

# Validation data for Anti-CoV2RBD-c2-hIgG1 (clone B38)

<https://www.invivogen.com/sars2-spike-b38-mab>

For research use only

Version 21B01-ED

Anti-CoV2RBD-c2-hIgG1 is a recombinant monoclonal antibody (mAb) featuring the variable region of the SARS-CoV-2 spike mAb (clone B38), which specifically targets the SARS-CoV-2 Spike receptor-binding domain (RBD), and the constant region of the human IgG1 (hIgG1) isotype. The binding of the antibody has been validated by ELISA using a coated Spike-RBD-His fusion peptide (Figure 1). Infection of InvivoGen's HEK-Blue™ hACE2 cells by SARS-CoV-2 Spike-(G614-variant)-pseudotyped lentiviral particles was successfully neutralized by Anti-CoV2RBD-c2-hIgG1 (clone B38) (Figure 2).

## Binding of Anti-CoV2RBD-c2-hIgG1 to SARS-CoV-2 spike RBD by ELISA

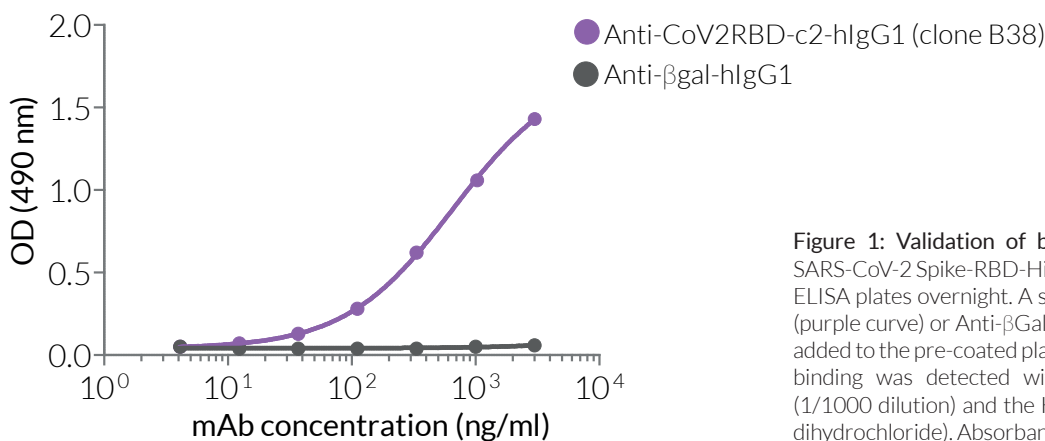


Figure 1: Validation of binding to SARS-CoV-2 RBD by ELISA. SARS-CoV-2 Spike-RBD-His fusion peptide (5 µg/ml) was coated onto ELISA plates overnight. A serial dilution of Anti-CoV2RBD-c2-hIgG1 (purple curve) or Anti-βGal-hIgG1 (control antibody; grey curve) was added to the pre-coated plate and incubated for 1 hour. Subsequently, binding was detected with a HRP-labelled anti-hIgG1 antibody (1/1000 dilution) and the HRP substrate OPD (o-phenylenediamine dihydrochloride). Absorbance was read at 490 nm.

## Anti-CoV2RBD-c2-hIgG1 inhibits infection of HEK293 cells by Spike-pseudotyped lentiviral particles

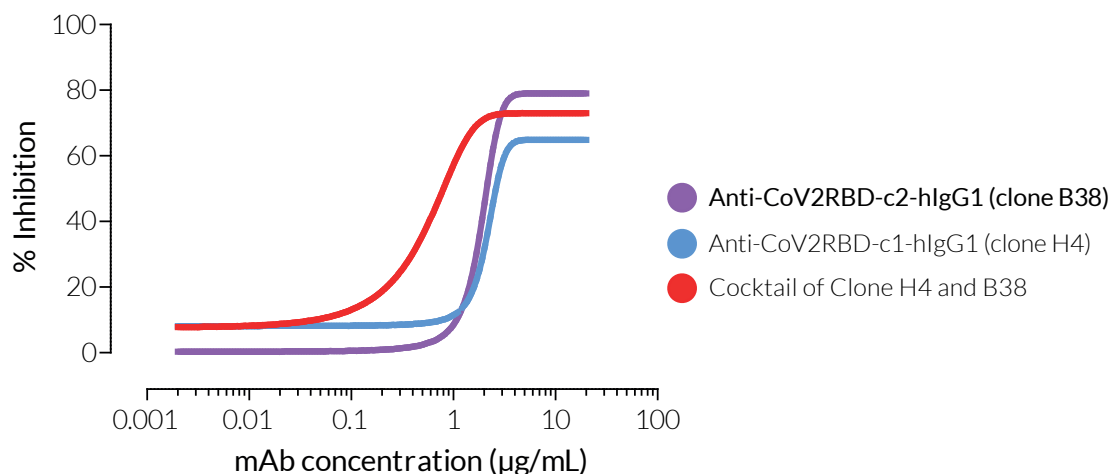


Figure 2: Inhibition of infection by SARS-CoV-2 spike (G614-variant) pseudotyped lentiviral particles. Anti-CoV2RBD-c2-hIgG1 (clone B38), Anti-CoV2RBD-c1-hIgG1 (clone H4), or a cocktail of both mAbs was incubated with SARS-CoV-2 spike-(G614-variant)-pseudotyped lentiviral particles for 1 hour at 37°C. Following this, HEK-Blue™ hACE2 cells were added and incubated for a further 72 hours. Infection by the pseudotyped particles (GFP-expression) was then measured using flow cytometry. Data are presented as % inhibition compared to the antibody-negative control.

### 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)