

Validation data for RAW-ASC KO-CASP11 cells

<https://www.invivogen.com/raw-kocasp11>

For research use only

Version 20G27-NJ

RAW-ASC KO-CASP11 cells were generated from the murine RAW-ASC macrophage cell line through the stable knockout of the murine caspase-11 (CASP11) gene. The KO status has been verified by PCR, Western blot (Figure 1A-B), and functional assays (Figure 2). While mature IL-1 β secretion in RAW-ASC KO-CASP11 cells is comparable to the parental cell line upon canonical inflammasome (NLRP3 and AIM2) activation (Figure 2A), this response is impaired upon activation of the non-canonical caspase-11 inflammasome using *E. coli* outer membrane vesicles (OMVs) or transfected LPS-EK (Figure 2B).

Validation of CASP11 Knockout (KO)

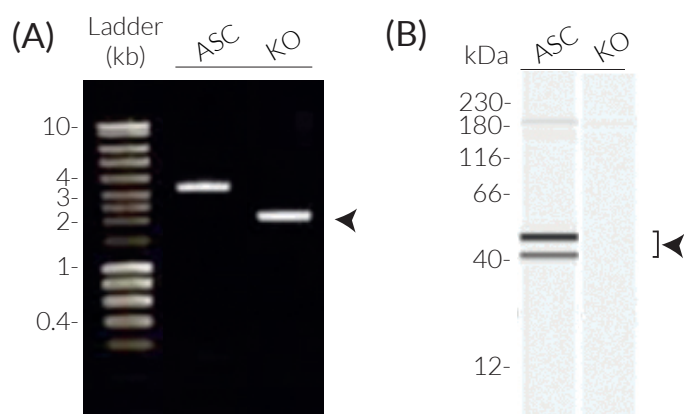


Figure 1: Validation of CASP11 KO in RAW-ASC KO-CASP11 cells by PCR and Western blot (WES™).

(A) The targeted CASP11 region in RAW-ASC (WT) and RAW-ASC KO-CASP11 (KO) was amplified by PCR. RAW-ASC KO-CASP11 feature a biallelic deletion (arrow). (B) Lysates from RAW-ASC (ASC) and RAW-ASC KO-CASP11 (KO) cells were analyzed using an anti-mouse CASP11 antibody, followed by a HRP-conjugated anti-rabbit secondary antibody. The arrow indicates the bands for the pro-caspase 11 protein isoforms (expected sizes ~38 and 43 kDa).

Functional validation of RAW-ASC KO-CASP11 cells

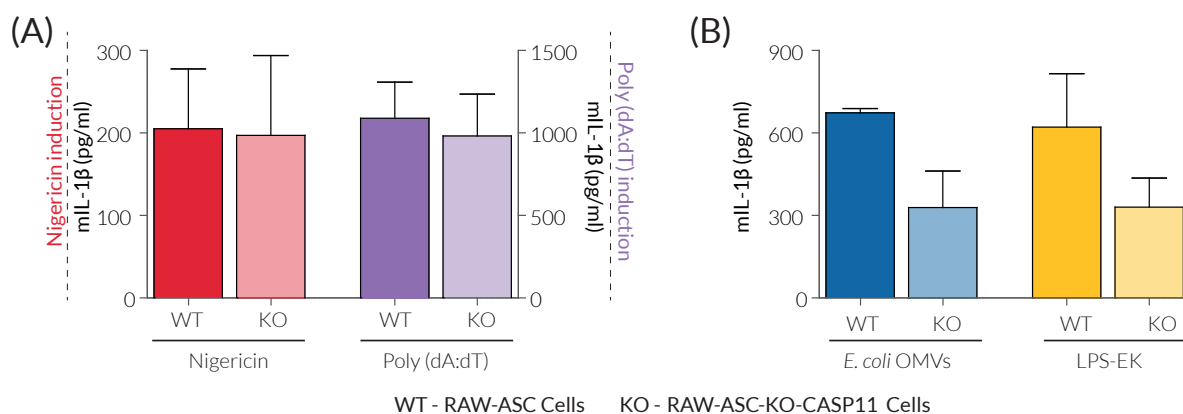


Figure 2: Secretion of mature IL-1 β by RAW-ASC KO-CASP11 cells and their parental RAW-ASC cells upon inflammasome activation.

~2x10⁵ RAW-ASC (WT) and RAW-ASC-KO-CASP11 (KO) cells were incubated for 3h at 37°C with Pam3CSK4 (100 ng/ml) (priming) and then incubated (activation) with (A) canonical inflammasome inducers Nigericin (5 μ M) or transfected Poly(dA:dT) (1 μ g/ml), and (B) non-canonical inflammasome inducers, *E. coli* outer membrane vesicles (OMVs) (100 μ g/ml) or transfected LPS-EK (5 μ g/ml). After 24h, the secretion of mature murine (m)IL-1 β was assessed in the culture supernatant using an ELISA assay.

Note: For non-canonical inflammasome activation, cells were pre-primed with recombinant murine IFN- γ (10 ng/ml) overnight at 37°C before priming.

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