

Cathepsin K Activity Assay Kit (Fluorometric) ab65303

[7 References](#) [2 Images](#)

Overview

Product name	Cathepsin K Activity Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Tissue Extracts, Cell Lysate
Assay type	Enzyme activity
Assay time	2h 00m
Product overview	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. Visit our FAQs page for tips and troubleshooting.

Notes	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. 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.
--------------	---

Platform	Microplate reader
-----------------	-------------------

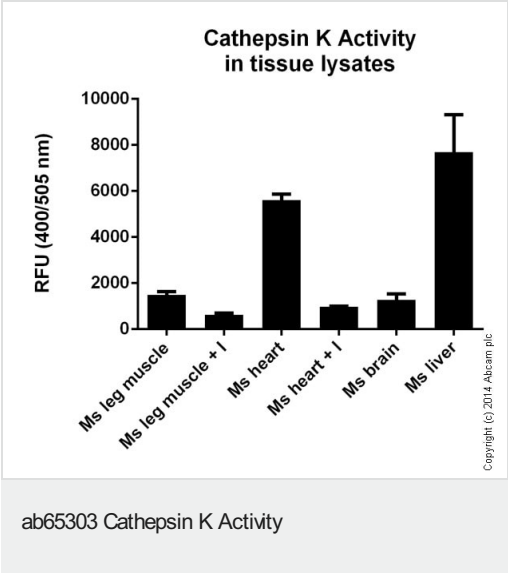
Properties

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

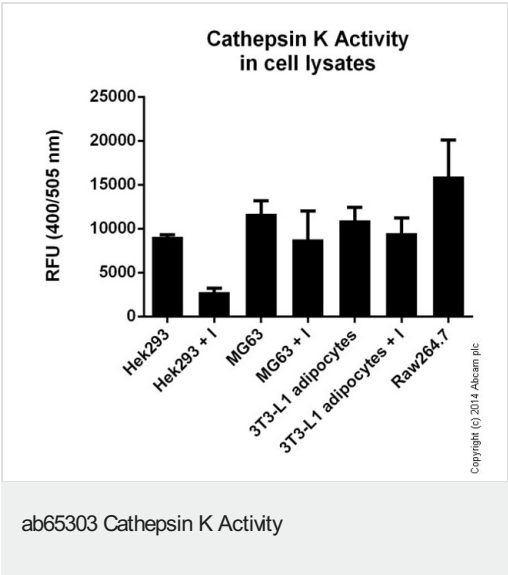
Components	100 tests
CK Cell Lysis Buffer	1 x 25ml
CK Inhibitor (1mM)	1 x 20µl
CK Reaction Buffer	1 x 5ml
CK Substrate Ac-LR-AFC (10mM)	1 x 200µl

Relevance	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.
Cellular localization	Lysosome.

Images



Cathepsin K Activity measured in tissue lysates (each with protein concentration of 4 mg/mL) after 1 hour of incubation with and without inhibitor (I).



Cathepsin K Activity measured in cell lysates (each with cells concentration of 4e7 cells/mL) after 1 hour of incubation with and without inhibitor (I).

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