



6405 Mira Mesa Blvd, Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Data sheet
Spike S1 RBD (SARS-CoV-2):
ACE2 Inhibitor Screening Assay Kit
Catalog #79931
Size: 96 reactions

DESCRIPTION: The pandemic coronavirus disease 2019 (COVID-19) is caused by Severe Acute Respiratory Syndrome Coronavirus 2 (SARS-CoV-2). As a first step of the viral replication strategy, the virus attaches to the host cell surface before entering the cell. The Spike protein receptor binding domain (RBD) recognizes and attaches to the Angiotensin-Converting Enzyme 2 (ACE2) receptor found on the surface of type I and II pneumocytes, endothelial cells, and ciliated bronchial epithelial cells. Drugs targeting the interaction between the Spike protein of SARS-CoV-2 and ACE2 may offer some protection against the viral infection.

The Spike S1 RBD (SARS-CoV-2): ACE2 Inhibitor Screening Assay Kit is designed for screening and profiling inhibitors of this interaction. This kit comes in a convenient 96-well format, with His-labeled ACE2, purified SARS-CoV-2 Spike S1 protein, HRP-labeled anti-His antibody, and assay buffer for 100 binding reactions. The key to this kit is the high sensitivity of detection of bound His-labeled ACE2 by HRP-labeled Anti-His. Only a few simple steps on a microtiter plate are required for the assay. First, SARS-CoV-2 Spike S1 is coated on a 96-well plate. Next, ACE2-His is incubated with SARS-CoV-2 Spike S1 on the plate. Finally, the plate is treated with Anti-His-HRP followed by addition of an HRP substrate to produce chemiluminescence, which then can be measured using a chemiluminescence reader.

COMPONENTS:

Catalog #	Component	Amount	Storage	
100684	SARS-CoV-2 Spike S1 Protein (RBD), mFc Tag	5 µg	-80°C	Avoid multiple freeze/thaw cycles!
11003	ACE2, His-Tag	5 µg	-80°C	
79311	3x Immuno Buffer 1	50 ml	-20°C	
79728	Blocking Buffer 2	50 ml	+4°C	
	Anti-His-HRP	10 µl	-80°C	
79670	ELISA ECL substrate A (transparent bottle)	6 ml	Room Temp	
	ELISA ECL substrate B (brown bottle)	6 ml	Room Temp	
79699	96-well white microplate	1	+4°C	

Note: As of October 1, 2020 this kit has been re-optimized. The old version can still be purchased upon special request.

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: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd, Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

APPLICATIONS: This kit is useful for screenings inhibitors of SARS-CoV-2 Spike binding to ACE2.

STABILITY: Up to 6 months from date of receipt, when stored as recommended.

REFERENCES:

Hoffmann, M. *et al.* 2020. *Cell*, **181**:1-10

Yan, R. *et al.* *Science*, In Press 4 March 2020. DOI: 10.1126/science.abb2762

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

PBS (Phosphate buffered saline)

Luminometer or microplate reader capable of reading chemiluminescence

Adjustable micropipettor and sterile tips

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Coating the plate with SARS-CoV-2 Spike S1:

- 1) Thaw **SARS-CoV-2 Spike S1** on ice. Upon first thaw, briefly spin tube containing **SARS-CoV-2 Spike S1** to recover the full contents of the tube. Aliquot into single use aliquots. Immediately store remaining **SARS-CoV-2 Spike S1** in aliquots at -80°C. Note: **SARS-CoV-2 Spike S1** is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles.
- 2) Dilute **SARS-CoV-2 Spike S1** to 1 µg/ml in PBS.
- 3) Add 50 µl of diluted **SARS-CoV-2 Spike S1** solution to each well and incubate overnight at 4°C.
- 4) Dilute **3x Immuno Buffer 1** to **1x Immuno Buffer 1** with water. Dilute only the amount required for the assay; store remaining 3x Immuno Buffer 1 undiluted.
- 5) Decant to remove supernatant. Wash the plate three times with 100 µl **1x Immuno Buffer 1**. Tap plate onto clean paper towels to remove liquid.
- 6) Block wells by adding 100 µl of **Blocking Buffer 2** to each well. Incubate for 1 hour at room temperature with slow shaking. Remove supernatant as described in step 5.

Step 1:

- 1) Add 20 µl of **1x Immuno Buffer 1** to each well.
- 2) Add 10 µl of inhibitor solution to each well designated "Test Inhibitor". For the "Positive Control" and "Blank", add 10 µl of the same solution without inhibitor (inhibitor buffer). Optionally, incubate at room temperature for one hour with slow shaking.

Note: It is recommendable to use PBS to dilute antibodies or other proteins acting as neutralization inhibitors. When using small molecules dissolved in DMSO, final DMSO

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd, Suite 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

concentration in the assay should be $\leq 1\%$. Inhibitor buffer should contain the same concentration of DMSO as the test inhibitor.

- 3) Thaw **ACE2-His** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Aliquot **ACE2-His** into single use aliquots. Immediately store remaining undiluted enzyme in aliquots at -80°C . Note: **ACE2-His** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.
- 4) Dilute **ACE2-His** to 2.5 ng/ μl (approximately 30 nM) in **1x Immuno Buffer 1**. Keep diluted protein on ice until use. Discard any unused diluted protein after use.
- 5) Add another 20 μl of **1x Immuno Buffer 1** to the wells designated "Blank".

	Blank	Positive Control	Test Inhibitor
1x Immuno Buffer 1	40 μl	20 μl	20 μl
Test Inhibitor	-	-	10 μl
5% DMSO in water (Inhibitor buffer)	10 μl	10 μl	-
ACE2-His (2.5 ng/ μl)	-	20 μl	20 μl
Total	50 μl	50 μl	50 μl

- 6) Initiate reaction by adding 20 μl of diluted **ACE2-His** (see Step 1-4) to wells labeled "Positive Control" and "Test Inhibitor". Incubate at room temperature for one hour with slow shaking.
- 7) Decant to remove supernatant. Wash the plate 3 times with 100 μl /well **1x Immuno Buffer 1**. Tap plate onto clean paper towels to remove liquid.
- 8) Block wells by adding 100 μl of **Blocking Buffer 2** to each well. Incubate for 10 minutes at room temperature. Remove supernatant as in Step 1-7.

Step 2:

- 1) Dilute **Anti-His-HRP** 1000-fold with **Blocking Buffer 2**.
- 2) Add 100 μl to each well. Incubate for 1 hour at room temperature with slow shaking.
- 3) Wash plate three times with **1x Immuno Buffer 1**. Tap plate onto clean paper towel to remove liquid.
- 4) Block wells by adding 100 μl of **Blocking Buffer 2** to each well. Incubate for 10 minutes at room temperature. Decant to remove supernatant. Tap plate onto clean paper towels to remove liquid.
- 5) Just before use, mix 50 μl **ELISA ECL Substrate A** and 50 μl **ELISA ECL Substrate B**, then add 100 μl to each well. Discard any unused chemiluminescent reagent after use.

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd, Suite 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

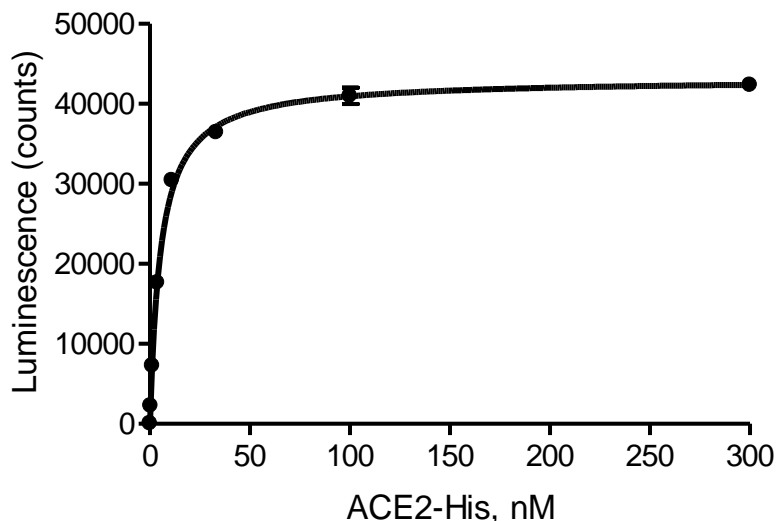
- 6) Immediately read sample in a luminometer or microtiter-plate capable of reading chemiluminescence. “Blank” value is subtracted from all readings.

Reading Chemiluminescence:

Chemiluminescence is the emission of light (luminescence) which results from a chemical reaction. The detection of chemiluminescence requires no wavelength selection because the method used is emission photometry and is not emission spectrophotometry.

To properly read chemiluminescence, make sure the plate reader is set for LUMINESCENCE mode. Typical integration time is 1 second; delay after plate movement is 100 milliseconds. Do not use a filter when measuring light emission. Typical settings for the Synergy 2 BioTek plate reader are: use the “hole” position on the filter wheel; Optics position: Top; Read type: endpoint. Sensitivity may be adjusted based on the luminescence of a control assay without enzyme (typically we set this value as 100).

