

**ab112153**

# **Protease Activity Assay Kit (Fluorometric – Red)**

## Instructions for Use

For detecting Protease activity in biological samples or to screen protease inhibitors using our proprietary red fluorescence probe

[View kit datasheet: www.abcam.com/ab112153](http://www.abcam.com/ab112153)

(use [www.abcam.cn/ab112153](http://www.abcam.cn/ab112153) for China, or [www.abcam.co.jp/ab112153](http://www.abcam.co.jp/ab112153) for Japan)

This product is for research use only and is not intended for diagnostic use.



# Table of Contents

---

|    |   |    |
|----|---|----|
| 1. | Introduction  | 3  |
| 2. | Protocol Summary  | 5  |
| 3. | Kit Contents  | 6  |
| 4. | Storage and Handling                                    | 6  |
| 5. | Assay Protocol  | 7  |
|    | Protocol I: Measurement of Protease Activity in Samples | 7  |
|    | Protocol II: Screening Protease Inhibitors              | 10 |
| 6. | Data Analysis   | 13 |
| 8. | Troubleshooting   | 16 |

# 1. 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.

ab112153 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.

ab112153 uses a red fluorescent casein conjugate that 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 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. Its signal can be easily read at Ex/Em = 540 /590 nm.

ab112153 has been used for screening protease inhibitors in a HTS mode

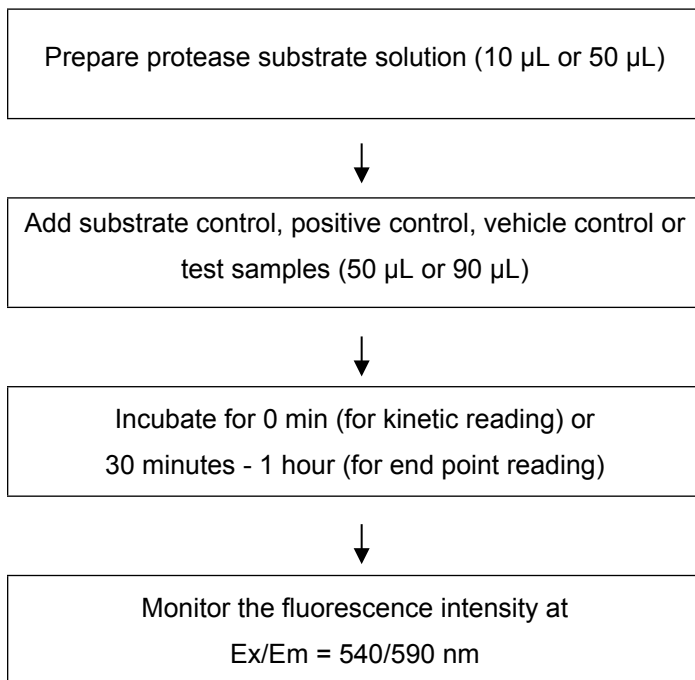
## Kit Key Features

- **Convenient Format:** Includes all the key assay components
- **Optimized Performance:** Optimized 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.

## 2. Protocol Summary

---

*Summary for One 96-well Plate (see each individual protocol for full details)*



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

### 3. Kit Contents

---

---

| <b>Components</b>                                 | <b>Amount</b> |
|---|---------------|
| Component A: Protease Substrate (Light Sensitive) | 300 $\mu$ L   |
| Component B: Trypsin 5U/ $\mu$ L                  | 100 $\mu$ L   |
| Component C: 2X Assay Buffer                      | 30 mL         |

---

### 4. Storage and Handling

---

Keep at -20°C. Avoid exposure to light.

Component C can be stored at 4°C for convenience

## 5. Assay Protocol

---

*Note: This protocol is for one 96 - well plate.*

Please choose Protocol I or II according to your needs.

### **Protocol I: Measurement of Protease Activity in Samples**

#### **A. Preparation of Working Solutions:**

1. Make protease substrate solution: Dilute Protease Substrate (Component A) at 1:100 in 2X assay buffer Component C). Use 50  $\mu\text{L}$ /well of protease substrate solution per assay for 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.*

2. Trypsin dilution: Dilute Trypsin (5 U/ $\mu\text{L}$ , Component B) at 1:50 in de-ionized water to get a concentration of 0.1 U/ $\mu\text{L}$ .



