



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

Data Sheet

3x ACE2-Spike TR-FRET Buffer

Catalog #: 79953

DESCRIPTION: The *3x ACE2-Spike TR-FRET Assay Buffer* is optimized for use with the Spike S1-Biotin (SARS-CoV-2): ACE2 TR-FRET Assay Kit (BPS Bioscience, #79949)

APPLICATIONS: Studying enzyme kinetics and ACE2 applications in drug discovery.

QUANTITY: 4 ml

STORAGE: -20°C. Avoid freeze/thaw cycles.

STABILITY: Stable for at least one year from date of receipt when stored as directed.

INSTRUCTIONS FOR USE: For a detailed protocol on the use of this product, please see the protocol for BPS Bioscience, #79949.

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
ACE2, His-Tag, Eu-labeled	100705	100 µg
ACE2, His-Avi-Tag, Biotin-labeled HiP™	100665	20 µg or 50 µg
ACE2, His-Tag	11003	20 µg or 100 µg
Spike S1, Fc fusion, Avi-tag, Biotin-Labeled (SARS-CoV-2)	100679	20 µg or 50 µg

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

Please visit our website at: www.bpsbioscience.com