

# Validation data for A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells

<https://www.invivogen.com/a549-dual-hace2tmprss2-cells>

For research use only

Version 21B08-NJ

A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells were generated from the A549-Dual™ KO-MDA5 lung carcinoma cell line through stable integration of the human ACE2 and TMPRSS2 genes. The A549 cell line is a well-characterized cellular model of the alveolar basal epithelial lung cells for the study of respiratory infections. ACE2 and TMPRSS2 overexpression in A549-Dual™ hACE2-TMPRSS2 and A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells has been verified by RT-qPCR (Figure 1), and ACE2 expression has been verified by cell surface staining (Figure 2). The biallelic deletion of the MDA5 gene has been verified by PCR (Figure 3). A549-Dual™ KO-MDA5 hACE2-TMPRSS2 and their parental cell line are highly permissive to infection with pseudotyped lentiviral particles expressing the SARS-CoV-2 Spike (G614) protein (Figure 4). A549-Dual™ KO-MDA5 hACE2-TMPRSS2 and their parental cell line display highly similar IRF and NF-κB activity in response to nucleic acid stimulation (Figure 5).

## Validation of ACE2 and TMPRSS2 overexpression by RT-qPCR      Validation of ACE2 surface expression by FACS

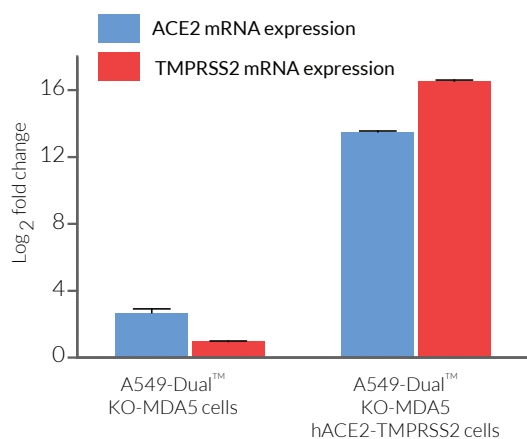


Figure 1: Human ACE2 and TMPRSS2 mRNA expression in A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells.

Total mRNA was extracted from ~1x10<sup>6</sup> A549-Dual™ hACE2-TMPRSS2 (parental) and A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells. ACE2 and TMPRSS2 mRNA were amplified using quantitative RT-qPCR. Data are represented as the log<sub>2</sub> fold change comparing ACE2 or TMPRSS2 expression to a housekeeping gene.

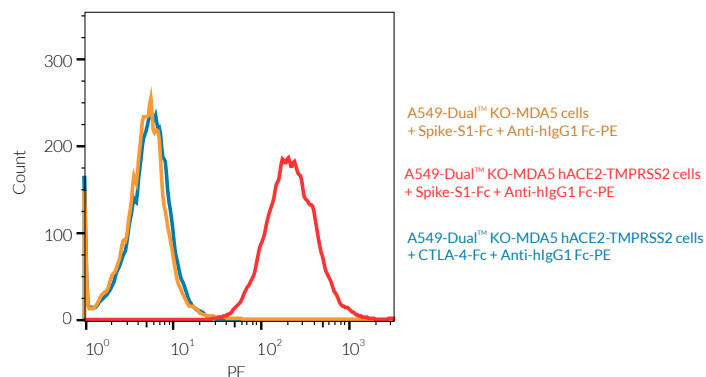


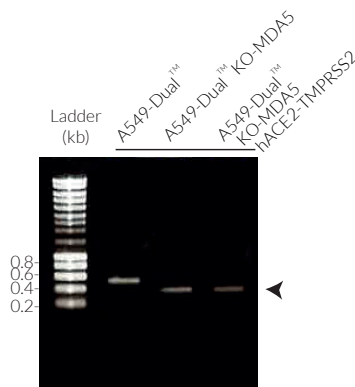
Figure 2: Surface expression of hACE2 by A549-Dual™ KO-MDA5 -hACE2-TMPRSS2 cells.

~2x10<sup>5</sup> A549-Dual™ KO-MDA5 and A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells were incubated with 1 μg of Spike-S1-Fc or CTLA-4-Fc fusion proteins for 1h at 4°C. Cells were then washed and incubated with 0.5 μg of a goat anti-hlgG1-Fc antibody coupled to PE for 1h at 4°C. Cell surface staining was analyzed by flow-cytometry.

## Validation of MDA5 knock-out by PCR

Figure 3: Validation of MDA5 KO in A549-Dual™ KO-MDA5-hACE2-TMPRSS2 cells.

The targeted MDA5 gene in A549-Dual™ (WT), A549-Dual™ KO-MDA5, and A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells was amplified by PCR. A549-Dual™ KO-MDA5, and A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells feature a biallelic deletion (arrow).



