

Protease Activity Assay Kit ab111750

4 References 2 图像

概述

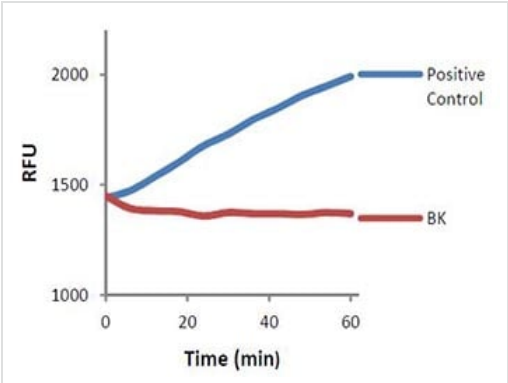
产品名称	Protease Activity Assay试剂盒
检测方法	Fluorescent
样品类型	Cell culture supernatant, Milk, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
检测类型	Enzyme activity
灵敏度	< 500 pg/well
检测时间	1h 00m
产品概述	<p>Abcam's Protease Activity Assay Kit is designed for the quantitative determination of proteases present in the protein sample. The assay uses fluorescein isothiocyanate (FITC)-labeled casein as a general protease substrate. The fluorescein label on the FITC-Casein is highly quenched. Upon digestion by proteases present in the sample the FITC-Casein substrate is cleaved into smaller peptides which abolishes the quenching of the fluorescence label. The fluorescence of the FITC-labeled peptide fragments is measured at Ex/Em = 485/530 nm. The kit is supplied with our Mass Spectrometry Grade (MSG), chemically stabilized Trypsin for use as a general protease control. However, other protease standard controls can also be used. This kit is easy to use and can detect < 500 pg/well Trypsin present in the sample</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K781 Protease Activity Fluorometric Assay Kit. K781-100 is the same size as the 100 test size of ab111750.</p> <p>Proteases are naturally present in all organisms. These enzymes are involved in a multitude of physiological reactions from simple digestion of food proteins to highly regulated cascades. Proteases can either break specific peptide bonds (limited proteolysis), depending on the amino acid sequence of a protein, or break down a complete peptide to amino acids (unlimited proteolysis). The activity can be a destructive change (abolishing a protein's function), an activation of a function (preform to mature form) or it can be a signal in a signaling pathway.</p>
平台	Microplate reader

性能

存放说明 Store at -20°C. Please refer to protocols.

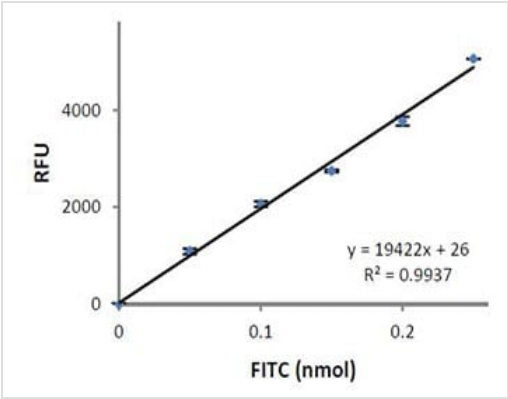
组件	100 tests
Assay Buffer I	1 x 25ml
Fluorescence Standard VI	1 x 200µl
Protease Positive Control	1 vial
Protease Substrate	1 vial

图片



Functional Studies - Protease Activity Assay Kit
(ab111750)

Positive control assayed using ab111750



Functional Studies - Protease Activity Assay Kit
(ab111750)

Example Standard curve obtained using ab111750

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.cn/abpromise> or contact our technical team.

Terms and conditions

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