

Validation data for HEK-Blue™ hACE2-TMPRSS2 cells

<https://www.invivogen.com/hek-blue-hace2-tmprss2-cells>

For research use only

Version 21F10-ED

HEK-Blue™ hACE2-TMPRSS2 cells were generated from the HEK-Blue™ hACE2 cells, which derive from the human embryonic kidney 293 (HEK-293) cell line. Overexpression of the SARS-CoV-2 host receptors ACE2 and TMPRSS2 in HEK-Blue™ hACE2-TMPRSS2 cells has been verified by RT-qPCR (Figure 1) and flow cytometry (Figure 2). In comparison to their parental cell line, these cells display increased sensitivity to pseudotyped lentiviral particles expressing the SARS-CoV-2 Spike protein (Figure 3). Additionally, since HEK-Blue™ hACE2-TMPRSS2 cells express an NF-κB-dependent SEAP reporter, they can be used as 'acceptor' cells in InvivoGen's Spike-ACE2-dependent cell fusion assay (Figure 4).

Validation of ACE2 and TMPRSS2 overexpression by RT-qPCR and flow cytometry

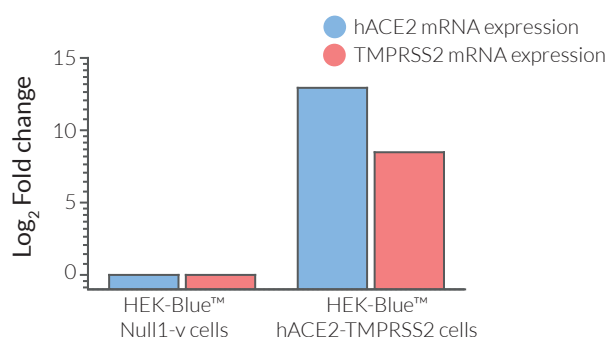


Figure 1: hACE2 and TMPRSS2 mRNA expression. Total mRNA was extracted from ~5x10⁵ HEK-Blue™ Null1-v and HEK-Blue™ hACE2-TMPRSS2 cells. ACE2 and TMPRSS2 mRNA were amplified using quantitative RT-qPCR. Data are represented as the log₂ fold change comparing hACE2 or TMPRSS2 relative expression between the cell lines.

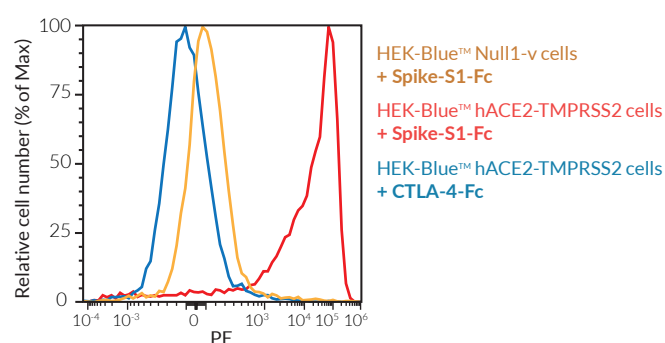


Figure 2: Surface expression of hACE2. ~5x10⁵ HEK-Blue™ Null1-v and HEK-Blue™ hACE2-TMPRSS2 cells were incubated with 1 μg of Spike-S1-Fc or CTLA-4-Fc fusion proteins for 1 hr at 4°C. Cells were then washed and incubated with 0.5 μg of a goat anti-IgG1-Fc antibody coupled to PE for 1 hr at 4°C. Cell surface staining was analyzed by flow cytometry.