- B.** Add reagents prepared in step 1 into a 96-well microplate according to Table 1 and Table 2:

|      |      |      |      |
|------|------|------|------|
| SC   | SC   | .... | .... |
| PC   | PC   | .... | .... |
| TS   | TS   | .... | .... |
| .... | .... |      |      |
| .... | .... |      |      |
| .... | .... |      |      |
| .... | .... |      |      |
| .... | .... |      |      |

**Table 1.** Layout of the substrate control, positive control, and test samples in a 96-well microplate. *Note: SC=Substrate Control, PC =Positive Control, TS=Test Samples.*

| Identifier               | Contents                    | Volume     |
|--------------------------|-----------------------------|------------|
| <b>Substrate Control</b> | De-ionized water:           | 50 $\mu$ L |
| <b>Positive Control</b>  | Trypsin dilution            | 50 $\mu$ L |
| <b>Test Sample</b>       | Protease-containing samples | 50 $\mu$ L |

**Table 2.** Reagent composition for each well. *Note: If less than 50  $\mu$ L of protease-containing biological sample is used, add ddH<sub>2</sub>O to make a total volume of 50  $\mu$ L.*

- C. Run the Enzyme Assay**

1. Add 50  $\mu\text{L}$  of protease substrate solution (from Step A.1) into each well of the assay plate. Mix the reagents well
2. Monitor the fluorescence increase with a fluorescence plate reader at Ex/Em = 540/590 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.

#### **D. Data Analysis**

Refer to the Data Analysis section.

## **Protocol II: Screening Protease Inhibitors (Purified Enzyme)**

### **A. Preparation of Working Solutions:**

1. Make 1X assay buffer: Add 5 mL de-ionized water into 5 mL of 2X Assay Buffer (Component C).
2. Make protease substrate solution: Dilute Protease Substrate (Component A) at 1: 20 in 1X assay buffer (from Step A.1). Use 10  $\mu$ L/well of protease substrate solution for 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*

3. Protease dilution: Dilute the protease in 1X assay buffer to a concentration of 500-1000 nM (For Trypsin 50-100 U/mL). Each well will need 10  $\mu$ L of protease dilution. Prepare an appropriate amount for all the test samples and extra for the positive control and vehicle control wells

- B.** Add reagents prepared in step 1 into a 96-well microplate according to Table 1 and Table 2:

|      |      |      |      |
|------|------|------|------|
| SC   | SC   | .... | .... |
| PC   | PC   | .... | .... |
| VC   | VC   | .... | .... |
| TS   | TS   |      |      |
| .... | .... |      |      |
| .... | .... |      |      |
| .... | .... |      |      |
| .... | .... |      |      |

**Table 1.** Layout of appropriate controls and test samples in a 96-well microplate.

*Note1: SC=Substrate Control, PC= Positive Control, VC=Vehicle Control, TS=Test Samples.*

*Note 2: It's recommended to test at least three different concentrations of each test compound. All the test samples should be done in duplicates or triplicates.*

| <b>Identifier</b>        | <b>Contents</b>  | <b>Total Volume</b> |
|--------------------------|--|---------------------|
| <b>Substrate Control</b> | 1X Assay Buffer  | 90 $\mu$ L          |
| <b>Positive Control</b>  | 1X assay buffer: 80 $\mu$ L<br>Protease dilution: 10 $\mu$ L                                 | 90 $\mu$ L          |
| <b>Vehicle Control</b>   | Vehicle*: X $\mu$ L<br>1X assay buffer: (80-X) $\mu$ L<br>Protease dilution: 10 $\mu$ L      | 90 $\mu$ L          |
| <b>Test Sample</b>       | Test compound: X $\mu$ L<br>1X assay buffer: (80-X) $\mu$ L<br>Protease dilution: 10 $\mu$ L | 90 $\mu$ L          |

**Table 2.** Reagent composition for each well.

*Note : \*For each volume of test compound added into a well, the same volume of solvent used to deliver test compound needs to be checked for the effect of vehicle on the activity of protease.*

### C. Run the Enzyme Reaction:

1. Add 10  $\mu\text{L}$  of protease substrate solution (from Step A.2) into the wells of positive control (PC), vehicle control (VC), and test sample (TS). Mix the reagents well.
2. Monitor the fluorescence intensity with a fluorescence plate reader at Ex/Em = 540 /590 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

### D. Data Analysis

Refer to the Data Analysis section.

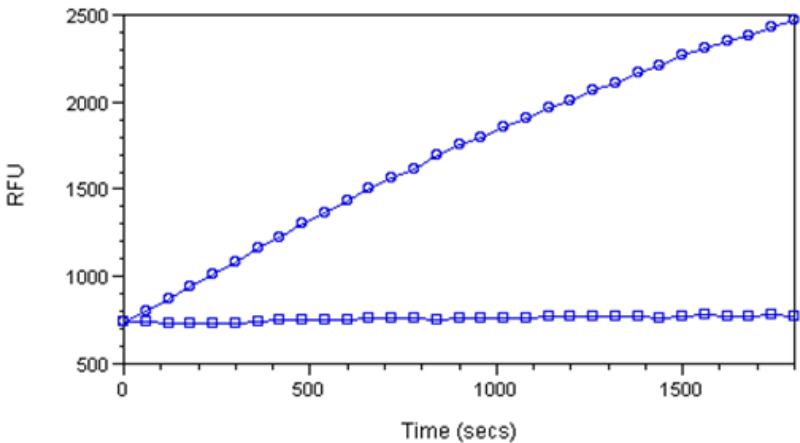
## 6. Data Analysis

---

The fluorescence in the substrate control wells is used as a control, and is subtracted from the values for other wells with the enzymatic reactions.

