

# Validation data for THP1-KO-CASP4 cells

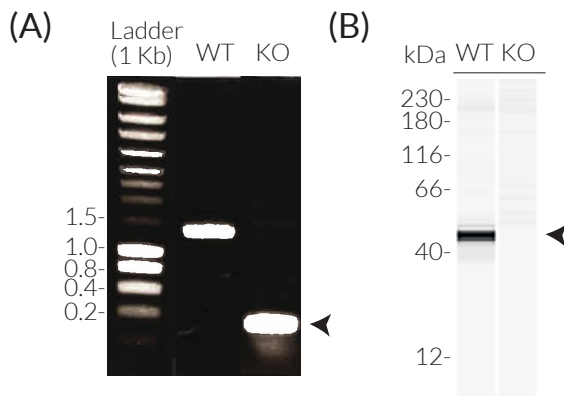
<https://www.invivogen.com/thp1-kocasp4>

For research use only

Version 20H01-NJ

THP1-KO-CASP4 cells were generated from the human monocytic THP1-Null2 (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). Although these cells are fully KO for the *CASP4* protein, they exhibit some IL-1 $\beta$  secretion after non-canonical inflammasome activation by transfected LPS (Figure 2) or *E. coli* outer membrane vesicles (OMVs) (data not shown). This remnant activity may be attributed to a partial rescue of the non-canonical inflammasome response by *CASP5* (functional homolog of *CASP4*). As expected, the mature IL-1 $\beta$  secretion is not affected upon canonical inflammasome activation (Figure 2).

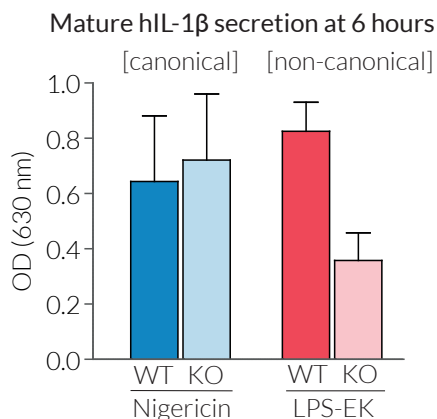
## Validation of *CASP4* Knockout (KO)



**Figure 1: Validation of *CASP4* knockout in THP1-KO-CASP4 cells.**

(A) The targeted *CASP4* region in THP1-Null2 (WT) and THP1-KO-CASP4 (KO) cells was amplified by PCR. THP1-KO-CASP4 cells feature a biallelic deletion (arrow). (B) Lysates from THP1-Null2 (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



WT - THP1-Null2 Cells    KO - THP1-KO-CASP4 Cells

**Figure 2: Altered mature IL-1 $\beta$  secretion by THP1-KO-CASP4 cells upon non-canonical inflammasome activation.**

$\sim 3 \times 10^5$  THP1-Null2 (WT) and THP1-KO-CASP4 cells (KO) were incubated for 3h at 37°C with LPS-EK (1  $\mu$ g/ml) (*priming*) and then incubated (*activation*) with Nigericin (5  $\mu$ M), or transfected LPS-EK (5  $\mu$ g/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 which express 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)



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

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)