

Amplite™ Universal Fluorimetric Protease Activity Assay Kit

Green Fluorescence

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 13500 (500 assays)	Store at -20 °C and keep from light Component C can be stored at 4 °C	Fluorescence microplate readers

Introduction

Protease assays are widely used for the investigation of protease inhibitors and detection of protease activities. Monitoring various protease activities has become a routine task for many biological laboratories. Some proteases have been identified as good drug development targets.

Our Amplite™ Universal Fluorimetric Protease Activity Assay Kit is an ideal choice to perform routine assays for the isolation of proteases, or for identifying the presence of contaminating proteases in protein samples. The kit uses a fluorescent casein conjugate which is proven to be a generic substrate for a broad spectrum of proteases (e.g. trypsin, chymotrypsin, thermolysin, proteinase K, protease XIV, and elastase). In the intact substrate, casein is heavily labeled with a green fluorescent dye, resulting in significant fluorescence quenching. Protease-catalyzed hydrolysis relieves its quenching effect, yielding brightly fluorescent dye-labeled short peptides. The increase in fluorescence intensity is directly proportional to protease activity. The assay can be performed in a convenient 96-well or 384-well microtiter plate format and readily adapted to automation. Its signal can be easily read with a fluorescence microplate reader at Ex/Em = 490/525 nm using FITC filter set.

Kit Key Features

Convenient Format :	Include all the key assay components.
Optimized Performance:	Optimal conditions for the detection of generic protease activity.
Continuous:	Easily adapted to automation without a separation step.
Convenient:	Formulated to have minimal hands-on time. No wash is required.
Non-Radioactive:	No special requirements for waste treatment.

Kit Components

Components	Amount
Component A: Protease Substrate	1 vial (300 µL), protected from light
Component B: Trypsin	1 vial (100 µL, 5 U/µL)
Component C: 2X Assay Buffer	1 bottle (30 mL)

Assay Protocols for One 96-well Plate

Please choose Protocol A or B according to your needs.

Protocol A: Measure protease activity in test samples

Brief Summary

Prepare protease substrate solution (50 µL) → Add substrate control, positive control or test samples (50 µL) → Incubate for 0 min (for kinetic reading) or 30 minutes - 1 hour (for end point reading) → Monitor fluorescence intensity at Ex/Em = 490/525 nm

Note: Thaw all the kit components at room temperature before starting the experiment.

1. Prepare working solutions:

- 1.1 Make protease substrate solution: Dilute Protease Substrate (Component A) at 1:100 in 2X assay buffer (Component C). Use 50 μL of protease substrate solution per assay in a 96-well plate.
*Note: The 2X Assay Buffer (Component C) is designed for detecting the activity of chymotrypsin, trypsin, thermolysin, proteinase K, protease XIV, and human leukocyte elastase. For other proteases, please refer to **Appendix I** for the appropriate assay buffer formula.*
- 1.2 Trypsin dilution: Dilute Trypsin (5 U/ μL , Component B) at 1:50 in de-ionized water to get a concentration of 0.1 U/ μL .

2. Add reagents prepared in step 1 into a 96-well microplate according to Table 1 and Table 2.**Table 1.** Layout of the substrate control, positive control, and test samples in a 96-well microplate

SC	SC										
PC	PC										
TS	TS										
....										

Note: SC=Substrate Control, PC =Positive Control, TS=Test Samples.

Table 2. Reagent composition for each well

Substrate Control	Positive Control	Test Sample
De-ionized water: 50 μL	Trypsin dilution: 50 μL	Protease-containing samples: 50 μL
Total volume: 50 μL	Total volume: 50 μL	Total volume: 50 μL

Note: If less than 50 μL of protease-containing biological sample is used, add ddH₂O to make a total volume of 50 μL .

3. Run the enzymatic reaction:

- 3.1 Add 50 μL of protease substrate solution (from Step 1.1) to all the wells in the assay plate. Mix the reagents well.
- 3.2 Monitor the fluorescence increase with a fluorescence plate reader at Ex/Em = 490/525 nm.
For kinetic reading: Immediately start measuring fluorescence intensity continuously and record data every 5 minutes for 30 minutes.
For end-point reading: Incubate the reaction at a desired temperature for 30 to 60 minutes, protected from light. Then measure the fluorescence intensity.

4. Data analysis:

Refer to **Data Analysis** section.

