

# Validation data for Anti-hACE2-mIgG2b

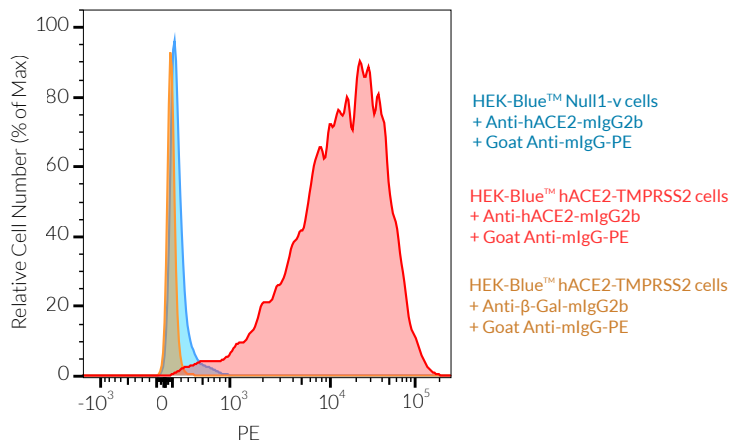
<https://www.invivogen.com/mabg-hace2>

For research use only

Version 21G20-NJ

Anti-hACE2-mIgG2b is a mouse monoclonal antibody (mAb) that targets the human ACE2 (angiotensin-I-converting enzyme-2) surface transmembrane protein. The cell surface staining capacity of Anti-hACE2-mIgG2b has been confirmed with HEK-Blue™ hACE2-TMPRSS2 cells stably overexpressing human ACE2 and TMPRSS2 (transmembrane protease serine 2) (Figure 1).

## Validation of Anti-hACE2-mIgG2b binding by FACS



**Figure 1:** Cell surface staining of human ACE2 using Anti-hACE2-mIgG2b mAb.

~5x10<sup>5</sup> HEK-Blue™ Null1-v (parental) or HEK-Blue™ hACE2-TMPRSS2 cells were incubated with 2 µg of Anti-hACE2-mIgG2b mAb or isotype control for 1h at 4°C. Cells were then washed and incubated with 0.25 µg of goat anti-mIgG 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 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)