

# Validation data for THP1-KO-CASP4 cells

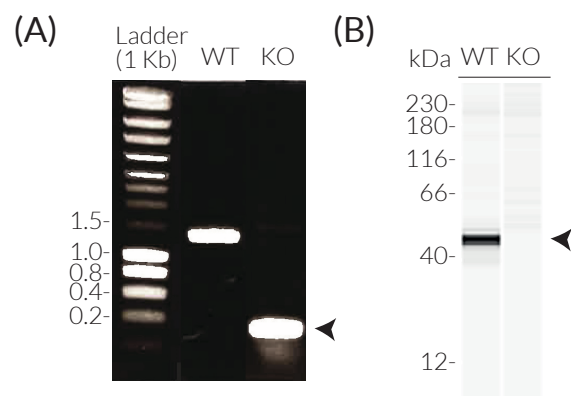
<https://www.invivogen.com/thp1-kocasp4>

For research use only

Version 19J17-ED

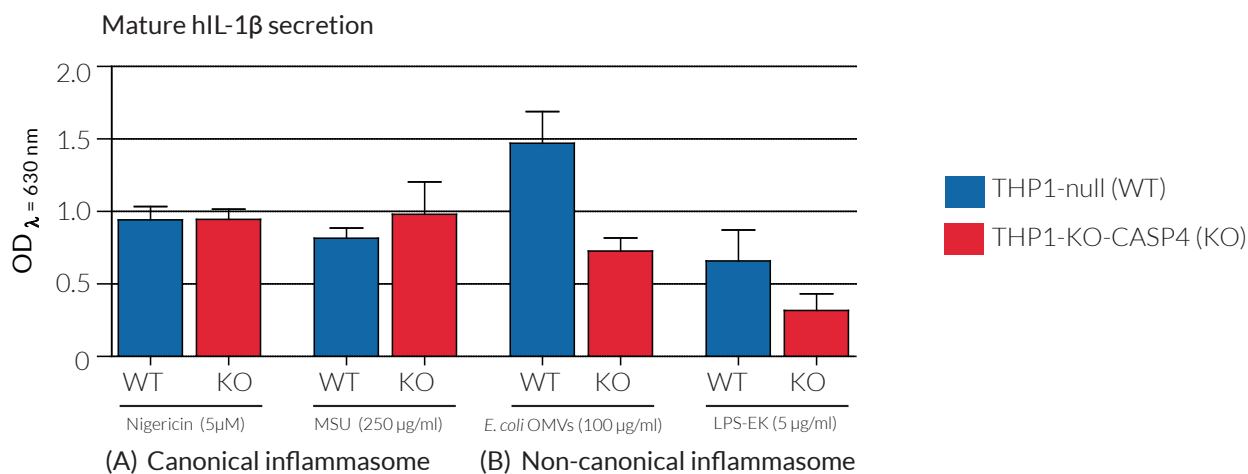
THP1-KO-CASP4 cells were generated from the human monocytic THP1 (WT) cell line through the stable knockout of the caspase-4 (CASP4) gene. The KO status has been verified by DNA sequencing, PCR (Figure 1A), and Western blot (Figure 1B). In THP1-KO-CASP4 cells, mature IL-1 $\beta$  secretion is not affected upon canonical inflammasome activation (Figure 2A), however, it is severely impaired after non-canonical inflammasome activation by *E. coli* outer membrane (OMVs) or transfected LPS (Figure 2B). Of note, due to the presence of CASP5 (functional homolog of CASP4), the secretion of IL-1 $\beta$  is not completely abolished.

## Validation of CASP4 Knockout (KO)



**Figure 1: Validation of CASP4 knockout in THP1-KO-CASP4 cells.** (A) The targeted CASP4 region in THP1-null (WT) and THP1-KO-CASP4 (KO) cells was amplified by PCR. THP1-KO-CASP4 cells feature a biallelic deletion (arrow). (B) Lysates from THP1-null (WT) and THP1-KO-CASP4 (KO) cells were analyzed by Western blot (Wes™) using an anti-human CASP4 antibody, followed by a HRP-conjugated anti-rabbit secondary antibody. The arrow indicates the expected band for the CASP4 protein (48 kDa).

## Functional validation of THP1-KO-CASP4 cells



**Figure 2: Secretion of mature IL-1 $\beta$  by THP1-KO-CASP4 cells upon inflammasome activation.**  $\sim 3 \times 10^5$  THP1-null (blue) and THP1-KO-CASP4 (red) cells were incubated for 3h at 37°C with LPS-EK (1 μg/ml) (priming) and then incubated (activation) with (A) canonical inflammasome inducers, Nigericin (5 μM) or MSU crystals (MSU; 250 μg/ml), and (B) non-canonical inflammasome inducers, *E. coli* outer membrane vesicles (OMVs) (100 μg/ml) or transfected with LPS-EK (5 μg/ml). After 6h, the secretion of mature human (h)IL-1 $\beta$  was assessed in the culture supernatant using HEK-Blue™ IL-1 $\beta$  sensor cells expressing an NF- $\kappa$ B SEAP reporter gene. QUANTI-Blue™ Solution was used to measure SEAP activity. Optical density (OD) was read at 630 nm.

### 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](mailto:info@invivogen.com)