

Validation data for A549-ASCoV2 cells

<https://www.invivogen.com/a549-ascg-cov>

For research use only

Version 22H04-NJ

A549-ASCoV2 cells are designed as a control cell line for A549-ASCoV2-NLRP1 cells to study the activation of the NLRP1 inflammasome in real time, in the context of SARS-CoV-2 infection. They are derived from A549-ASC cells, a human A549 lung carcinoma epithelial cell line featuring an NF- κ B-inducible ASC::GFP reporter gene. Additionally, A549-ASCoV2 cells stably express the SARS-CoV-2 receptors genes, human ACE2 and *TMPRSS2*. The expression of *hTMPRSS2* was confirmed using quantitative RT-PCR (Figure 1) and of hACE2 using flow cytometry (Figure 2). Western blot analysis confirmed stable ASC::GFP expression and absence of endogenous NLRP1 expression (Figure 3). The formation of ASC specks upon NLRP1 inflammasome activation by SARS-CoV-2 viral infection was monitored using fluorescence microscopy (Figure 4), and subsequent pyroptotic cell death using the LDH-release assay (Figure 5). As expected, the lack of endogenous NLRP1 expression in A549-ASCoV2 cells precludes the formation of single ASC specks and cell death.

Validation of *TMPRSS2* overexpression

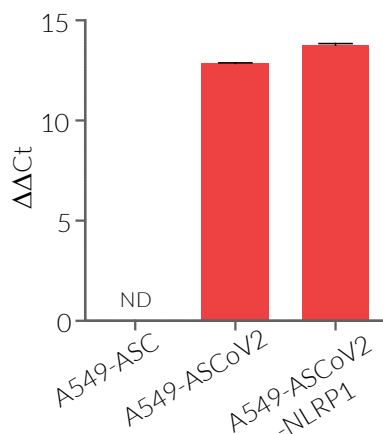


Figure 1: Human *TMPRSS2* mRNA expression in A549-ASC, A549-ASCoV2, A549-ASCoV2-NLRP1 cells. Total mRNA was extracted from $\sim 1 \times 10^6$ cells for each cell line. *TMPRSS2* mRNA was amplified using quantitative RT-PCR. Data are represented as $\Delta\Delta CT$ comparing *TMPRSS2* expression to a housekeeping gene. ND: not detected.

Validation of ACE2 surface expression

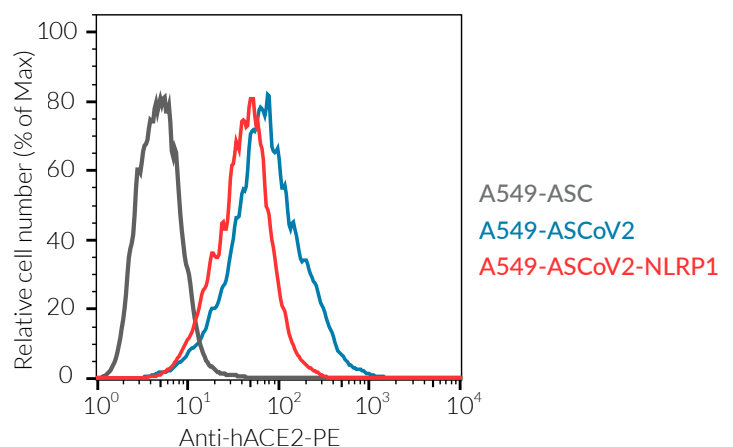


Figure 2: Surface expression of hACE2 by A549-ASC, A549-ASCoV2, A549-ASCoV2-NLRP1 cells. Each cell line was incubated with Anti-ACE2-PE antibody for 1 hour at 4°C. Cell surface staining was analyzed by flow cytometry.

Validation of NLRP1 and ASC::GFP expression

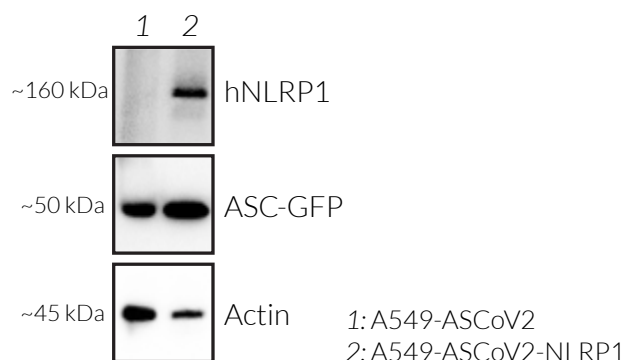


Figure 3: Validation of the expression of ASC::GFP and human NLRP1. Lysates from A549-ASCoV2 (1) and A549-ASCoV2-NLRP1 (2) cells were analyzed by western blot using an anti-human NLRP1 and an anti-human ASC antibody, followed by HRP conjugated secondary antibody. Actin was used as a loading control.

