

# Validation data for THP1-NLRC4 cells

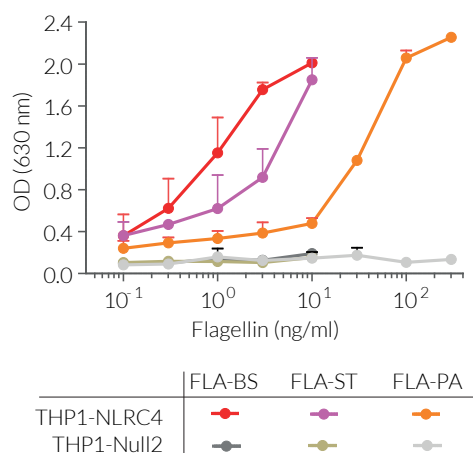
<https://www.invivogen.com/thp1-nlrc4>

For research use only

Version 21B25-NJ

THP1-NLRC4 cells were generated from the human monocytic THP1-Null2 cell line. Substantial Flagellin amounts are required for NLRC4 inflammasome activation in THP-1 cells. THP1-NLRC4 cells stably overexpress NLRC4 and naturally express TLR5. NLRC4 overexpression increases the cell line sensitivity to Flagellins from various Gram-negative bacteria, without impacting the inflammasome response to highly potent NLRC4 inducers, such as Needle-Tox (LFn-Needle from *B. thailandensis* combined with the anthrax protective antigen).

## Inflammasome response in THP1-NLRC4 cells using different Flagellins

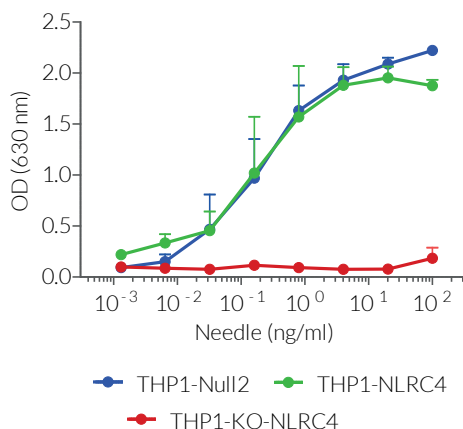


**Figure 1: Secretion of mature IL-1 $\beta$  by THP1-NLRC4 cells and their parental THP1-Null2 cells upon inflammasome activation with Flagellins.**

$\sim 3 \times 10^5$  THP1-Null2 and THP1-NLRC4 cells were incubated for 3h at 37°C with LPS-EK (1  $\mu\text{g/ml}$ ) (*priming*) and then incubated (*activation*) with increasing concentrations of Flagellin-BS (from *B. subtilis*), Flagellin-ST (from *S. typhimurium*), or Flagellin-PA (from *P. aeruginosa*). After 24h, 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-inducible SEAP reporter gene. QUANTI-Blue™ Solution was used to measure SEAP activity. Optical density (OD) was read at 630 nm.

*Note: Ultrapure Flagellin grades were used.*

## Inflammasome response in THP1-NLRC4 cells using Needle-Tox



**Figure 2: Secretion of mature IL-1 $\beta$  by THP1-NLRC4, THP1-KO-NLRC4, and their parental THP1-Null2 cells upon inflammasome activation with Needle-Tox.**

$\sim 3 \times 10^5$  THP1-Null2, THP1-NLRC4, and THP1-KO-NLRC4 cells were incubated for 3h at 37°C with LPS-EK (1  $\mu\text{g/ml}$ ) (*priming*) and then incubated (*activation*) with increasing concentrations of Needle-Tox. After 24h, 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-inducible 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, #171E) (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 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)