

Validation data for RAW-ASC cells

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

For research use only

Version 20G07-ED

RAW-ASC cells were generated by stable transfection of the murine ASC gene into the murine RAW 264.7 macrophage cell line which is otherwise ASC-deficient. The expression of ASC in RAW-ASC cells was verified by Western blot (Figure 1) and functional assays (Figure 2). RAW-ASC cells produce and secrete mature IL-1 β upon activation of canonical (NLRP3 and AIM2) (Figure 2A) and non-canonical (caspase-11) inflammasomes (Figure 2B).

Validation of RAW-ASC cells

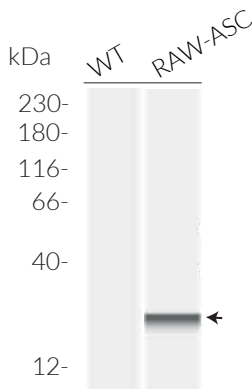


Figure 1: Validation of ASC expression in RAW-ASC cells by Western blot (WES™). Lysates from RAW 264.7 (WT) and RAW-ASC cells were analyzed using an anti-mouse ASC antibody, followed by a HRP-conjugated anti-rabbit secondary antibody. The arrow indicates the expected band for the ASC protein (27 kDa).

Functional validation of RAW-ASC cells

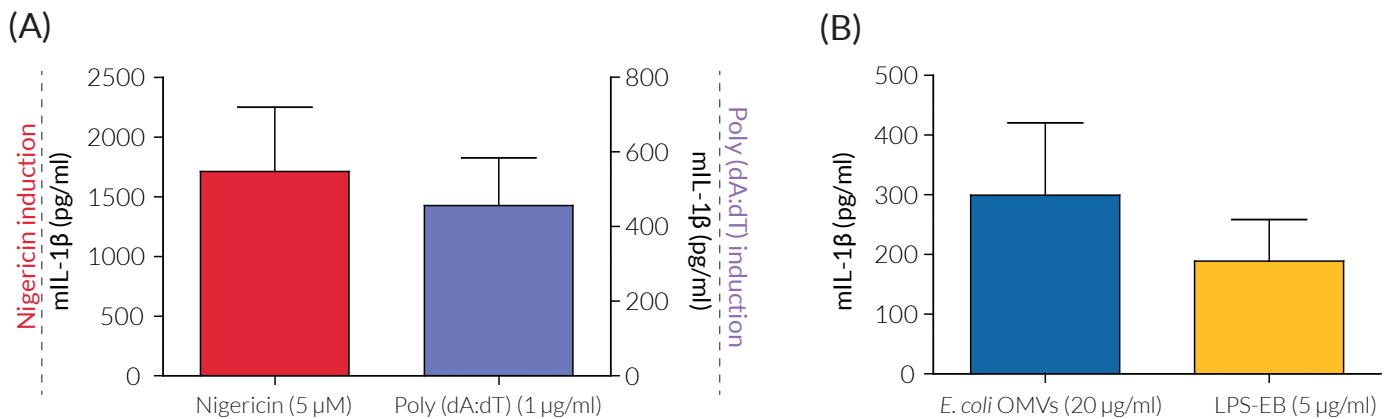


Figure 2: Secretion of mature IL-1 β by RAW-ASC cells upon inflammasome activation. $\sim 1 \times 10^6$ RAW-ASC cells were incubated for 3h at 37°C with Pam3CSK4 (100 ng/ml) (priming) and then incubated (activation) with (A) canonical inflammasome inducers Nigericin (5 μ M) or transfected Poly (dA:dT) (1 μ g/ml), and (B) non-canonical inflammasome inducers, *E. coli* outer membrane vesicles (OMVs) (20 μ g/ml) or transfected LPS-EB (5 μ g/ml). After 6h, the secretion of mature 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