

Protease Activity Assay Kit (Fluorometric - Green)
ab112152

2 References 1 Image

Overview

Product name	Protease Activity Assay Kit (Fluorometric - Green)
Detection method	Fluorescent
Sample type	Serum, Plasma, Cell culture extracts, Other biological fluids
Assay type	Enzyme activity
Assay time	1h 00m
Product overview	<p>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.</p> <p>ab112152 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. ab112152 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.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
Platform	Microplate reader

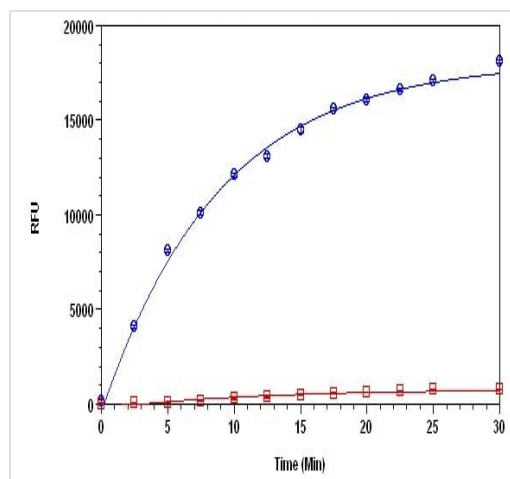
Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	500 tests
2X Assay Buffer	1 x 30ml

Components	500 tests
Protease Substrate	1 x 300µl
Trypsin	1 x 100µl

Images



Functional Studies - Protease Activity Assay Kit
(Fluorometric - Green) (ab112152)

Trypsin protease activity was analyzed by using ab112152. 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 microplate reader with a filter set of Ex/Em = 490/525 nm. Samples were done in triplicates.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors