



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

Data Sheet

3CL Protease Assay Buffer

Catalog #: 79956

Description: The *3CL Protease Assay Buffer* is optimized for use with the 3CL Protease (SARS-CoV-2) Assay Kit (BPS Bioscience #79955) and the 3CL Protease (SARS-CoV) Assay Kit (BPS Bioscience #78015).

Applications: Study enzyme kinetics, and screen small molecular inhibitors of 3CL SARS-CoV-2 protease for HTS applications in drug discovery.

Supplied As: Liquid

Format: Aqueous buffer solution

Storage/Stability: Store at -20°C. Avoid freeze/thaw cycles. Stable at least 6 months from date of receipt, when stored as directed.

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)