



6042 Cornerstone Court W, Ste 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**  
***ACE2 Inhibitor Screening Assay Kit***  
**Catalog #79923**  
**Size: 96 reactions**

**BACKGROUND:** Angiotensin converting enzyme 2 (ACE2) is an exopeptidase that catalyzes the conversion of angiotensin II to angiotensin 1-7 and L-phenylalanine. Angiotensin II is part of the classical renin angiotensin system (RAS), a hormone system that regulates fluid balance, blood pressure and maintains vascular tone. ACE2 has been also proved to be the receptor for the human respiratory coronavirus NL63, the SARS-coronavirus (SARS-CoV) and the novel coronavirus 2019-nCoV/SARS-CoV-2.

**DESCRIPTION:** The *ACE2 Inhibitor Screening Assay Kit* is designed to measure the exopeptidase activity of ACE2 for screening and profiling applications. The ACE2 assay kit comes in a convenient 96-well format, with purified ACE2, its substrate, and ACE2 buffer for 96 reactions.

**COMPONENTS:**

Catalog #	Component	Amount	Storage	
11003	ACE2	2 µg	-80 °C	Avoid multiple freeze/thaw cycles!
	ACE2 Fluorogenic Substrate	2.5 ml	-80 °C	
	ACE2 Buffer	3 ml	-20 °C	
79685	96-well black microplate	1	Room Temp	

**MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:**

Adjustable micropipettor and sterile tips  
Fluorescent microplate reader

**APPLICATIONS:** Useful for studying enzyme kinetics and screening small molecule inhibitors and antibodies for drug discovery and HTS applications.

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

**REFERENCES:**

Danilczyk, U., *et al. Circulation Research*. 2006. **98(4)**: 463-471.  
Ge, X. Y., *et al. Nature*. 2013. **503(7477)**: 535-538.

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)



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

#### ASSAY PROTOCOL:

*All samples and controls should be tested in duplicate.*

1. Thaw **ACE2** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Note: **ACE2** is sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.
2. Prepare **Enzyme solution** (0.5 ng/μl ACE2) by diluting ACE2 in **ACE2 Buffer**.
3. Add 20 μl of **Enzyme solution** (0.5 ng/μl ACE2) to each well designated "Positive Control" and "Test Inhibitor," and 20 μl of **ACE2 buffer** to each well designated "Blank."
4. Add 5 μl of **Test Inhibitor** solution to each well designated "Test Inhibitor." For the wells labeled "Positive Control" and "Blank," add 5 μl of 10% DMSO in water (Inhibitor buffer). Note: Keep the final concentration of DMSO in the reaction ≤1%.
5. Add 25 μl of **ACE2 Fluorogenic Substrate** to all wells. Protect from light and incubate reaction at room temperature for 60 minutes.

	Positive Control	Test Inhibitor	Blank
Enzyme solution (0.5 ng/μl ACE2)	20 μl	20 μl	-
ACE2 Buffer	-	-	20 μl
Test inhibitor	-	5 μl	-
10% DMSO in water (Inhibitor buffer)	5 μl	-	5 μl
ACE2 Fluorogenic Substrate	25 μl	25 μl	25 μl
<b>Total</b>	<b>50 μl</b>	<b>50 μl</b>	<b>50 μl</b>

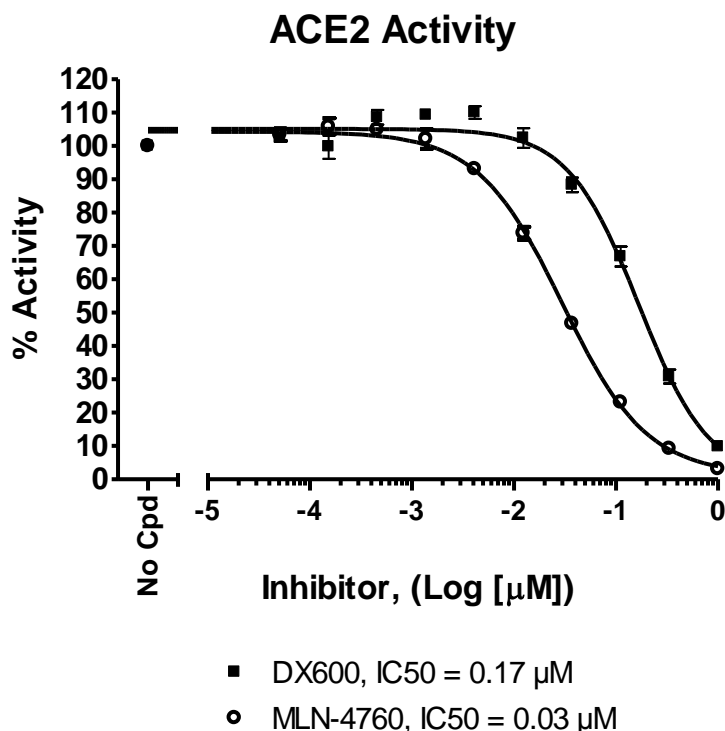
6. Read fluorescence intensity of the samples ( $\lambda_{\text{excitation}} = 555 \text{ nm}$ ;  $\lambda_{\text{emission}} = 585 \text{ nm}$ ) in an appropriate microplate reader.

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)

**Example of assay results:**



ACE2 inhibition by DX600 (Cayman, #22186) and MLN-4760 (Sigma, #5306160001) measured using the *ACE2 Inhibitor Screening Assay Kit*, BPS Bioscience Catalog #79923. Fluorescence was measured using a Bio-Tek microplate reader. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**RELATED PRODUCTS:**

<u>Product</u>	<u>Catalog#</u>	<u>Size</u>
ACE2	11003	10 μg
PrCP, His-tag	80380	10 μg
Renin, Human, His-tag	80200	100 μg
Renin, Rat, His-tag	80201	100 μg
Renin Assay Kit	80211	96 rxns.

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)