

Product datasheet

Protease Activity Assay Kit (Fluorometric - Red)
ab112153

1 Image

Overview

Product name	Protease Activity Assay Kit (Fluorometric - Red)
Detection method	Fluorescent
Sample type	Plasma, Cell culture extracts, Other biological fluids, Whole Blood
Assay type	Enzyme activity (quantitative)
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>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.</p> <p>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.</p> <p>ab112153 has been used for screening protease inhibitors in a HTS mode.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
Platform	Microplate reader

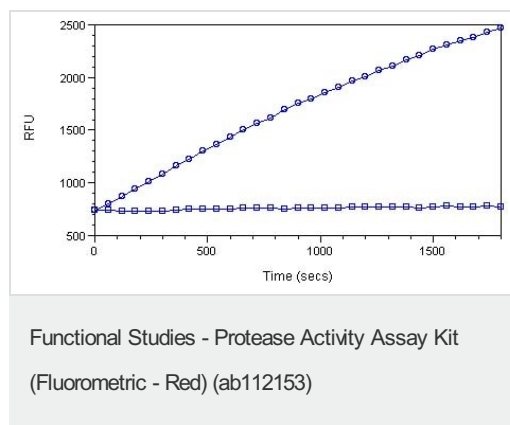
Properties

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

Components	500 tests
Assay Buffer	1 x 30ml

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

Images



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

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