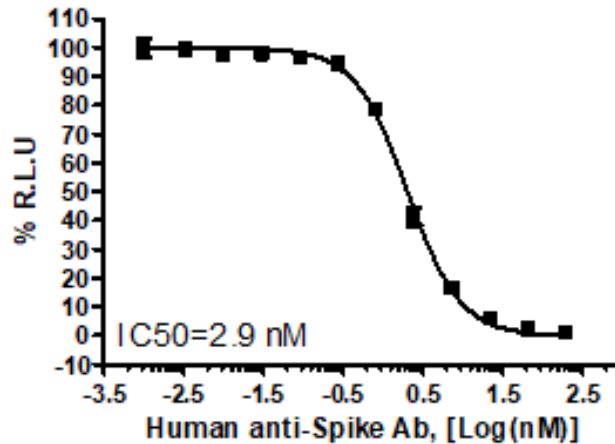
Example of assay results:



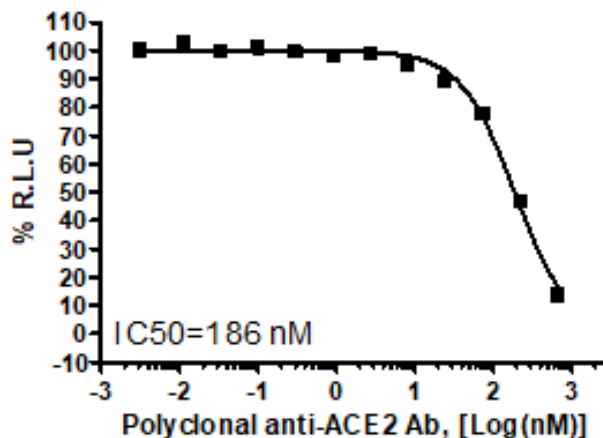
ACE2-His (BPS Bioscience #11003) binding to immobilized SARS-CoV-2 Spike S1 (RBD) (BPS Bioscience #100684) using the *Spike S1 RBD (SARS-CoV-2): ACE2 Inhibitor Screening Assay Kit (BPS Bioscience #79931)*. Luminescence was measured using a BioTek microplate reader. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at support@bpsbioscience.com.

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: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



Human anti-Spike Ab (BPS Bioscience, #100793) at the indicated concentrations was added to the SARS-CoV-2 Spike S1 (RBD) (BPS Bioscience #100684) coated plate and incubated for 1 hour at room temperature. Then, ACE2-His (BPS Bioscience #11003) was added to the wells and followed the Spike S1 RBD (SARS-CoV-2): ACE2 Inhibitor Screening Assay Kit (BPS Bioscience #79931) protocol. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at support@bpsbioscience.com.



ACE2-His (BPS Bioscience #11003) was incubated with various concentrations of the polyclonal anti-ACE2 Ab (R&D Systems, AF933) for 1 hour at room temperature and added to the SARS-CoV-2 Spike S1 (RBD) (BPS Bioscience #100684) coated plate. Then followed the Spike S1 RBD (SARS-CoV-2): ACE2 Inhibitor Screening Assay Kit (BPS Bioscience #79931) protocol. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at support@bpsbioscience.com.

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: support@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd, Suite 100
San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog#</u>	<u>Size</u>
ACE2-His	71158	100 µg
ACE2 Inhibitor Screening Assay Kit	79923	96 reactions
Immuno Buffer 1	79311	50 ml
Blocking Buffer 2	79728	50 ml
ELISA ECL Substrate	79760-1	200 ml

TROUBLESHOOTING GUIDE

Problem	Possible cause	Solution
Luminescence signal of positive control reaction is weak	SARS-CoV-2 Spike ACE2-His has lost activity	Proteins lose activity upon repeated freeze/thaw cycles. Use fresh ACE2-His (BPS Bioscience #11003) and fresh SARS-CoV-2 Spike (BPS Bioscience #TBD). Store proteins in single-use aliquots. Increase time of enzyme incubation. Increase enzyme concentration.
	Incorrect settings on instruments	Refer to instrument instructions for settings to increase sensitivity of light detection.
	Chemiluminescent reagents mixed too soon	Chemiluminescent solution should be used within 15 minutes of mixing. Ensure both reagents are properly mixed.
Luminescent signal is erratic or varies widely among wells	Inaccurate pipetting/technique	Run duplicates of all reactions. Use a multichannel pipettor. Use master mixes to minimize errors.
	Bubbles in wells	Pipette slowly to avoid bubble formation. Tap plate lightly to disperse bubbles; be careful not to splash between wells.
Background (signal to noise ratio) is high	Insufficient washes	Increase number of washes. Increase wash volume.
	Sample solvent is inhibiting the enzyme	Run negative control assay including solvent. Maintain DMSO level at <1% Increase time of enzyme incubation.
	Results are outside the linear range of the assay	Use different concentrations of ACE2-His (BPS Bioscience #11003) to create a standard curve

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com