ng/ml human (h)TNF- α (NF- κ B-SEAP positive control), 100 ng/ml LPS-EK Ultrapure (TLR4 agonist), 10 μ g/ml CLO75, 10 μ g/ml R848 (both TLR7/8 agonists), 1 μ g/ml VacV70 complexed with LyoVec™ (CDS agonist), 1 μ g/ml 3p-hpRNA complexed with LyoVec™ (RIG-I agonist) and 30 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™. The IRF induction of each ligand is expressed relative to that of hIFN- β at 1×10^3 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