

Validation data for THP1-KO-GSDMD cells

<https://www.invivogen.com/thp1-kogsdmd>

For research use only

Version 20G27-NJ

THP1-KO-GSDMD cells were generated from the human monocytic THP1-Null2 cell line through the stable knockout of the gasdermin D (*GSDMD*) gene. The KO status has been verified by PCR (Figure 1A), Western blot (Figure 1B), and functional assays (Figure 2). THP1-KO-GSDMD cells exhibit an impaired IL-1 β secretion upon canonical and non-canonical inflammasome activation with Nigericin and *E. coli* outer membrane vesicles (OMVs), respectively (Figure 2A). Although these cells are fully KO for the GSDMD protein, mature IL-1 β secretion is detected with the AIM2 inducer Poly(dA:dT) (Figure 2A), and with NLRP3 inducers such as MSU crystals or Alum Hydroxide (at 24 hours, data not shown). This observation illustrates the reported IL-1 β secretion upon GSDMD-independent cell death. Importantly, pyroptosis is abrogated in THP1-KO-GSDMD cells upon NLRP3 activation with Nigericin (Figure 2B).

Validation of GSDMD Knockout (KO)

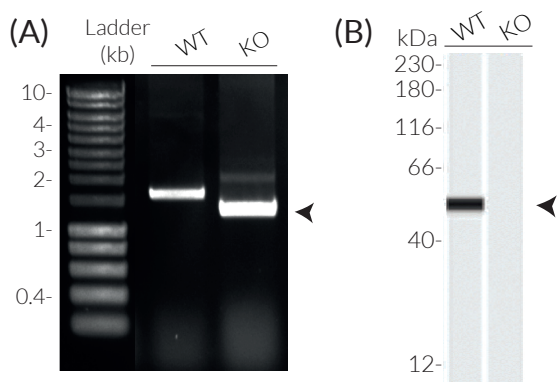


Figure 1: Validation of GSDMD KO in THP1-KO-GSDMD cells

(A) The targeted GSDMD region in THP1-Null2 (WT) and THP1-KO-GSDMD (KO) cells was amplified by PCR. THP1-KO-GSDMD cells feature a biallelic deletion (arrow). (B) Lysates from THP1-Null2 (WT) and THP1-KO-GSDMD (KO) cells were analyzed by Western blot (Wes™) using an anti-human GSDMD antibody and a HRP-conjugated anti-mouse secondary antibody. The arrow indicates the expected band for the GSDMD protein (48 kDa).

Functional validation of THP1-KO-GSDMD cells

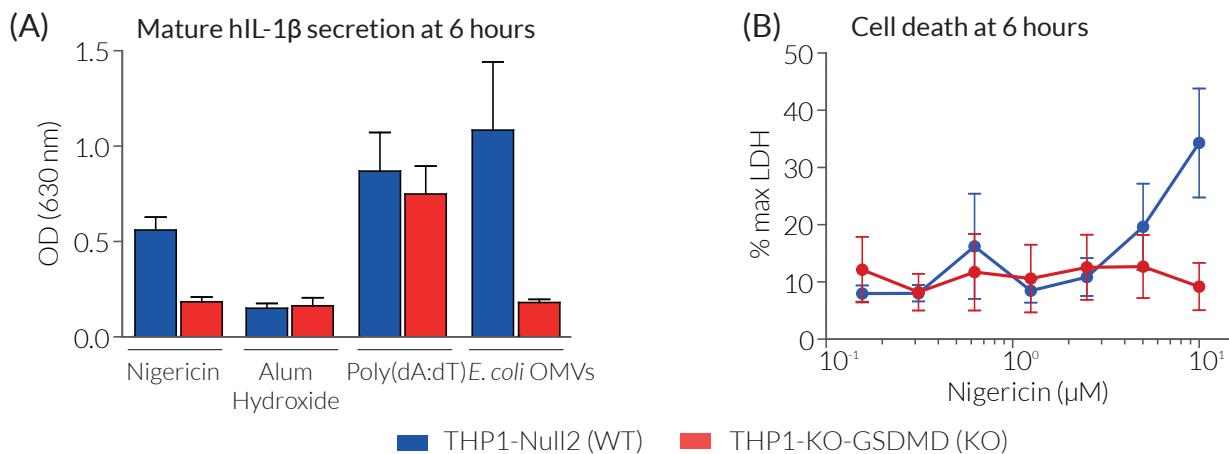


Figure 2: Altered mature IL-1 β secretion and pyroptosis by THP1-KO-GSDMD cells upon inflammasome activation.

~3x10⁵ THP1-Null2 (blue) and THP1-KO-GSDMD 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), 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 6h, (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