Protocol B: Screening protease inhibitors using a purified enzyme**Brief Summary**

Prepare protease substrate solution (10 μL) → Add substrate control, positive control, vehicle control or test samples (90 μL) → Incubate for 0 minute (for kinetic reading) or 30 minutes - 1 hour (for end point reading) → Monitor fluorescence intensity at Ex/Em = 490/525 nm

1. Prepare working solutions:

- 1.1 Make 1X assay buffer: Add 5 mL de-ionized water into 5 mL of 2X Assay Buffer (Component C).
- 1.2 Make protease substrate solution: Dilute Protease Substrate (Component A) at 1: 20 in 1X assay buffer (from Step 1.1). Use 10 μL /well of protease substrate solution for a 96-well plate.

- A variety of data analyses can be done, e.g., determining inhibition %, IC_{50} , K_m , K_i , etc.

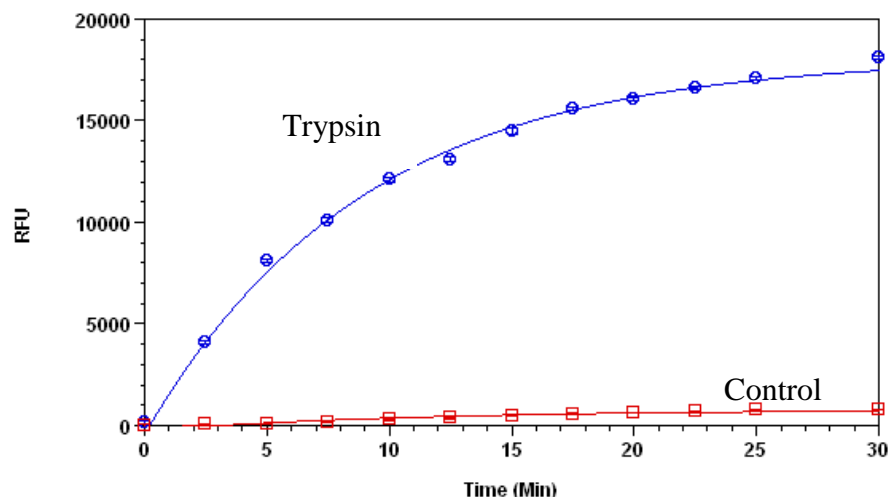


Figure 1. Trypsin protease activity was analyzed by using Amplitude™ Universal Fluorimetric kit. Protease substrate was incubated with 1 unit trypsin in the kit assay buffer. The control wells had protease substrate only (without trypsin). The fluorescence signal was measured starting from time 0 when trypsin was added using a NOVOSTar microplate reader (BMG Labtech) with a filter set of Ex/Em = 490/525 nm. Samples were done in triplicates.

Appendix I

Protease	1X Assay Buffer*
Cathepsin D	20 mM Sodium Citrate, pH 3.0
Papain	20 mM sodium acetate, 20 mM cysteine, 2 mM EDTA, pH 6.5
PAE	20 mM sodium phosphate, pH 8.0
Pepsin	10 mM HCl, pH 2.0
Porcine pancreas elastase	10 mM Tris-HCl, pH 8.8
Subtilisin	20 mM potassium phosphate buffer, pH 7.6, 150 mM NaCl

Note: * For protocol A, 2X assay buffer is needed. For protocol B, 1X assay buffer is needed.

References

1. Vineyard D, Zhang X, Lee I. (2006) Transient kinetic experiments demonstrate the existence of a unique catalytic enzyme form in the peptide-stimulated ATPase mechanism of Escherichia coli Lon protease. *Biochemistry*, 45, 11432.
2. Yadav SC, Pande M, Jagannadham MV. (2006) Highly stable glycosylated serine protease from the medicinal plant Euphorbia milii. *Phytochemistry*, 67, 1414.
3. Frohbieter KA, Ismail B, Nielsen SS, Hayes KD. (2005) Effects of Pseudomonas fluorescens M3/6 bacterial protease on plasmin system and plasminogen activation. *J Dairy Sci*, 88, 3392.
4. Chauhan V, Sheikh AM, Chauhan A, Spivack WD, Fenko MD, Malik MN. (2005) Fibrillar amyloid beta-protein inhibits the activity of high molecular weight brain protease and trypsin. *J Alzheimers Dis*, 7, 37.
5. Lee EH, Kim CS, Cho JB, Ahn KJ, Kim KH. (2003) Measurement of protease activity of live Uronema marinum (Ciliata: Scuticociliatida) by fluorescence polarization. *Dis Aquat Organ*, 54, 85.
6. Cilenti L, Lee Y, Hess S, Srinivasula S, Park KM, Junqueira D, Davis H, Bonventre JV, Alnemri ES, Zervos AS. (2003) Characterization of a novel and specific inhibitor for the pro-apoptotic protease Omi/HtrA2. *J Biol Chem*, 278, 11489.

Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of the kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at info@aatbio.com if you have any questions.