

Description

The Spike S1 RBD (B.1.1.7 Variant) (N501Y) (SARS-CoV-2):ACE2 Inhibitor Screening Chemiluminescence Assay Kit is designed for screening and profiling inhibitors of the interaction of ACE2 with the B.1.1.7 Variant (N501Y) of the SARS-CoV-2 Spike S1 RBD protein. This kit comes in a convenient 96-well format, with purified SARS-CoV-2 Spike S1 RBD B.1.1.7 Variant (N501Y) and ACE2-Biotin proteins, streptavidin-HRP, chemiluminescence HRP substrate, and assay buffer for 100 binding reactions. The key to this kit is the high sensitivity of detection of ACE2-Biotin protein by Streptavidin-HRP. Only a few simple steps on a microtiter plate are required for the assay. First, Spike S1 RBD protein B.1.1.7 Variant is coated on a 96-well transparent plate. Next, ACE2-Biotin is incubated with Spike S1 RBD B.1.1.7 Variant on the plate. Finally, the plate is treated with streptavidin-HRP followed by addition of an HRP substrate to produce chemiluminescence, which can then be measured using a luminometer or microplate reader capable of reading chemiluminescence.

Background

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 Receptor Binding Domain (RBD) of Spike protein 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. It has been widely suggested that active as well as passive immunizations targeting the interaction between the Spike protein of SARS-CoV-2 and ACE2 offer promising protection against the viral infection. However recent reports showed that a mutant strain first identified in the UK (B1.1.7) exhibits higher transmissibility and infectivity. Investigations on the effects of mutations on viral replication and pathogenesis will be critical for developing effective strategies for vaccines and antibody therapies against COVID-19.

Applications

This kit is useful for screening for inhibitors of ACE2 binding to SARS-CoV-2 Spike S1 RBD B.1.1.7 Variant (N501Y).

Supplied Materials

Catalog #	Name	Amount	Storage
	Spike S1 RBD (N501Y), His-tag (SARS-CoV-2)	5 µg	-80°C
100665	ACE2, His-Avi-Tag, Biotin-labeled HiP™	5 µg	-80°C
79742	Streptavidin-HRP	15 µl	+4°C
79311	3x Immuno Buffer 1	50 ml	-20°C
79728	Blocking Buffer 2	50 ml	+4°C
79670	ELISA ECL Substrate A (translucent bottle)	6 ml	Room Temp
79670	ELISA ECL Substrate B (brown bottle)	6 ml	Room Temp
79699	White 96-well microplate	1	Room Temp

Materials Required but Not Supplied

PBS (Phosphate buffered saline)
Rotating or rocker platform
Microplate reader capable of reading chemiluminescence

Storage Conditions



Upon arrival, store individual kit components at the recommended temperatures. Avoid freeze/thaw cycles of frozen components.

Stability

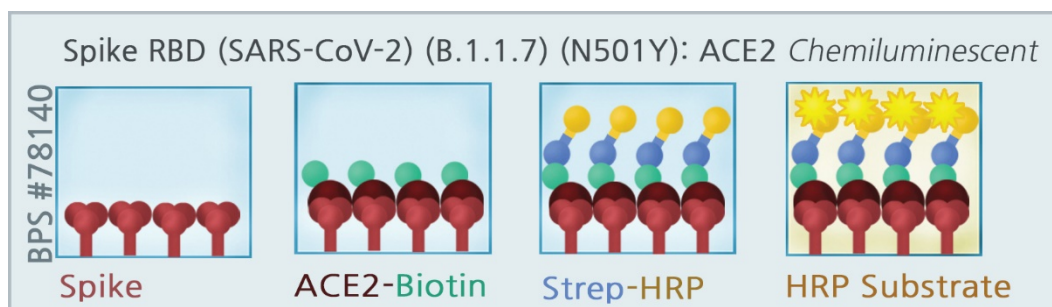
This assay kit will perform optimally for up to 6 months from date of receipt when the materials are stored as directed.

Safety



This product is for research purposes only and not for human or therapeutic use. This product should be considered hazardous and is harmful by inhalation, in contact with skin, eyes, clothing, and if swallowed. If contact occurs, wash thoroughly.

Assay Principle



Assay Protocol

All samples and controls should be tested in duplicate. We recommend preincubating antibodies or protein inhibitors with the target protein. For small molecule inhibitors, pre-incubation may also be beneficial, depending on the experimental conditions.


Coating the plate with Spike S1 RBD B.1.1.7 Variant (N501Y)

1. Thaw Spike S1 RBD B.1.1.7 Variant (N501Y) on ice. Upon first thaw, briefly spin tube containing Spike S1 RBD B.1.1.7 Variant (N501Y) to recover the full contents of the tube. Aliquot into single use aliquots. Immediately store remaining Spike S1 RBD B.1.1.7 Variant (N501Y) in aliquots at -80°C . Note: Spike S1 RBD B.1.1.7 Variant (N501Y) is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles.
2. Dilute Spike RBD to $1\ \mu\text{g}/\text{ml}$ in PBS.
3. Add $50\ \mu\text{l}$ of diluted Spike S1 RBD B.1.1.7 Variant (N501Y) solution to each well of the microtiter plate 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 ($\sim 60\ \text{ml}$ of 1x Immuno Buffer is needed for a 96-well plate).
5. Decant to remove supernatant. Wash the plate three times with $100\ \mu\text{l}$ 1x Immuno Buffer 1. Tap plate onto clean paper towels to remove liquid.
6. Block wells by adding $100\ \mu\text{l}$ of Blocking Buffer 2 to each well. Incubate for 1 hour at room temperature with slow shaking. Remove supernatant as described in step 5.

