

Validation data for *E. coli* OMVs

<https://www.invivogen.com/ecoli-omvs>

For research use only

Version 19A17-ED

Outer membrane vesicles (OMVs) are secreted, small, non-replicative, immunogenic spherical bodies that contain many pathogen-associated molecular patterns (PAMPs) from the parent bacterium such as DNA, RNA, peptidoglycan, lipoproteins, toxins, and lipopolysaccharides (LPS). Invivogen has isolated OMVs with a diameter between 40-60 nm from late-stationary phase cultures of the commonly used *Escherichia coli* BL21 strain (Figure 1). In particular, LPS one of the most abundant constituents of OMVs, is recognized by and activates caspase-11 (human homologs caspase 4/5) forming a non-canonical inflammasome, which in turn can also activate the canonical inflammasome pathway. These events are characterized by the release of the alarmin HMGB1 upon membrane rupture (Figure 2) and the release of pro-inflammatory cytokines such as IL-1 β (Figure 3). Concomitantly, the activation of the caspase 11-4/5 inflammasome complex by *E. coli* OMVs leads to pyroptotic cell death through the action of gasdermin D (GSDMD) (Figure 4).

Characterization of *E. coli* OMVs

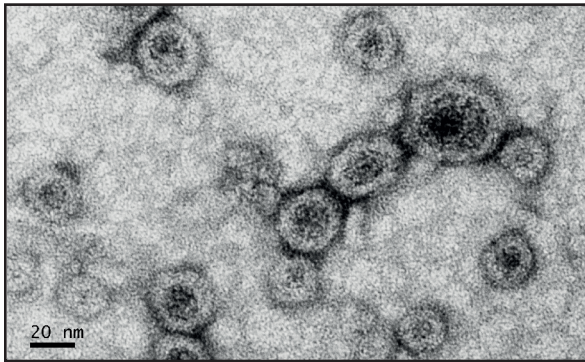


Figure 1: *E. coli* OMVs were prepared for electron microscopy using conventional negative staining procedures. *E. coli* OMVs (20 μ l) were absorbed onto Formvar carbon-coated grids for 2 min, blotted, and then negatively stained with 1% uranyl acetate for 1 min. Grids were then examined with a transmission electron microscope (Jeol JEM-1400, JEOL Inc, Peabody, MA, USA) at 80 kV. Images were acquired with a digital camera (Gatan Orius, Gatan Inc, Pleasanton, CA, USA) using a range of magnifications (20 nm magnification seen to the left).

Evaluation of the efficiency of *E. coli* OMVs on a pyroptosis reporter cell line

THP1-HMGB1-Lucia™ cells are designed to study inflammasome-mediated cell death, characterized by the release of cytoplasmic components upon cell membrane rupture. THP1-HMGB1-Lucia™ cells stably express a fusion protein, in which the C-terminus of the alarmin HMGB1 (High Mobility Group Box-1) is fused to Lucia luciferase. Levels of the fusion protein in the supernatant can be readily monitored using the detection reagent QUANTI-Luc™. When these cells are activated with *E. coli* OMVs there is a dose-dependent response seen when measuring luciferase activity. The optimal concentration of *E. coli* OMVs is approximately 20 μ g/ml.

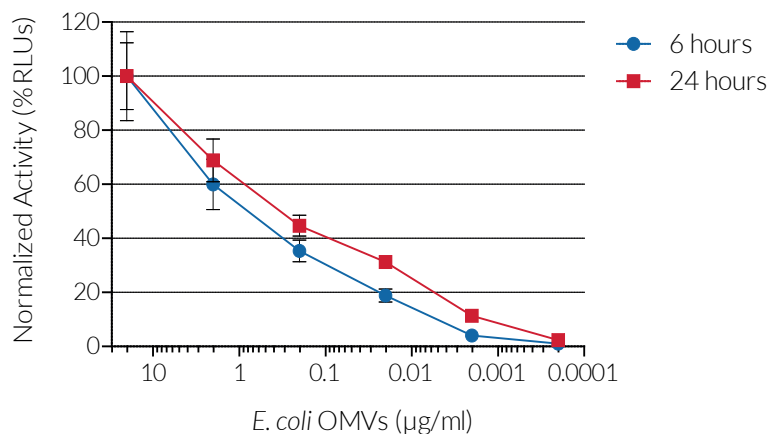


Figure 2: THP1-HMGB1-Lucia™ cells were incubated with 0.2 ng/ml-20 μ g/ml of *E. coli* OMVs for 6 (blue) and 24 hrs (red). After the incubation, Lucia luciferase activity was determined by measuring relative light units (RLUs) in a luminometer using the QUANTI-Luc™ detection reagent. Results are presented as normalized activity (percentage % RLUs).

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

 **InvivoGen**
www.invivogen.com

Evaluation of *in vitro* induction of the caspase 11-4/5 inflammasome

After endocytosis of OMVs into the cytosol, LPS on the surface of the OMVs activates the caspase 11-4/5 inflammasome leading to pyroptotic cell death and the release of cytokines such as IL-1 β . Treatment of murine bone-marrow derived macrophages (BMDMs) with *E. coli* OMVs shows a clear dose-dependent response for the release of the biological marker for cellular cytotoxicity lactate dehydrogenase (LDH) (Figure 3A) and the pro-inflammatory cytokine IL-1 β (Figure 3B). Additionally, knockout (KO)-caspase11 murine BMDMs show very minimal release of LDH and IL-1 β , highlighting that caspase 11-4/5 inflammasome -induced cell death by OMVs is highly specific.