### TECHNICAL SUPPORT

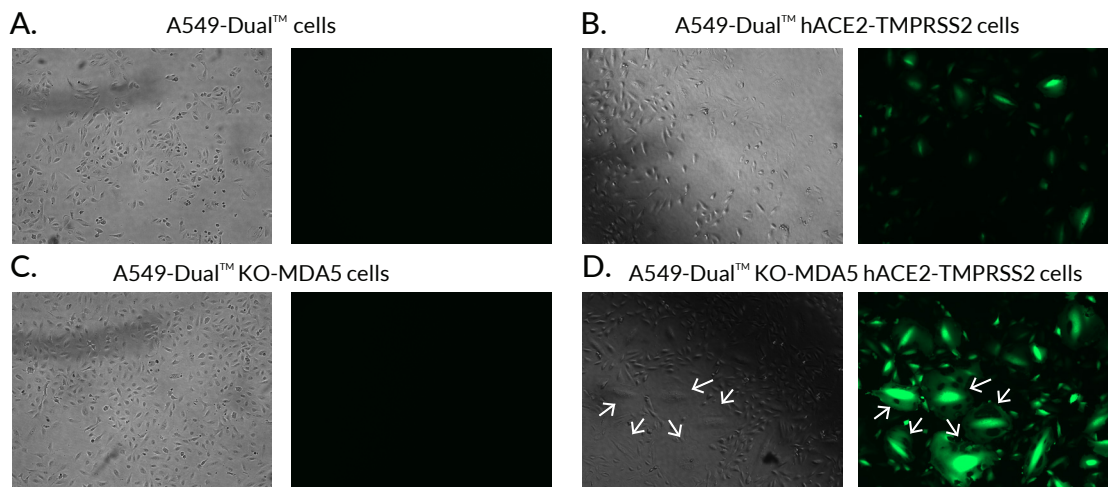
InvivoGen USA (Toll-Free): 888-457-5873  
 InvivoGen USA (International): +1 (858) 457-5873  
 InvivoGen Europe: +33 (0) 5-62-71-69-39  
 InvivoGen Hong Kong: +852 3-622-34-80  
 E-mail: [info@invivogen.com](mailto:info@invivogen.com)



Any questions about our cell lines?  
 Visit our FAQ page.

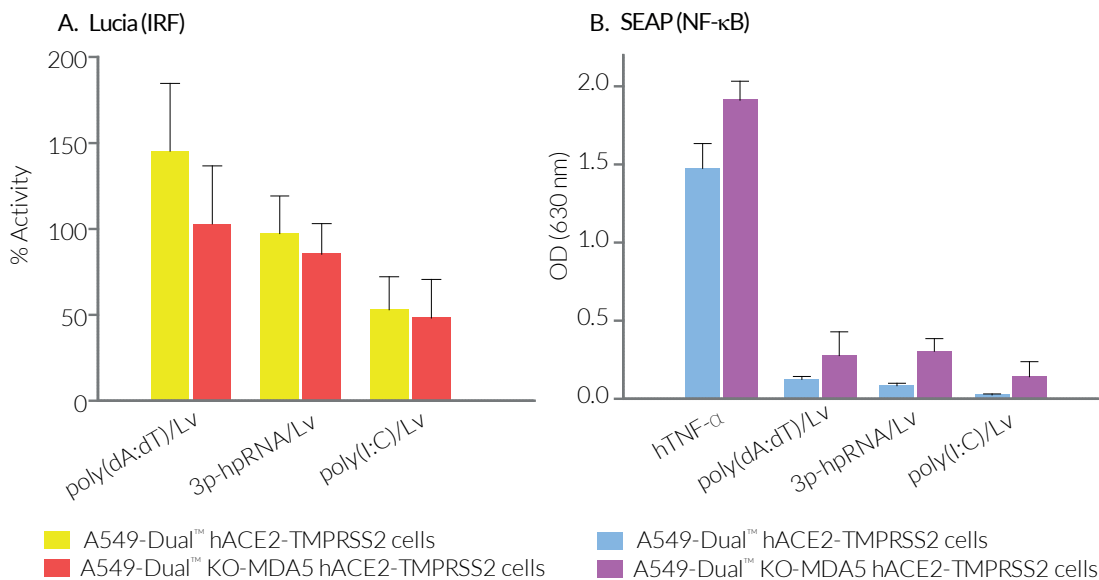


## Infection of A549-Dual™ -derived cells by SARS-CoV-2 Spike (G614) pseudotyped lentiviral particles



**Figure 4: Infection of A549-Dual™-derived cells by Spike (G614) pseudotyped lentiviral particles.**  $\sim 1 \times 10^4$  (A) A549-Dual™, (B) A549-Dual™ hACE2-TMPRSS2, (C) A549-Dual™ KO-MDA5, and (D) A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells were cultured in the presence of SARS-CoV-2 Spike G614-variant pseudotyped GFP lentiviral particles. After 72h, the transduction efficiency of the Spike pseudotyped GFP particles was evaluated by fluorescence microscopy. Syncytia are indicated with white arrows.

## Functional validation of IRF and NF-κB reporter systems in A549-Dual™ KO-MDA5 ACE2-TMPRSS2 cells



**Figure 5: Activation of A549-Dual™-derived cells.**  $5 \times 10^5$  A549-Dual™ hACE2-TMPRSS2 or A549-Dual™ KO-MDA5 hACE2-TMPRSS2 cells were incubated with 100 ng/ml poly(dA:dT) complexed with LyoVec™, 100 ng/ml 3p-hpRNA complexed with LyoVec™, 100 ng/ml poly(I:C) HMW complexed with LyoVec™, or 10ng/ml hTNF-α. After overnight incubation, IRF responses were assessed by measuring the Lucia luciferase bioluminescent activity in the supernatant using QUANTI-Luc™. Activity normalized on response upon incubation with  $10^3$  U/ml hIFNβ is shown (A). The NF-κB activity in A549-Dual™-derived cells was assessed by measuring the SEAP activity in the supernatant using QUANTI-Blue™ Solution. Reading of optical density (OD) at 630 nm is shown (B).

### 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](mailto:info@invivogen.com)



Any questions about our cell lines?  
 Visit our FAQ page.