Component	Blank	Positive Control	Test Sample
Neutralizing antibody (or Protein inhibitor) in Blocking Buffer 2	-	-	25 μ l
Blocking Buffer 2	25 μ l	25 μ l	-
Biotin-labeled ACE2 (1-2 ng/ μ l)	-	25 μ l	25 μ l
Blocking Buffer 2	25 μ l	-	-
Total	50 μ l	50 μ l	50 μ l

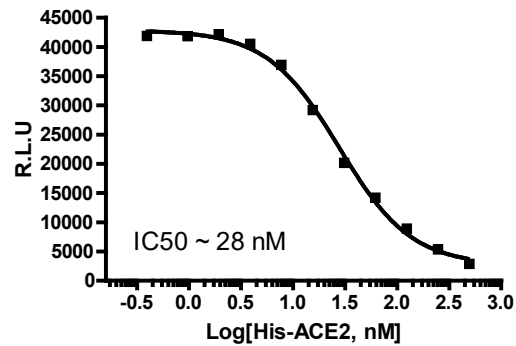
For the SARS-CoV2 Spike S1 RBD B.1.1.7 Variant (N501Y) neutralizing antibody (or protein inhibitor)

1. Prepare serial dilutions of the test neutralizing antibody or protein inhibitor in Blocking Buffer 2.
2. Add 25 μ l of the serially diluted neutralizing antibody or protein inhibitor to each well designated "Test Sample." For the wells designated "Blank" and "Positive Control," add 25 μ l of Blocking Buffer 2.
3. Incubate the plate for 30 minutes at room temperature (neutralizing antibody or protein inhibitor can be preincubated up to 60 minutes).
4. Thaw the biotin-labeled ACE2 on ice. After thawing, briefly spin the tube containing biotin-labeled ACE2 to recover the full contents of the tube. Aliquot into single use aliquots. Immediately store remaining biotin-labeled ACE2 in aliquots at -80°C. Note: Biotin labeled ACE2 is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles.
5. Dilute biotin-labeled ACE2 to 1.5 ng/ μ l in Blocking Buffer 2.
6. Add 25 μ l of the diluted biotin-labeled ACE2 to the wells designated "Test Sample" and "Positive Control"
7. For the wells designated "Blank", add 25 μ l Blocking Buffer 2.
8. Incubate the plate for 1 hour at room temperature with slow shaking.
9. After 1 hour, decant the solution and wash the plate three times with 100 μ l 1x Immuno Buffer 1. Tap plate onto clean paper towels to remove liquid.
10. Dilute Streptavidin-HRP 1000-fold with Blocking Buffer 2.
11. Add 100 μ l to each well. Incubate for 30 minutes at room temperature with slow shaking.
12. After 30 minutes, decant the solution and wash plate three times with 1x Immuno Buffer 1. Tap plate onto clean paper towel to remove liquid.
13. Just before use, mix 50 μ l ELISA ECL Substrate A and 50 μ l ELISA ECL Substrate B, then add 100 μ l to each well. Discard any unused chemiluminescent reagent after use.
14. 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 Results



His-ACE2 protein (BPS Bioscience, #11003) competes and blocks the binding of biotin-labeled ACE2 to SARS-CoV-2 Spike S1 RBD B.1.1.7 Variant (N501Y). Experiment was performed using the Spike S1 RBD (B.1.1.7 Variant) (N501Y) (SARS-CoV-2) :ACE2 Inhibitor Screening Chemiluminescence Assay Kit, BPS Bioscience, #78140. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at support@bpsbioscience.com

General considerations

“Blank” Control: The “Blank” control is important to determine the background absorbance in the assay.

Visit bpsbioscience.com/assay-kits-faq for detailed troubleshooting instructions. For all further questions, please email support@bpsbioscience.com

References

1. Wang P. *et al.*, Increased Resistance of SARS-CoV-2 Variants B.1.351 and B.1.1.7 to Antibody Neutralization. *bioRxiv* 2021 Jan 26; 2021.01.25.428137
2. Shen X., *et al.*, SARS-CoV-2 Variant B.1.1.7 is susceptible to neutralizing antibodies elicited by ancestral Spike vaccines. *bioRxiv*. 2021 Jan 29; 2021.01.27.428516
3. Hoffman M. *et al.*, SARS-CoV-2 Cell Entry Depends on ACE2 and TMPRSS2 and Is Blocked by a Clinically Proven Protease Inhibitor. *Cell* 2020; 181:1-10

Related Products

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
Spike S1 RBD (B.1.1.7 Variant) (N501Y) (SARS-CoV-2): ACE2 Inhibitor Screening Colorimetric Assay Kit	78133	96 reactions
Spike S1 RBD (SARS-CoV-2): ACE2 Inhibitor Screening Assay Kit	79931	96 reactions
ACE2 : Spike S1 RBD (SARS-CoV-2) Inhibitor Screening Assay Kit	79936	96 reactions
ACE2 : Spike S1 RBD (SARS-CoV-2) Inhibitor Screening Colorimetric Assay Kit	78031	96 reactions
SARS-CoV-1 Spike Trimer (S1+S2) : ACE2 Inhibitor Screening Colorimetric Assay Kit	78012	96 reactions
ACE2, His-Avi-Tag	11003-1	20 µg
ACE2, His-Avi-Tag, Biotin-labeled HiP™	100665-1	20 µg
Spike (B.1.1.7 Variant) (SARS-CoV-2) Pseudotyped Lentivirus (Luc Reporter)	78112-2	500 µl x 2
Spike (SARS-CoV-2) Pseudotyped Lentivirus (Luc Reporter)	79942-2	500 µl x 2
ACE2: Spike S1 RBD (SARS-CoV-2) Inhibitor Screening Assay Kit	79936	96 reactions