

Data Sheet

3CL Protease, MBP-tagged (SARS-CoV-2) Assay Kit

Catalog #79955-1
Size: 96 reactions

Datasheet updated December 2021

BACKGROUND: Coronaviruses (CoVs) cause respiratory and intestinal infections in humans and animals. The 3CL protease, also known as Main Protease (Mpro), plays a vital role in processing the polyproteins that are translated from the viral RNA. Protease inhibitors that can block viral replication are promising potential drug candidates for the treatment of patients suffering from COVID-19 infection.

DESCRIPTION: The *3CL Protease Assay Kit* is designed to measure 3CL Protease activity for screening and profiling applications, in a homogeneous assay with no time-consuming washing steps. The kit comes in a convenient 96-well format, with purified 3CL Protease MBP-tag (BPS Bioscience, #100707), fluorogenic substrate, and 3CL Protease assay buffer for 100 enzyme reactions. 3CL inhibitor GC376 is also included as control.

The 3CL Protease Substrate is an internally quenched 14-mer fluorogenic peptide (DABCYL-KTSAVLQSGFRKME-EDANS). When the donor (EDANS) and acceptor (DABCYL) fluorophores are in close proximity, the energy emitted from EDANS is quenched by DABCYL (intact substrate). Upon proteolysis by 3CL, the peptide substrate is cleaved between glutamine and serine by the 3CL protease to generate the highly fluorescent peptide fragment (SGFRKME-EDANS). The fluorescence intensity increases proportionally to the activity of 3CL. More information on the substrate, including MW and structure, can be found on our website (BPS Bioscience, #79952).

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

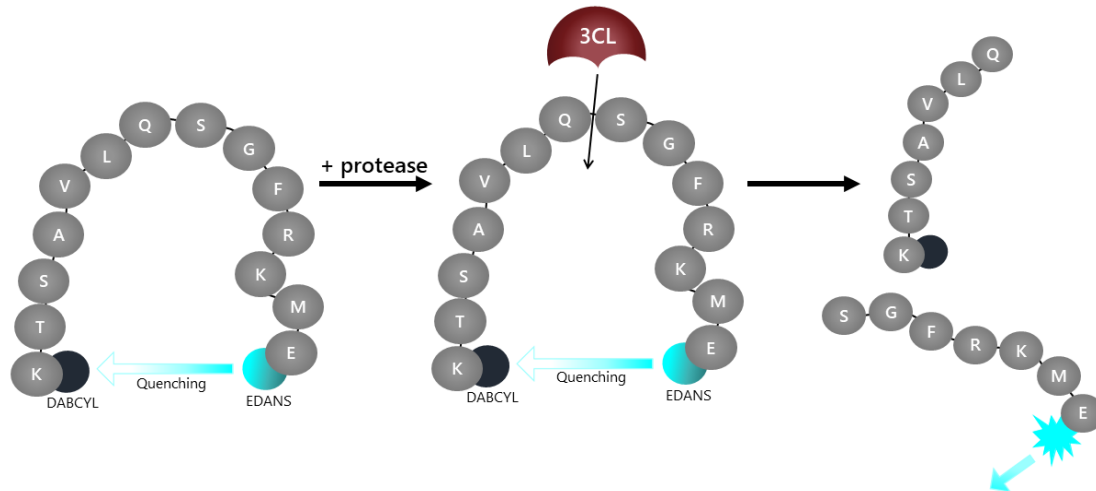


Figure 1: Illustration of the principle behind the 3CL protease assay. The 3CL Protease Substrate is an internally quenched 14-mer fluorogenic peptide (DABCYL-KTSAVLQSGFRKME-EDANS). When the donor (EDANS) and acceptor (DABCYL) fluorophores are in close proximity the energy emitted from EDANS is quenched by DABCYL (intact substrate). Upon proteolysis by 3CL, the peptide substrate is cleaved between the glutamine and serine residues to generate the highly fluorescent peptide fragment (SGFRKME-EDANS). The fluorescence intensity increases proportionally to the activity of 3CL. More information on the substrate, including MW and structure, can be found on our website (BPS Bioscience #79952).

