

Validation data for THP1-KO-NLRC4 cells

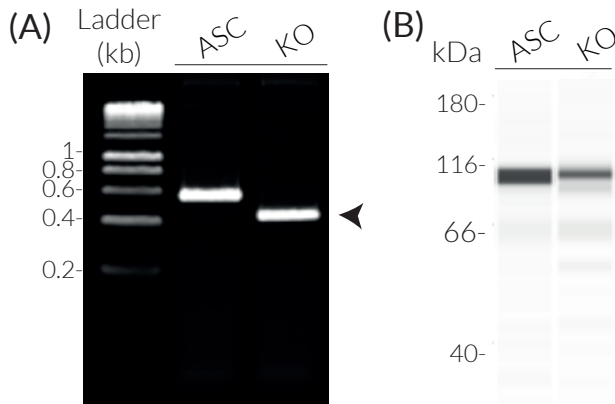
<https://www.invivogen.com/thp1-konlrc4>

For research use only

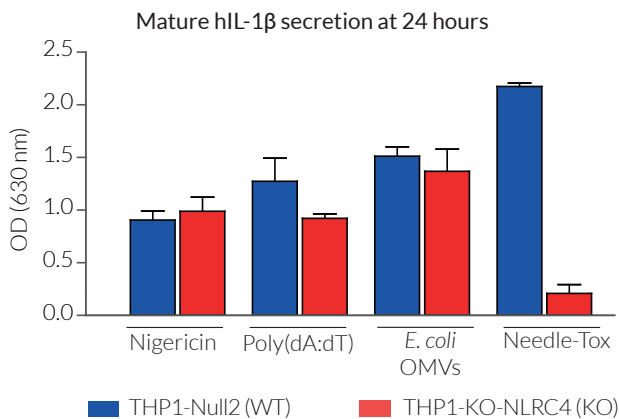
Version 21A11-NJ

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)



Functional validation of THP1-KO-NLRC4 cells



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 3622-3480
E-mail: info@invivogen.com



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

InvivoGen
www.invivogen.com