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 Asia: +852 3622-3480
E-mail: info@invivogen.com

Monitoring of ASC speck formation upon NLRP1 inflammasome activation

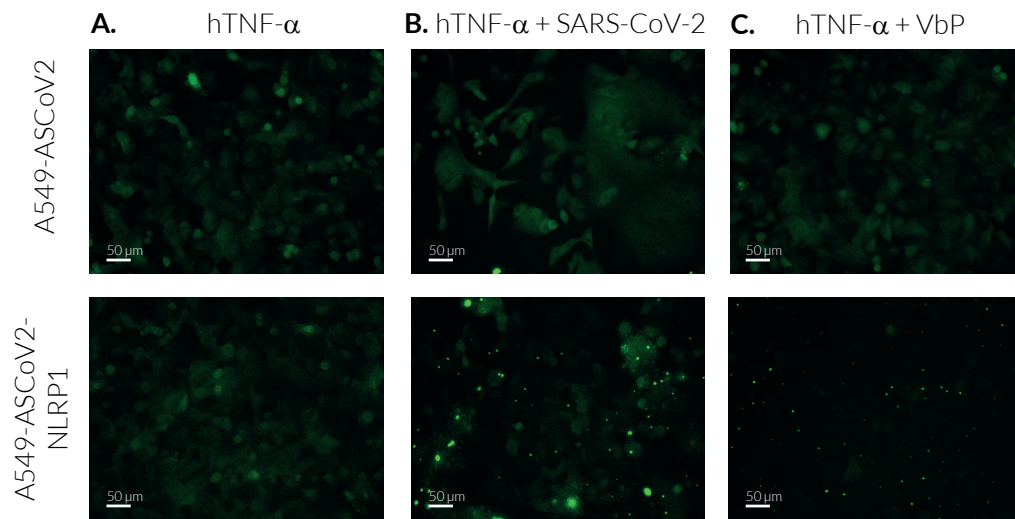


Figure 4: SARS-CoV-2 virus and Val-boroPro induce ASC speck formation in A549-ASCoV2-NLRP1 cells. A549-ASCoV2 cells and A549-ASCoV2-NLRP1 cells were incubated with 0.5 ng/ml human TNF- α overnight at 37°C, 5% CO₂ (A-C). The following day, the cells were incubated in test medium containing SARS-CoV-2 particles (MOI of 0.1) for 1 hour, then for 24 hours in fresh test medium (B). Alternatively, cells were incubated in test medium containing 10 μ M Val-boroPro (VbP) for 24 hours (C). The ASC::GFP expression and ASC speck formation were monitored using fluorescence microscopy. Scale bar: 50 μ m.

Monitoring of pyroptotic cell death upon NLRP1 inflammasome activation

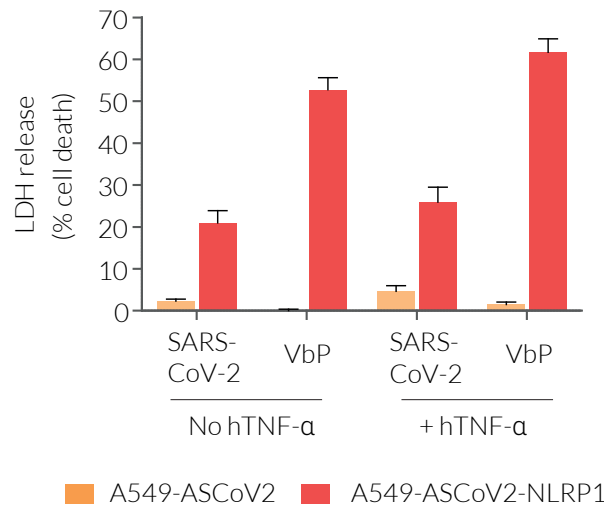


Figure 5: SARS-CoV-2 virus and Val-boroPro induce pyroptotic cell death in A549-ASCoV2-NLRP1 cells. A549-ASCoV2 cells and A549-ASCoV2-NLRP1 cells were incubated with/without 0.5 ng/ml human TNF- α overnight at 37°C, 5% CO₂. The following day, the cells were incubated in test medium containing SARS-CoV-2 particles (MOI of 0.1) for 1 hour, then for 24 hours in fresh test medium. Alternatively, cells were incubated in test medium containing 10 μ M Val-boroPro (VbP) for 24 hours. Cell death was assessed using the lactate dehydrogenase (LDH) assay. Data is shown as percentage of cell death (mean \pm SEM).

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 Asia: +852 3622-3480
 E-mail: info@invivogen.com