- Plot data in the format of relative fluorescence unit (RFU) versus time for each sample (as shown in Figure 1).

- Determine the range of initial time points during which the reaction is linear. 10-15% conversion appears to be the optimal range.
- Obtain the initial reaction velocity ( $V_0$ ) in RFU/min. Determine the slope of the linear portion of the data plot.
- A variety of data analyses can be done, e.g., inhibition percentage,  $IC_{50}$ ,  $K_m$ ,  $K_i$ , etc.



**Figure 1.** Trypsin protease activity was analyzed using ab112153. Protease substrate was incubated with 3 units 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 microplate reader with a filter set of Ex/Em = 540/590nm. Samples were done in triplicates.

## 7. Appendix I

---

| <b>Protease</b>           | <b>1X Assay Buffer*</b>                                 |
|---------------------------|---|
| 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 I, 2X assay buffer is needed. For Protocol II, 1X assay buffer is needed.*



## 8. Troubleshooting

| <b>Problem</b>     | <b>Reason</b>                            | <b>Solution</b>   |
|--------------------|--|---|
| Assay not working  | Assay buffer at wrong temperature        | Assay buffer must not be chilled - needs to be at RT  |
|                    | Protocol step missed                     | Re-read and follow the protocol exactly   |
|                    | Plate read at incorrect wavelength       | Ensure you are using appropriate reader and filter settings (refer to datasheet)  |
|                    | Unsuitable microtiter plate for assay    | Fluorescence: Black plates (clear bottoms);<br>Luminescence: White plates;<br>Colorimetry: Clear plates.<br>If critical, datasheet will indicate whether to use flat- or U-shaped wells |
| Unexpected results | Measured at wrong wavelength             | Use appropriate reader and filter settings described in datasheet   |
|                    | Samples contain impeding substances      | Troubleshoot and also consider deproteinizing samples   |
|                    | Unsuitable sample type                   | Use recommended samples types as listed on the datasheet  |
|                    | Sample readings are outside linear range | Concentrate/ dilute samples to be in linear range   |

| <b>Problem</b>                                  | <b>Reason</b>   | <b>Solution</b>   |
|---|---|---|
| Samples with inconsistent readings              | Unsuitable sample type                                | Refer to datasheet for details about incompatible samples   |
|   | Samples prepared in the wrong buffer                  | Use the assay buffer provided (or refer to datasheet for instructions)                                    |
|   | Samples not deproteinized (if indicated on datasheet) | Use the <b>10kDa spin column (ab93349)</b> or <b>Deproteinizing sample preparation kit (ab93299)</b>      |
|   | Cell/ tissue samples not sufficiently homogenized     | Increase sonication time/ number of strokes with the Dounce homogenizer                                   |
|   | Too many freeze-thaw cycles                           | Aliquot samples to reduce the number of freeze-thaw cycles  |
|   | Samples contain impeding substances                   | Troubleshoot and also consider deproteinizing samples   |
|   | Samples are too old or incorrectly stored             | Use freshly made samples and store at recommended temperature until use                                   |
| Lower/ Higher readings in samples and standards | Not fully thawed kit components                       | Wait for components to thaw completely and gently mix prior use   |
|   | Out-of-date kit or incorrectly stored reagents        | Always check expiry date and store kit components as recommended on the datasheet                         |
|   | Reagents sitting for extended periods on ice          | Try to prepare a fresh reaction mix prior to each use   |
|   | Incorrect incubation time/ temperature                | Refer to datasheet for recommended incubation time and/ or temperature                                    |
|   | Incorrect amounts used                                | Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume) |

|                              |  |  |
|------------------------------|--|--|
| Standard curve is not linear | Not fully thawed kit components                        | Wait for components to thaw completely and gently mix prior use  |
|                              | Pipetting errors when setting up the standard curve    | Try not to pipette too small volumes   |
|                              | Incorrect pipetting when preparing the reaction mix    | Always prepare a master mix  |
|                              | Air bubbles in wells                                   | Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates |
|                              | Concentration of standard stock incorrect              | Recheck datasheet for recommended concentrations of standard stocks  |
|                              | Errors in standard curve calculations                  | Refer to datasheet and re-check the calculations   |
|                              | Use of other reagents than those provided with the kit | Use fresh components from the same kit   |

**For further technical questions please do not hesitate to contact us by email ([technical@abcam.com](mailto:technical@abcam.com)) or phone (select “contact us” on [www.abcam.com](http://www.abcam.com) for the phone number for your region).**

**For all technical and commercial enquires  
please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)