COMPONENTS:

Catalog #	Component	Amount	Storage	
100707	Recombinant 3CL Protease, MBP-tag*	20 µg	-80°C	Avoid freeze/thaw cycles!
79952	3CL Protease Substrate (10 mM)	25 µl	-80°C	
79956	3CL Protease Assay Buffer	25 ml	-20°C	
78013	GC376 (MW=507.5 Dalton)	50 µg	-20°C	
	0.5 M DTT	200 µl	-20°C	
79685	Black, low binding microtiter plate	1	Room Temperature	
	Plate sealing film	1		

**The exact concentration of protein is lot-specific and will be indicated on the tube containing the protein.*

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

3CL inhibitor GC376 is provided as a technical control for 3CL inhibition. More information on GC376, including MW and molecular structure, can be found on our website (BPS Bioscience, #78013).

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Fluorescent microplate reader capable of reading exc/em=360 nm/460 nm

APPLICATIONS: Great for studying enzyme kinetics and for high throughput (HTS) applications.

STABILITY: At least one year from date of receipt when stored as directed.

REFERENCE(S):

1. Jared S. Morse, *et al.*, 2020 *Chem.Bio.Chem.* **21**:730 – 738.
2. Zhang, L., *et al.* 2020, *Science* **368 (6489)**: 409-412.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

- 1) Just before use, dilute **0.5 M DTT** 500 times into the **3CL Protease Assay Buffer** to obtain a DTT concentration of 1 mM. For example, add 10 µl of **0.5 M DTT** to 5 ml of assay buffer. Prepare enough DTT-containing buffer as required for the assay. Store the remaining assay buffer at -20°C.
- 2) Thaw the **3CL Protease** on ice. Briefly spin the tube containing the enzyme to recover the full content of the tube. The kit provides enough enzyme for one plate. Note: **3CL Protease** is sensitive to freeze/thaw cycles. Do not re-use the diluted enzyme.
- 3) Dilute **3CL Protease** in the **Assay buffer** (containing 1 mM DTT) at 3-5 ng/µl (90-150 ng per reaction).

Note: The exact concentration and volume of enzyme is lot-specific and will be indicated on the tube. Calculate required dilution from the information in the tube. It may be desirable to dilute the enzyme serially to avoid using large amounts of protease assay buffer for the dilution.

- 4) Add 30 µl of **diluted 3CL Protease** to the wells designated as “Positive Control”, “Inhibitor Control” (i.e. GC376) and “Test Inhibitor”. Add 30 µl of **Assay buffer** (containing 1 mM DTT) to the “Blank” wells.
- 5) Dilute the 50 µg **GC376** in 200 µl of Assay Buffer to obtain a 500 µM solution. Add 10 µl of **GC376** (500 µM) to the wells labeled “Inhibitor Control”. Aliquot and store the remaining solution at -80°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

Note: adding 10 μ l of GC376 (500 μ M solution) to the assay in a final volume of 50 μ l results in a final concentration of 100 μ M.

6) Prepare the Test Inhibitor.

The final concentration of DMSO in the assay should not exceed 1%. If the Test Inhibitor is dissolved in DMSO, make a 100-fold higher concentration of the compound than the highest desired concentration. Then make a 20-fold dilution in 1x **Assay Buffer** (containing DTT). At this step the compound concentration is 5-fold higher than the final concentration.

If the test inhibitor is dissolved in water, make a 5-fold higher concentration of the test compound than the final desired concentration in the **Assay buffer** (containing DTT).

Add 10 μ l of Test Inhibitor to each well designated “Test Inhibitor”.

7) Diluent Solution (no inhibitor): add 10 μ l of assay buffer (if the test compound is water soluble and was diluted in assay buffer) or 10 μ l of 5% DMSO diluted in assay buffer (if DMSO was used to dissolve the test inhibitor) to “Blank” and “Positive Control” wells.

Component	Positive Control	Test Inhibitor	Inhibitor Control	Blank
3CL Protease (3-5 ng/ μ l)	30 μ l	30 μ l	30 μ l	–
Assay Buffer (with DTT)	–	–	–	30 μ l
GC376 (500 μ M)	–	–	10 μ l	–
Test Inhibitor	–	10 μ l	–	–
Diluent Solution (no inhibitor)	10 μ l	–	–	10 μ l
Total	40 μl	40 μl	40 μl	40 μl

8) Preincubate for 30 min at room temperature with slow shaking.

