

Validation data for LFn-Needle

<https://www.invivogen.com/lfn-needle>

For research use only

Version 21A04-NJ

LFn-Needle is a model of NLRC4/NAIP inflammasome agonist. Needle is a component of the type III secretion systems (T3SS) of intracellular bacteria described as an NLRC4/NAIP ligand. It is fused to the amino-terminal domain of *B. anthracis* lethal factor (LFn). This fusion system, when co-administered with the anthrax toxin's protective antigen (PA), allows intracellular delivery of the bacterial ligand. The combination of LFn-Needle with the anthrax protective antigen (PA) is named Needle-Tox. LFn-Needle specificity for NLRC4 has been verified using THP1-KO-NLRC4 cells (Figure 1). It induces NLRC4-dependent IL-1 β secretion and pyroptosis in a dose-dependent manner (Figure 2).

Functional validation of LFn-Needle using THP1-derived cells

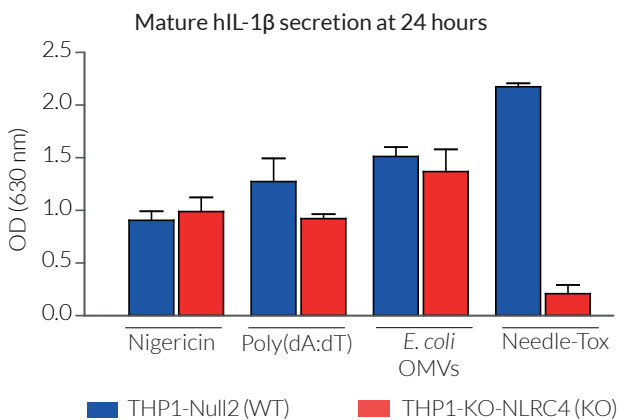


Figure 1: Secretion of mature IL-1 β by THP1-KO-NLRC4 cells and their parental THP1-Null2 cells upon inflammasome activation.

$\sim 3 \times 10^5$ 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.

Human NLRC4 inflammasome response to LFn-Needle

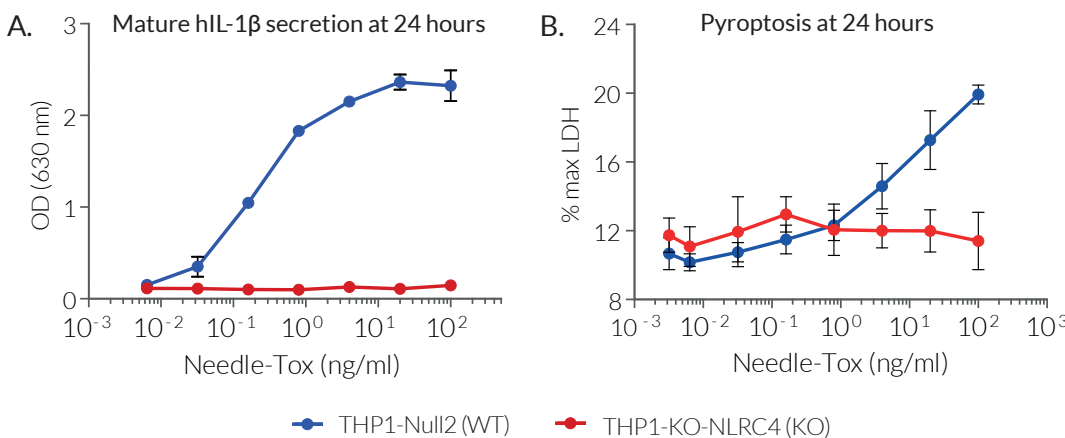


Figure 2: Secretion of mature IL-1 β and pyroptosis upon THP1-derived cell activation using LFn-Needle.

$\sim 3 \times 10^5$ THP1-Null2 (blue) and THP1-KO-NLRC4 cells (red) were incubated for 3h at 37°C with LPS-EK (1 μ g/ml) (priming) and then incubated (activation) with Needle-Tox (0.006-100 ng/ml). After 24h, mature human (h) IL-1 β was assessed in the culture supernatant using HEK-Blue™ IL-1 β sensor (A), and cell death was assessed using the lactate dehydrogenase (LDH) assay (B).

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