

Protease Activity Assay Kit (Fluorometric - Green) ab112152

画像数 1

製品の概要

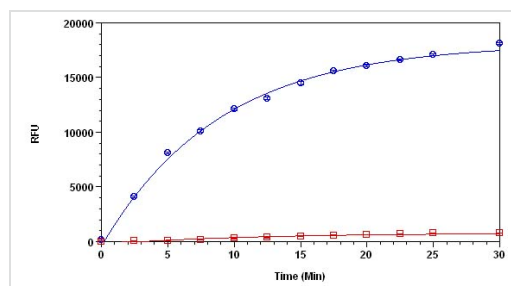
製品名	Protease Activity Assay Kit (Fluorometric - Green)
サンプルの種類	Plasma, Cell culture extracts, Other biological fluids, Whole Blood
アッセイタイプ	Quantitative
全工程の試験時間	1h 00m
製品の概要	<p>Protease assays are widely used for the investigation of protease inhibitors and detection of protease activities. Monitoring 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>

Visit our [FAQs page](#) for tips and troubleshooting.

製品の特性

保存方法 Please refer to protocols.

内容	500 tests
2X Assay Buffer	1 x 30ml
Protease Substrate	1 x 300µl
Trypsin	1 x 100µl



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 AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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