

Validation data for THP1-KO-ASC cells

<https://www.invivogen.com/thp1-ko-kd-asc>

For research use only

Version 20F24-NJ

THP1-KO-ASC cells were generated from the human monocytic THP1-Null2 (WT) cell line through the stable knockout of the ASC gene. The KO status has been verified by DNA sequencing, PCR, Western blot (Figure 1), and functional assays (Figure 2). Mature IL-1 β secretion is abolished in THP1-KO-ASC cells (KO) after activation with inducers of the ASC-dependent NLRP3 (i. e. Nigericin, MSU crystals, or Alum Hydroxide) and AIM2 (i.e. Poly (dA:dT)) canonical inflammasomes, when compared to the parental (WT) cell line (Figure 2A). The NLRP3-caspase-1 dependent maturation of IL-1 β is abolished upon indirect activation of NLRP3 by induction of the non-canonical inflammasome by *E. coli* outer membrane vesicles (OMVs) (Figure 2A). Furthermore, pyroptosis is abrogated in THP1-KO-ASC cells as shown by an LDH assay (Figure 2B).

Validation of ASC Knockout (KO)

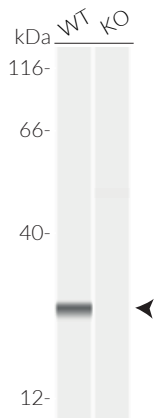
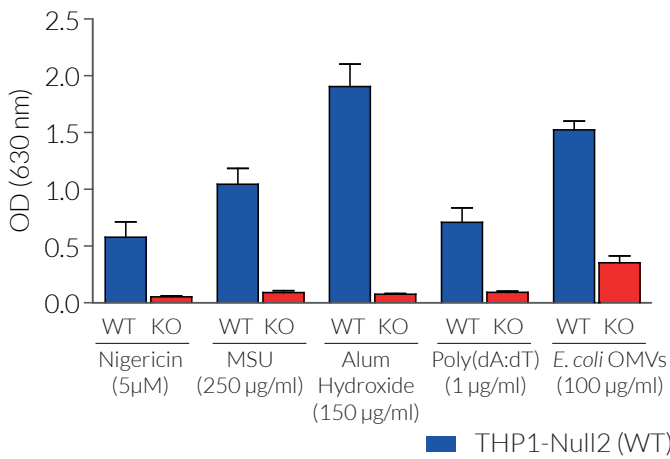


Figure 1: Validation of ASC KO in THP1-KO-ASC cells by Western blot (WES™).

Lysates from THP1-Null2 (WT) and THP1-KO-ASC (KO) cells were analyzed using an anti-human ASC antibody, followed by a HRP-conjugated anti-rabbit secondary antibody. The arrow indicates the expected band for the ASC protein (27 kDa).

Functional validation of THP1-KO-ASC cells

(A) Mature hIL-1 β secretion at 24 hours



(B) Cell death at 6 hours

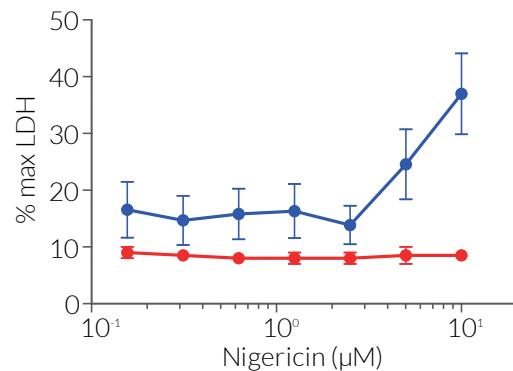


Figure 2: Absence of mature IL-1 β secretion and pyroptosis by THP1-KO-ASC cells upon inflammasome activation.

$\sim 3 \times 10^5$ THP1-Null2 (blue) and THP1-KO-ASC cells (red) were incubated for 3h at 37°C with LPS-EK (1 μ g/ml) (priming) and then incubated (activation) with Nigericin (A - 5 μ M; B - 0.15-10 μ M), MSU crystals (MSU; 250 μ g/ml), Alum Hydroxide (150 μ g/ml), transfected Poly (dA:dT) (1 μ g/ml), or *E. coli* outer membrane vesicles (OMVs) (100 μ g/ml). After 24h, (A) 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. (B) After a 6 hour incubation with Nigericin, cell death was assessed using the lactate dehydrogenase (LDH) assay.

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



Any questions about our cell lines?
[Visit our FAQ page.](#)

 **InvivoGen**
www.invivogen.com