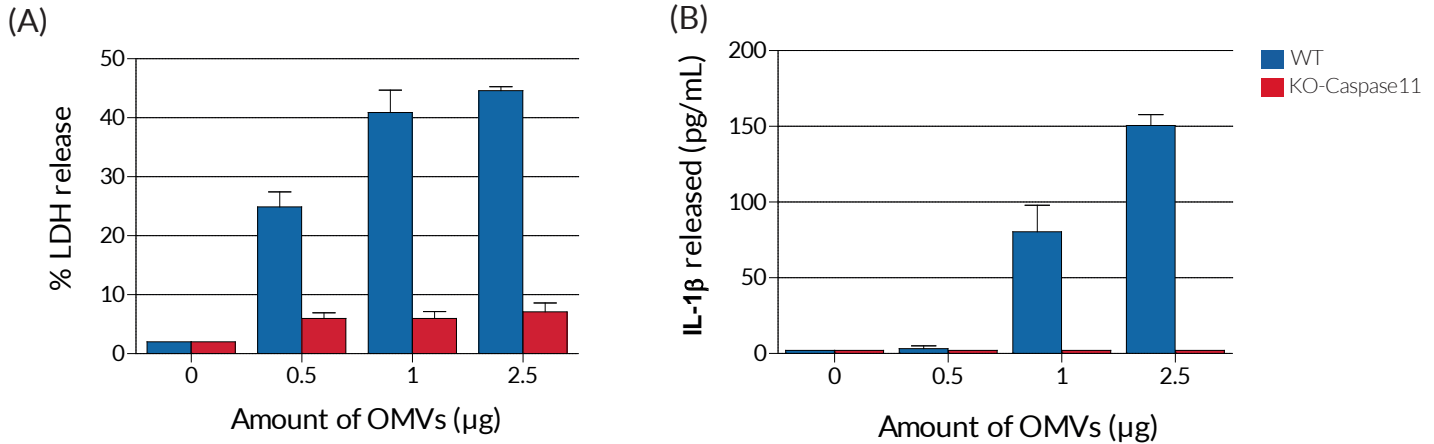


Figure 3: Bone marrow-derived macrophages (BMDMs) from wild type (WT) (blue) and caspase-11 knockout (KO) (red) mice were treated with a range of different concentrations (from 0.5 to 2.5 μg/250 000 cells) of *E. coli* OMVs (prepared in PBS) for 16-20 hours. After the incubation, the cell supernatants were used for both a (A) lactate dehydrogenase (LDH) assay and an (B) ELISA for measuring the release of IL-1 β .

Evaluation of *in vivo* induction of the caspase 11-4/5 inflammasome

E. coli OMVs can also be used for *in vivo* studies that have been designed to mimic pathogenic bacterial infection of a host. These studies are essential to elucidate the role of the caspase 11-4/5 inflammasome in the outcome of the infection. We recommend to use *E. coli* OMVs InvivoFit™ grade, as these have been thoroughly tested to ensure sterility. Treatment of mice intraperitoneally with *E. coli* OMVs InvivoFit™ shows a dose-dependent response. Additionally, this response is completely dependent on the presence of gasdermin-D, a key player in pyroptotic cell death.

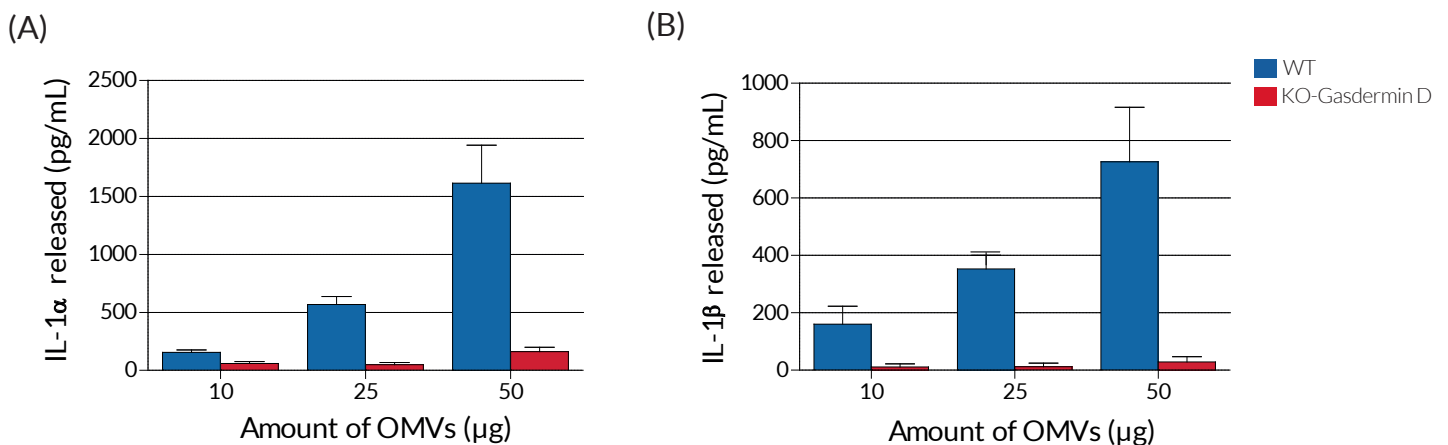


Figure 4: Wild type (WT) (blue) and Gasdermin-D knock out (KO) (red) mice (20-25g) were injected with 150 μg of poly IC (LMW) for the priming stage of inflammasome activation, and then left for 6 hours. Following this, a range of different concentrations from 10-50 μg of *E. coli* OMVs (prepared in PBS -/-) were injected intraperitoneally into the mice. After an additional 5 hours, peritoneal (A) IL-1 α and (B) IL-1 β cytokine levels were assayed using ELISA kits.

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