



6042 Cornerstone Court West, Suite B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### Anti-Spike S1 Antibody (SARS-CoV-2)

SARS-CoV-2, Recombinant, Monoclonal  
**Catalog #:** 100715  
**Lot #:** 200415      **Conc:** 0.10 mg/ml

**Host Species | Isotype:**

Human | Human IgG1

**Formulated in:** PBS and 0.02% sodium azide

**Stability:** At least 6 months at -80°C. *Avoid freeze/thaw cycles.*

**References:**

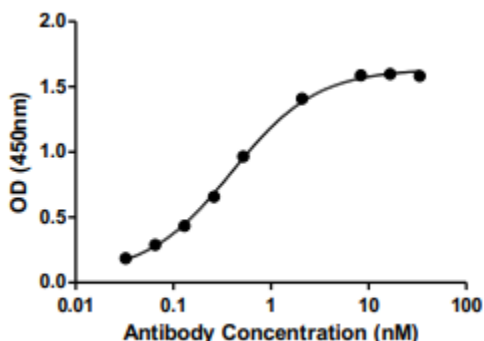
1. Zhou P., *et al.*, *Nature*. 2020; **579**: 270-289.
2. Xiao X., *et al.*, *Cell Mol Life Sci*. 2004; **61(19-20)**: 2428-2430.

**Description:** Recombinant human monoclonal antibody recognizing the SARS-CoV-2 Spike RBD glycoprotein. This antibody cross-reacts with the Spike protein from the SARS-CoV virus.

**Purification:** Purified using protein A resin.

**Applications:** Useful in ELISA (0.1 ug/ml) and FACS (1 ug/ml) assays for detecting Spike RBD. This antibody recognizes a conformation epitope and is not suitable for denaturing Western blotting applications.

### Quality Assurance



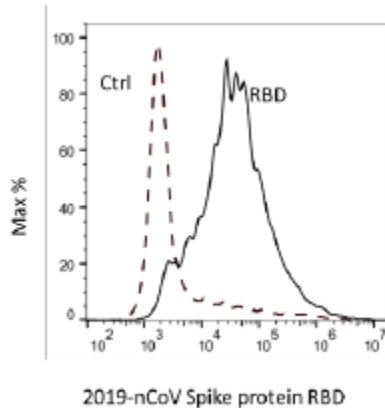
Binding of Anti-Spike Antibody (SARS-CoV-2) to Spike S1 RBD, His-tag (SARS-CoV-2) (BPS Bioscience #100687)

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

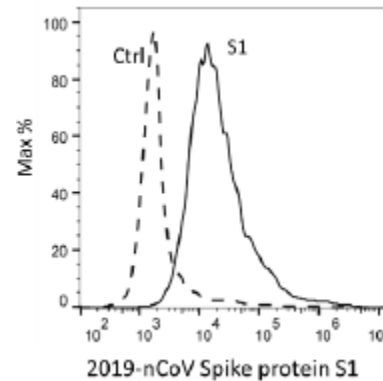
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

## FACS Assay

**Figure A**



**Figure B**



FACS assay showed Spike S1 RBD, Fc fusion (BPS Bioscience #100699) and Spike S1 RBD, His-tag (SARS-CoV-2) (BPS Bioscience #100687) can bind to ACE2 overexpressing Cells (Figure A and B, respectively). ACE2 overexpressing cells were stained with Spike S1 RBD, Fc fusion or Spike S1 RBD, followed by Anti-Spike Antibody (SARS-CoV-2) and fluorescence-conjugated secondary antibody.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)