

Cathepsin K Activity Assay Kit (Fluorometric) ab65303

[7 References](#) [2 图像](#)

概述

产品名称	Cathepsin K Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Tissue Extracts, Cell Lysate
检测类型	Enzyme activity
检测时间	2h 00m
产品概述	<p>Cathepsin K Activity Assay Kit (Fluorometric) (ab65303) is a fluorescence-based assay that utilizes the preferred cathepsin K substrate sequence LR labeled with AFC (amino-4-trifluoromethyl coumarin). Cell lysates or other samples that contain cathepsin K will cleave the synthetic substrate LR-AFC to release free AFC. The released AFC can easily be quantified using a fluorometer or fluorescence plate reader.</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 K141 Cathepsin K Activity Fluorometric Assay Kit. K141-100 is the same size as the 100 test size of ab65303.</p> <p>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.</p>
----	--

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

性能

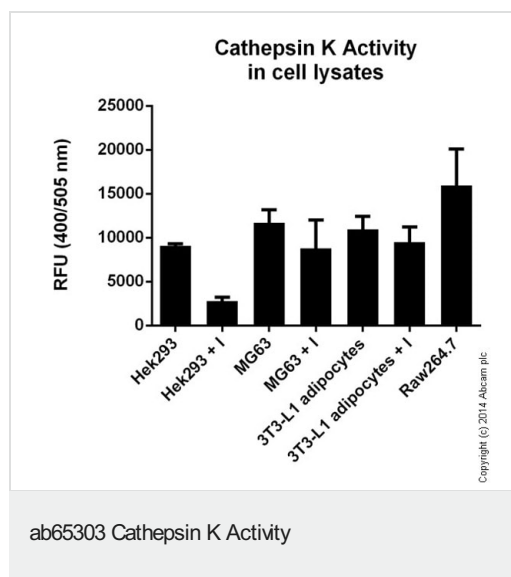
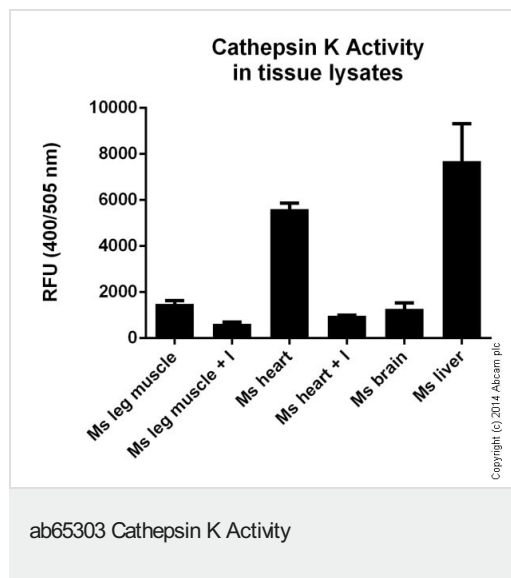
存放说明	Store at -20°C. Please refer to protocols.
------	--

组件	100 tests
CK Cell Lysis Buffer	1 x 25ml
CK Inhibitor (1mM)	1 x 20µl
CK Reaction Buffer	1 x 5ml
CK Substrate Ac-LR-AFC (10mM)	1 x 200µl

相关性 Cathepsin K is closely involved in osteoclastic bone resorption and may participate partially in the disorder of bone remodeling. Displays potent endoprotease activity against fibrinogen at acid pH. May play an important role in extracellular matrix degradation.

细胞定位 Lysosome.

图片



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