

# Validation data for 293-SARS2-S-dfur Cells

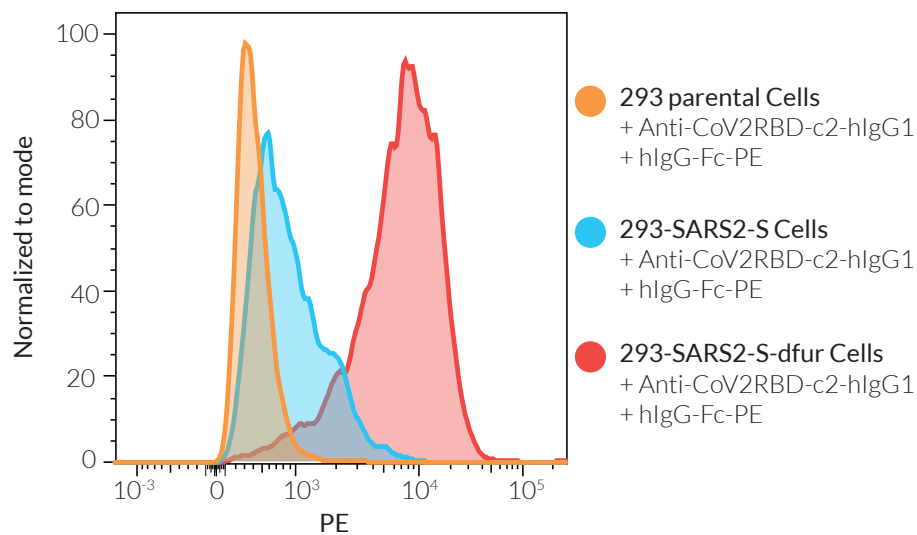
<https://www.invivogen.com/293-sars2-spike>

For research use only

Version 21E03-ED

293-SARS2-S-dfur cells were generated from the human embryonic kidney (HEK)-293 cell line, by stable transfection of InvivoGen's pUNO1-Spike-dfur. This plasmid encodes the original Wuhan-Hu-1 SARS-CoV-2 Spike (D614) protein with an inactivated furin cleavage site (dfur) facilitating improved surface expression for detection by flow cytometry. Therefore, there is strong binding of the Anti-Spike mAb, Anti-CoV2RBD-c2-hIgG1, to 293-SARS2-S-dfur cells when compared to 293-SARS2-S (with a functional furin site) (Figure 1).

## Detection of Spike-binding antibodies by flow cytometry



**Figure 1: Detection of SARS-CoV-2 Spike-binding antibodies by flow cytometry.** 293XL/null (orange), 293-SARS2-S (blue), and 293-SARS2-S-dfur (red) cells were incubated with Anti-CoV2RBD-c2-hIgG1 (clone B38) for 1h at 4°C. Cells were then washed and incubated with goat anti-human hIgG-FC coupled to PE for 1h at 4°C. Cell surface staining was analyzed 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 3-622-34-80  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)



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