

Cathepsin H Activity Assay Kit (Fluorometric) ab65305

1 图像

概述

产品名称	Cathepsin H Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Tissue Extracts, Cell Lysate
检测类型	Quantitative
检测时间	2h 00m
产品概述	Cathepsin H Activity Assay Kit (Fluorometric) (ab65305) is a fluorescence-based assay that utilizes the preferred cathepsin-B substrate Arginine labeled with AFC (amino-4-trifluoromethyl coumarin). Cell lysates or other samples that contain cathepsin-B will cleave the synthetic substrate R-AFC to release free AFC. The released AFC can easily be quantified using a fluorometer or fluorescence plate reader. The released AFC can easily be quantified using a fluorometer or fluorescence plate reader. The Cathepsin H assay is simple, straightforward, and can be adapted to 96-well plate assays. Assay conditions have been optimized to obtain the maximal activity. Visit our FAQs page for tips and troubleshooting.

说明	This product is manufactured by BioVision, an Abcam company and was previously called K145 Cathepsin H Activity Fluorometric Assay Kit. K145-100 is the same size as the 100 test size of ab65305. Apoptosis can be mediated by mechanisms other than the traditional caspase-mediated cleavage cascade. There is growing recognition that alternative proteolytic enzymes such as the lysosomal cathepsin proteases may initiate or propagate proapoptotic signals. Cathepsins are lysosomal enzymes that are also used as sensitive markers in various toxicological investigations.
----	---

平台	Microplate reader
----	-------------------

性能

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

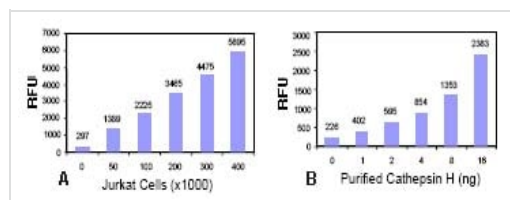
组件	100 tests
CH Cell Lysis Buffer	1 x 25ml
CH Inhibitor (1 mM)	1 x 20µl

组件	100 tests
CH Reaction Buffer	1 x 5ml
CH Substrate R-AFC (10 mM)	1 x 200µl

相关性 The cathepsin family of proteolytic enzymes contains several diverse classes of proteases. The cysteine protease class comprises cathepsins B, L, H, K, S, and O. The aspartyl protease class is composed of cathepsins D and E. Cathepsin G is in the serine protease class. Most cathepsins are lysosomal and each is involved in various cellular events such as peptide biosynthesis and protein degradation. The propeptide region of cathepsin H is highly homologous to other cathepsins. This region is thought to play a role in the processing of cysteine proteases. Cathepsin H is thought to be involved in several pathological conditions, including neuro degenerative disorders and cancer.

细胞定位 Lysosome

图片



Cathepsin H assays were performed using various numbers of Jurkat cells (A) or various amounts of purified human liver cathepsin H (B), as indicated. Results were analyzed using a fluorescence plate reader (Ex/Em = 400/505 nm) as described in the kit instructions.

Functional Studies - Cathepsin H Activity
Fluorometric Assay Kit (ab65305)

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