



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

3CL Protease Fluorogenic Standard

Catalog #: 78006

Description: Fluorogenic Standard for the determination of 3CL protease (Main protease, Mpro) Activity. Corresponds to the Nsp5 cleavage sequence of SARS-CoV and SARS-Cov-2.

Application: Useful as a fluorogenic standard for the 3CL Protease Assay Kit (BPS Bioscience, #79955)

Sequence: H-SGFRKME-EDANS-NH₂

Counter Ion: Trifluoroacetate

Theoretical Mass: 1101.3

Observed Mass: 1100.8

Purity: ≥95%

Concentration: 1 mg at 5 mM

Supplied as: DMSO solution

Storage: Stable at +4°C for up to 6 months. For longer storage, store in single use aliquots at -20°C. Avoid freeze/thaw cycles. Protect from light and moisture. Aqueous solutions are stable for one day.

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