Infection by SARS-CoV-2 Spike pseudotyped lentiviral particles

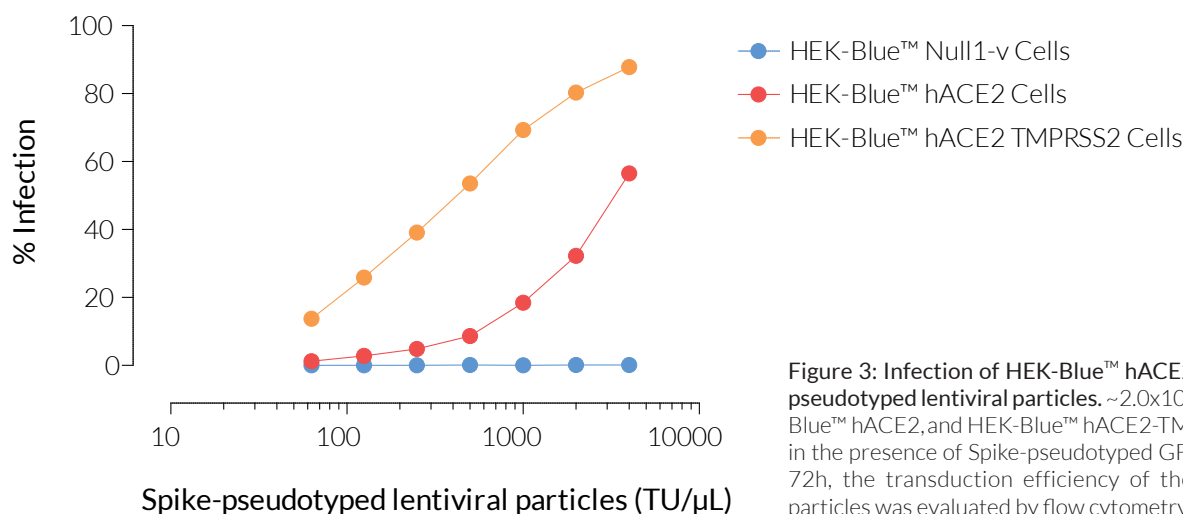


Figure 3: Infection of HEK-Blue™ hACE2-TMPRSS2 cells by Spike pseudotyped lentiviral particles. ~2.0x10⁴ HEK-Blue™ Null1-v, HEK-Blue™ hACE2, and HEK-Blue™ hACE2-TMPRSS2 cells were cultured in the presence of Spike-pseudotyped GFP lentiviral particles. After 72h, the transduction efficiency of the Spike pseudotyped GFP particles was evaluated by flow cytometry.

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



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

 **InvivoGen**
www.invivogen.com

Assessing cell fusion with InvivoGen's COVID-19 cell lines

To generate the 'donor cell line', 293-hMyD88 cells were transiently transfected with InvivoGen's pUNO1-Spike expression plasmid, which encodes the Wuhan-Hu-1 Spike with a functional furin cleavage site to facilitate cell fusion. Upon co-culture of a dilution series of the 'donor cell line' with HEK-Blue™ hACE2-TMPRSS2 cells ('acceptor' cells), fusion was triggered. Following this, MyD88 activated a signalling cascade in the HEK-Blue™ hACE2-TMPRSS2 cells, ultimately leading to readily assessable SEAP production.

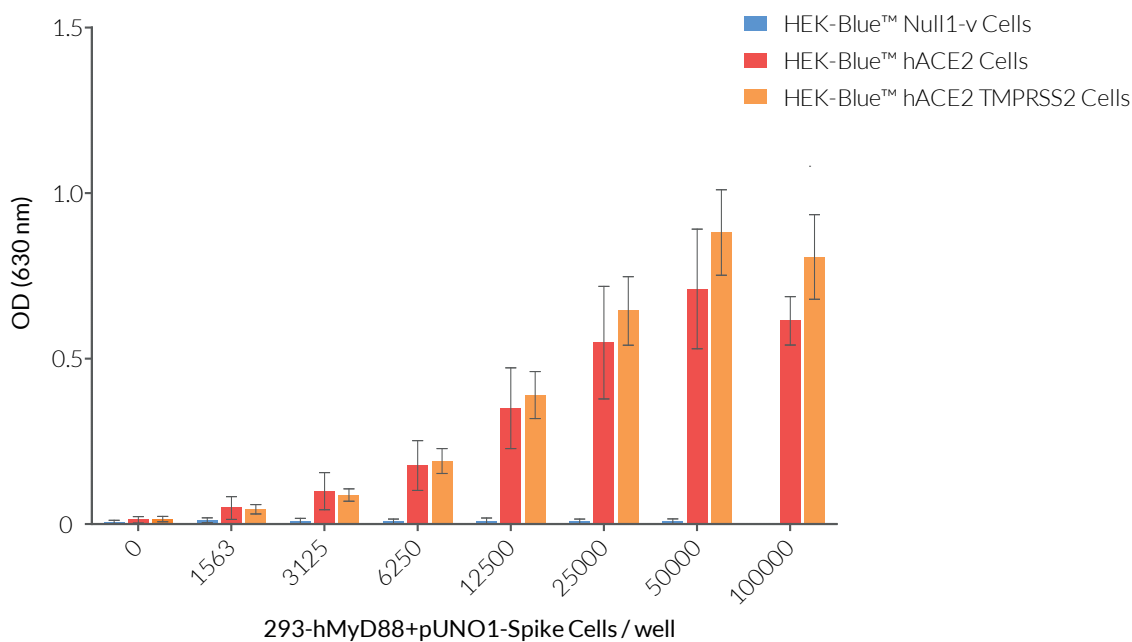


Figure 4: Assessing cell fusion with 293-hMyD88 cells. 293-hMyD88 cells were transiently transfected with a SARS-CoV-2 Spike expression plasmid (pUNO1-Spike) using LyoVec™. After 24 hours, the cells were washed, and a dilution series of the 'donor' 293-hMyD88-Spike cells was co-cultured with either 2.0×10^4 HEK-Blue™ Null1-v, HEK-Blue™ hACE2, or HEK-Blue™ hACE2-TMPRSS2 cells. After overnight incubation, cell fusion was assessed by measuring the activity of SEAP in the supernatant using QUANTI-Blue™ Solution, a SEAP detection reagent. Data are presented as OD_{630nm} ± 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



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