9) Dilute the 25 μ l of **3CL Protease substrate** (10mM) in 1 ml of the **Assay Buffer** containing DTT, to make a 250 μ M solution. The final concentration of the **3CL Protease substrate** in a 50 μ l reaction is 50 μ M.

10) Start the reaction by adding 10 μ l of the substrate solution to all the wells, including the “Blank”. This brings the final reaction volume to 50 μ l.

Seal the plate with the plate sealer and incubate overnight at room temperature.

11) Measure the fluorescence intensity in a microtiter plate-reading fluorimeter capable of excitation at a wavelength 360 nm and detection of emission at a wavelength

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**

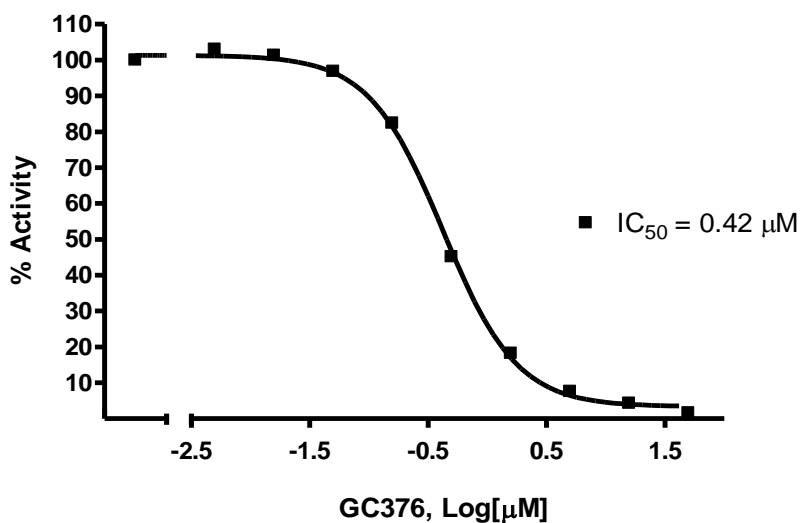
Or you can Email us at: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

460 nm. The fluorescence intensity can also be measured kinetically. The “Blank” value is subtracted from all other values.

EXAMPLE OF ASSAY RESULTS:

3CL Protease SARS-CoV-2 Activity



Inhibition of 3CL Protease enzyme activity by increasing concentrations of GC376, measured using the *Fluorogenic 3CL Protease Assay Kit* (BPS Bioscience #79955). Fluorescence intensity was measured using a Tecan fluorescent microplate reader. Results are expressed as percent of activity relative to the no-inhibitor control (set at 100%).

Data shown is representative. For lot-specific information, please contact BPS Bioscience, Inc. at support@bpsbioscience.com

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



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

RELATED PRODUCTS

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
3CL Protease, Untagged (SARS-CoV-2) Assay Kit	78042	96 reactions
3CL Protease, Untagged (SARS-CoV-2) Assay Kit	78042	384 reactions
3CL Protease (MERS-CoV) Inhibitor Screening Assay Kit	78278	96 reactions
3CL Protease (SARS-CoV-1) Assay Kit	78015	96 reactions
3CL Protease (Mpro), MBP-tag (SARS-CoV-2)	100707-1	100 µg
3CL Protease (Mpro), MBP-tag (SARS-CoV-2)	100707-2	1 mg
PLPro, His-tag (SARS-CoV-2)	100735	20 µg/50 µg
PLPro, His-tag (SARS-CoV)	81091	25 µg
SARS-CoV-2 Spike:ACE2 Inhibitor Screening Kit	79931	96 reactions
ACE2:SARS-CoV-2 Spike Inhibitor Screening Kit	79936	96 reactions
ACE2:SARS-CoV-2 Spike S1-Biotin Inhibitor Screening Kit	79945	96 reactions
SARS-CoV-2 Spike S1-Biotin:ACE2 TR-FRET Kit	79949	96 reactions

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