

Cathepsin B Inhibitor Screening Kit (Fluorometric) ab185438

4 References 5 图像

概述

产品名称	Cathepsin B抑制剂Screening试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Inhibitor compounds
检测类型	Semi-quantitative
产品概述	Abcam's Cathepsin B Inhibitor Screening Kit (ab185438) utilizes the ability of Cathepsin B to cleave the synthetic AFC based peptide substrate to release AFC, which can be easily quantified using a fluorometer or fluorescence microplate reader. In the presence of a Cathepsin B-specific inhibitor, the cleavage of the substrate is reduced/abolished resulting in decrease or total loss of the AFC fluorescence. This high-throughput adaptable assay kit is simple, sensitive, and rapid tool to screen the potential inhibitors of Cathepsin B.

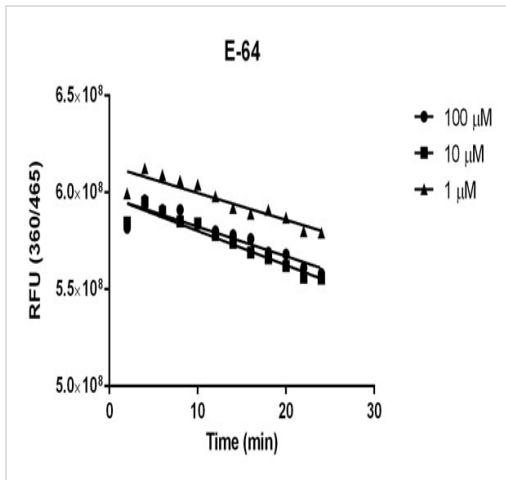
说明	<p>Visit our FAQs page for tips and troubleshooting.</p> <p>This product is manufactured by BioVision, an Abcam company and was previously called K147 Cathepsin B (CTSB) Inhibitor Screening Kit (Fluorometric). K147-100 is the same size as the 100 test size of ab185438.</p> <p>Cathepsin B (CTSB, EC 3.4.22.1) is a lysosomal cysteine proteinase that is suggested to participate in intracellular degradation and turnover of proteins. It has also been implicated in tumor invasion and metastasis.</p>
----	---

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

性能

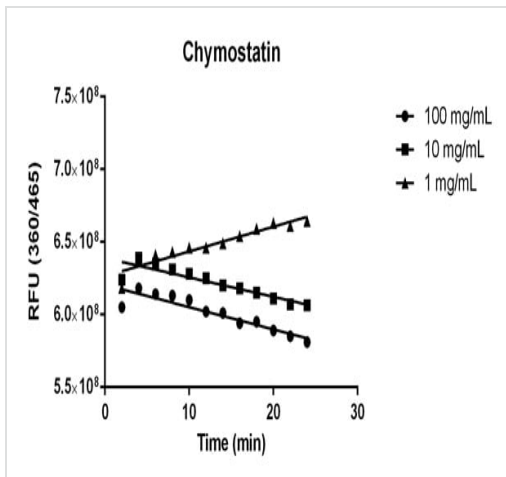
存放说明	Store at -20°C. Please refer to protocols.
------	--

组件	100 tests	100 tests
Cathepsin B Assay Buffer	1 x 15ml	1 x 15ml
Cathepsin Inhibitor	1 x 20µl	1 x 20µl



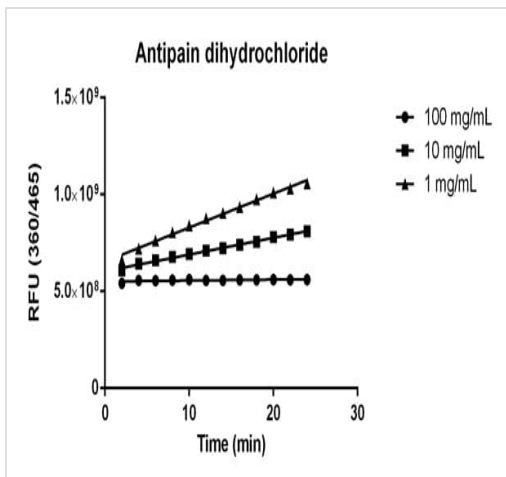
Inhibition of Cathepsin B activity by the Cathepsin B inhibitor E-64

Inhibition of Cathepsin B activity by the Cathepsin B inhibitor E-64



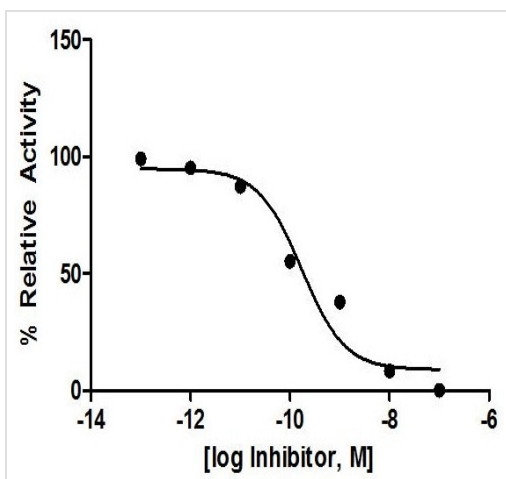
Inhibition of Cathepsin B activity by the Cathepsin B inhibitor Chymostatin

Inhibition of Cathepsin B activity by the Cathepsin B inhibitor Chymostatin



Inhibition of Cathepsin B activity by the Cathepsin B inhibitor
Antipain dihydrochloride

Inhibition of Cathepsin B activity by the Cathepsin B
inhibitor Antipain Dihydrochloride



Inhibition of Cathepsin B activity by Cathepsin B Inhibitor Screening
Kit was performed following kit protocol. This is example data only.

Inhibition of Cathepsin B activity by the Cathepsin B
Inhibitor [F-F-FMK]

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