



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

Data Sheet
Papain-like Protease Assay Buffer
Catalog #: 78039

Description: The *Papain-like Protease (PLPro) Assay Buffer* is optimized for use with the Papain-like Protease (SARS-CoV-2): Protease Activity Assay Kit (BPS Bioscience #79995).

Applications: Study enzyme kinetics, and screen small molecular inhibitors of Papain-like Protease (PLPro) (SARS-CoV-2) 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.

Instructions for Use: Add 0.5 M DTT to PLPro Assay Buffer so final DTT concentration is 1 mM. For example, add 10 µl of 0.5 M DTT to 5 ml assay buffer. (DTT should be added just before use. Prepare only enough DTT-containing buffer as required for the assay. Store the remaining assay buffer at -20°C).

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