

# Validation data for A549-ASCoV2-NLRP1 cells

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

For research use only

Version 22H02-NJ

A549-ASCoV2-NLRP1 cells are designed to study NLRP1 inflammasome activation upon SARS-CoV-2 infection. They are derived from A549-ASC cells, a human A549 lung carcinoma epithelial cell line featuring an NF- $\kappa$ B-inducible ASC::GFP reporter gene. Additionally, A549-ASCoV2-NLRP1 cells stably express the human *NLRP1*, as well as 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 *NLRP1* and ASC::GFP 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).

## Validation of *TMPRSS2* overexpression

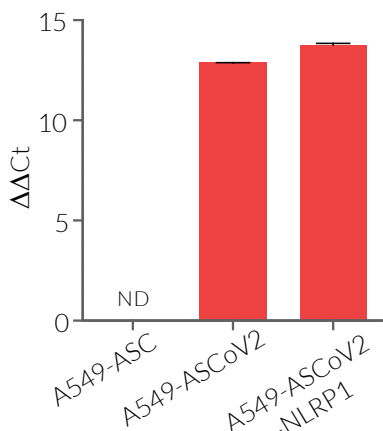


Figure 1: Human *TMPRSS2* mRNA expression in A549-ASC, A549-ASCoV2, A549-ASCoV2-NLRP1 cells. Total mRNA was extracted from  $\sim 1 \times 10^6$  cells for each cell line. *TMPRSS2* mRNA was amplified using quantitative RT-PCR. Data are represented as  $\Delta\Delta C_t$  comparing *TMPRSS2* expression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

## Validation of ACE2 surface expression

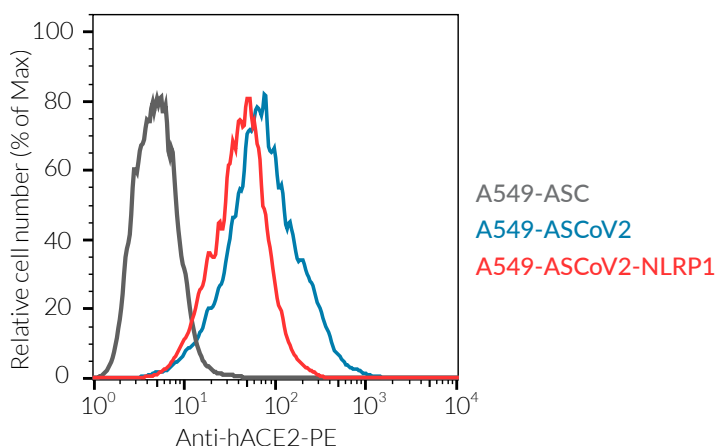


Figure 2: Surface expression of *hACE2* by A549-ASC, A549-ASCoV2, A549-ASCoV2-NLRP1 cells. Each cell line was incubated with Anti-ACE2-PE antibody for 1 hour at 4°C. Cell surface staining was analyzed by flow cytometry.

## Validation of NLRP1 and ASC::GFP expression

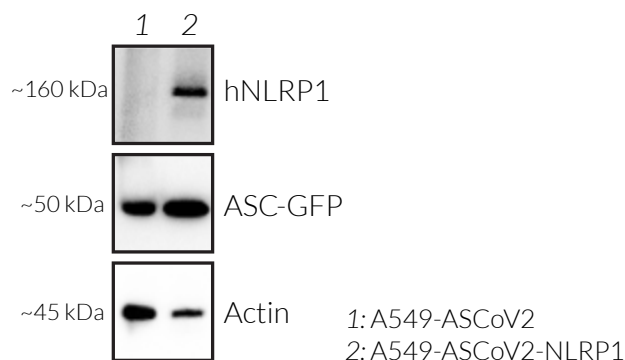
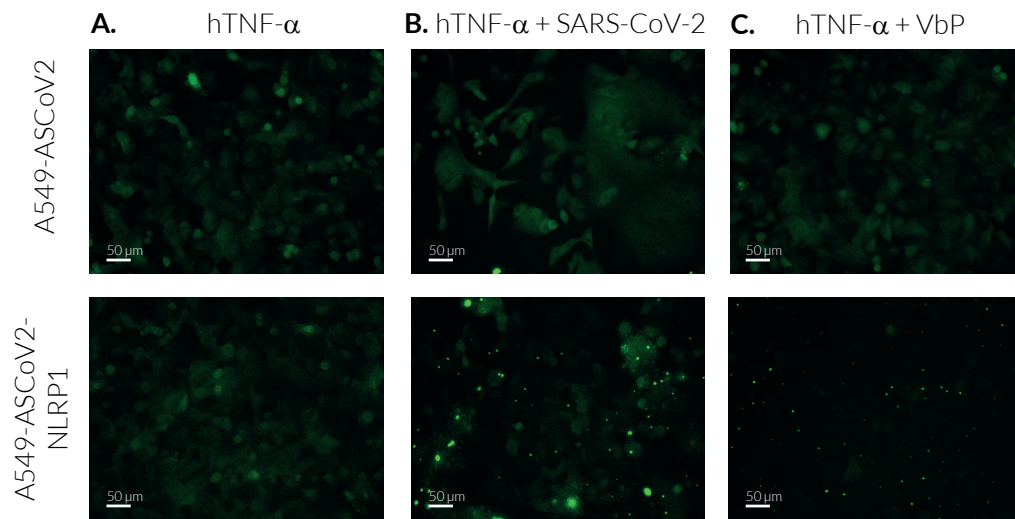


Figure 3: Validation of the expression of ASC::GFP and human *NLRP1*. Lysates from A549-ASCoV2 (1) and A549-ASCoV2-NLRP1 (2) cells were analyzed by western blot using an anti-human *NLRP1* and an anti-human ASC antibody, followed by HRP conjugated secondary antibody. Actin was used as a loading control.

## TECHNICAL SUPPORT

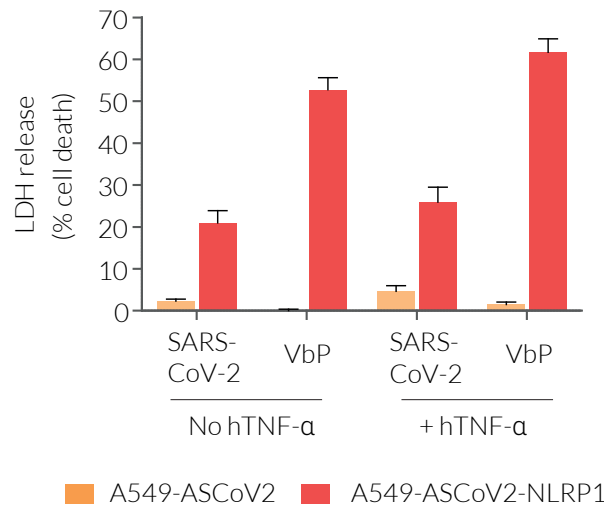
InvivoGen USA (Toll-Free): 888-457-5873  
InvivoGen USA (International): +1 (858) 457-5873  
InvivoGen Europe: +33 (0) 5-62-71-69-39  
InvivoGen Asia: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)

### 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- $\alpha$  overnight at 37°C, 5% CO<sub>2</sub> (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  $\mu$ M Val-boroPro (VbP) for 24 hours (C). The ASC::GFP expression and ASC speck formation were monitored using fluorescence microscopy. Scale bar: 50  $\mu$ 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- $\alpha$  overnight at 37°C, 5% CO<sub>2</sub>. 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  $\mu$ 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](mailto:info@invivogen.com)