

# Validation data for Anti-Spike-RBD mIgG2a (clone CR3022)

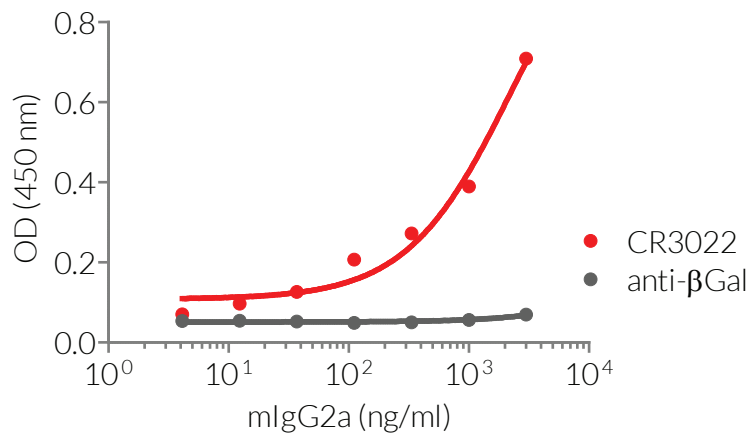
<https://www.invivogen.com/sars2-spike-cr3022-mab-isotypes>

For research use only

Version 20G03-NJ

Anti-Spike-RBD mIgG2a is a recombinant monoclonal antibody (mAb) featuring the variable region of the CR3022 human mAb that specifically targets the SARS-CoV and SARS-CoV-2 Spike receptor-binding domain (RBD) and a mouse IgG2a constant region. The binding of the antibody has been validated by ELISA, using coated Spike-RBD-His fusion peptide and an Anti-mIgG2-HRP secondary antibody. Anti-βGal mIgG2a has been used as a negative control (**Figure 1**).

Binding of Anti-Spike-RBD mIgG2a to SARS-CoV-2 RBD



**Figure 1:** SARS-CoV-2 Spike-RBD-His fusion peptide (5 µg/ml) was coated on ELISA plates overnight. A 3-fold serial dilution of the Anti-Spike-RBD mIgG2a (red curve) or of the Anti-βGal mIgG2a control antibody (grey curve) was performed for the capture step. A HRP-labelled anti-mIgG2a antibody (1/1000 dilution) and the HRP substrate OPD (o-phenylenediamine dihydrochloride) were used for the detection step. Absorbance was read at 490 nm.

## 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)