

Validation data for THP1-KO-NLRP3 cells

<https://www.invivogen.com/thp1-ko-kd-nlrp3>

For research use only

Version 20F22-NJ

THP1-KO-NLRP3 cells were generated from the human monocytic THP1-Null2 cell line through the deletion of the N-terminal region of the *NLRP3* gene. Only an inactive *NLRP3* C-terminal fragment is expressed. The KO status has been verified by DNA sequencing, PCR (Figure 1A), Western blot (Figure 1B, C), and functional assays (Figure 2). Mature IL-1 β secretion is abolished in THP1-KO-NLRP3 cells (KO) after activation with inducers of the *NLRP3* inflammasome (i.e. Nigericin, MSU crystals, or Alum Hydroxide) when compared to the parental (WT) cell line (Figure 2A). The *NLRP3*-caspase-1 dependent maturation of IL-1 β is abolished upon indirect activation of *NLRP3* by inducers of the AIM2 inflammasome (i.e. Poly (dA:dT)) and non-canonical inflammasome (i.e. *E. coli* outer membrane vesicles (OMVs)) (Figure 2A). Furthermore, pyroptosis is abrogated in THP1-KO-NLRP3 cells upon *NLRP3* activation (Figure 2B).

Validation of *NLRP3* Knockout (KO)

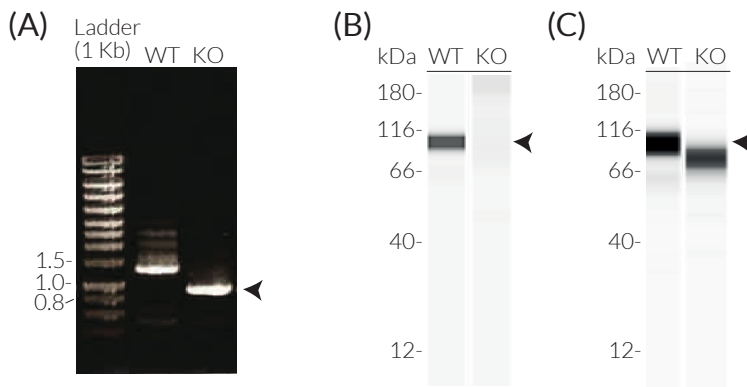


Figure 1: Validation of *NLRP3* KO in THP1-KO-NLRP3 cells. (A) The targeted *NLRP3* region in THP1-Null2 (WT) and THP1-KO-NLRP3 (KO) cells was amplified by PCR. THP1-KO-NLRP3 cells feature a biallelic deletion (arrow). (B and C) Lysates from THP1-Null2 (WT) and THP1-KO-NLRP3 (KO) cells were analyzed by Western blot (Wes™) using an anti-human *NLRP3* antibody, targeting the N-terminal region of *NLRP3* (B), or the C-terminal region of *NLRP3* (C), and followed by a HRP-conjugated anti-mouse or anti-rabbit secondary antibody. The arrow indicates the expected band for the *NLRP3* protein (110 kDa).

Functional validation of THP1-KO-NLRP3 cells

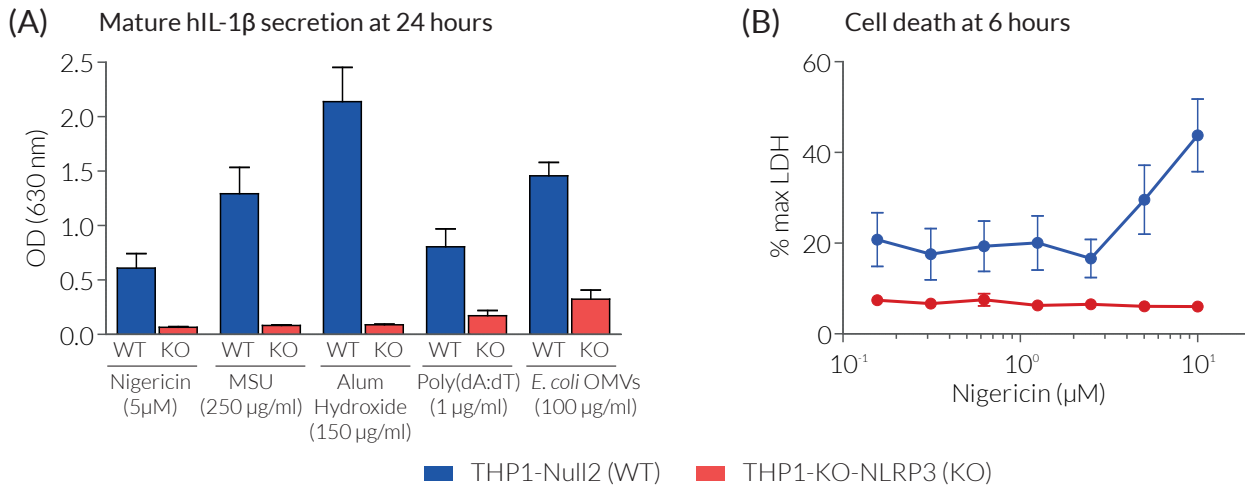


Figure 2: Absence of mature IL-1 β secretion and pyroptosis by THP1-KO-NLRP3 cells upon inflammasome activation.

$\sim 3 \times 10^5$ THP1-Null2 (blue) and THP1-KO-NLRP3 cells (red) were incubated for 3h at 37°C with LPS-EK (1 μ g/ml) (priming) and then incubated (activation) with Nigericin (A: 5 μ M; B: 0.15-10 μ M), MSU crystals (MSU; 250 μ g/ml), Alum Hydroxide (150 μ g/ml), transfected Poly(dA:dT) (1 μ g/ml), or *E. coli* outer membrane vesicles (OMVs) (100 μ g/ml). After 24h, (A) the secretion of mature human (h)IL-1 β was assessed in the culture supernatant using HEK-Blue™ IL-1 β sensor cells which express an NF- κ B SEAP reporter gene. QUANTI-Blue™ Solution was used to measure SEAP activity. Optical density (OD) was read at 630 nm. (B) After a 6 hour incubation with Nigericin, cell death was assessed using the lactate dehydrogenase (LDH) assay.

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 3-622-34-80
E-mail: info@invivogen.com

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

 **InvivoGen**
www.invivogen.com