

Validation data for THP1-KO-NLRC4 cells

<https://www.invivogen.com/thp1-konlrc4>

For research use only

Version 22J13-NJ

THP1-KO-NLRC4 cells were generated from the human monocytic THP1-Null2 cell line through a deletion encompassing the nucleotide-binding domain (NBD) in the gene encoding NLRC4. NBD is essential for the NLRC4 polymerization and inflammasome assembly. Thus, THP1-KO-NLRC4 cells express an inactive NLRC4 protein, unable to form a functional inflammasome. The KO status has been verified by PCR, Western blot (Figure 1), and functional assays (Figure 2). These cells do not produce mature IL-1 β upon NLRC4/NAIP inflammasome activation using Needle-Tox (LFn-Needle from *B. thailandensis* combined to the anthrax protective antigen). However, they retain the ability to respond to other canonical inflammasome inducers, as well as to non-canonical inflammasome inducers (Figure 2).

Validation of NLRC4 Knockout (KO)

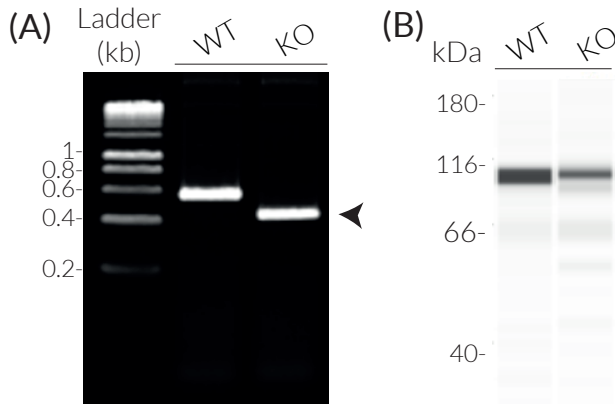


Figure 1: Validation of NLRC4 KO in THP1-KO-NLRC4 cells by PCR and Western blot (WES™).

(A) The targeted NLRC4 region in THP1-Null2 (WT) and THP1-KO-NLRC4 (KO) cells was amplified by PCR. THP1-KO-NLRC4 cells feature a biallelic deletion (arrow).

(B) Lysates from THP1-Null2 (WT) and THP1-KO-NLRC4 (KO) cells were analyzed by Western blot (Wes™) using an NLRC4 antibody (Abcam, ref 201792), followed by a HRP-conjugated anti-rabbit secondary antibody. The arrow indicates the band for the NLRC4 protein (expected size ~115 kDa). Note: THP1-KO-NLRC4 cells feature a deletion encompassing the NLRC4 nucleotide binding domain, which explains the presence of a band in the KO lane.

Functional validation of THP1-KO-NLRC4 cells

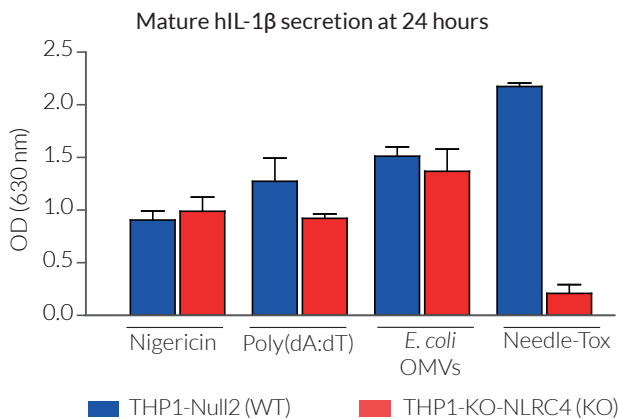


Figure 2: Secretion of mature IL-1 β by THP1-KO-NLRC4 cells and their parental THP1-Null2 cells upon inflammasome activation.

~3x10⁵ THP1-Null2 (WT) and THP1-KO-NLRC4 (KO) cells were incubated for 3h at 37°C with LPS-EK (1 μ g/ml) (priming) and then incubated (activation) with inflammasome inducers: Nigericin (5 μ M), transfected Poly (dA:dT) (1 μ g/ml), *E. coli* outer membrane vesicles (OMVs) (100 μ g/ml), or Needle-Tox (4 ng/ml). After 24h, 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.

Note: Needle-Tox is a combination of LFn-Needle (4 ng/ml) with the anthrax protective antigen (PA) (20 ng/ml). PA allows LFn-Needle translocation into the cytosol.

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



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

 **InvivoGen**
www.invivogen.com