

Caspase-1 Inhibitor Assay Kit ab102489

1 References 1 图像

概述

产品名称	Caspase-1抑制剂Assay试剂盒
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Quantitative
检测时间	1h 30m
产品概述	<p>Abcam's Caspase 1 Inhibitor Drug Detection Kit provides an effective means for screening caspase inhibitors using fluorometric methods. The assay utilizes synthetic peptide substrate YVAD-AFC (AFC, 7-amino-4-trifluoromethyl coumarin). Active caspase-1 cleaves the synthetic substrate to release free AFC which can then be quantified by fluorometry. Compounds to be screened can directly be added to the reaction and the level of inhibition of caspase-1 activity can be determined by comparison of the fluorescence intensity in samples with and without the testing inhibitors.</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K151 Caspase-1 Inhibitor Drug Screening Kit (Fluorometric). K151-100 is the same size as the 100 test size of ab102489.</p> <p>Caspases have been shown to play a crucial role in apoptosis induced by various deleterious and physiologic stimuli. Inhibition of caspases can delay apoptosis, implicating a potential role in drug screening efforts.</p>
平台	Microplate reader

性能

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

组件	100 tests
2X Reaction Buffer I	1 x 10ml
Active Caspase 1	1 x 100 units
DTT II	1 x 100µl

组件	100 tests
YVAD-AFC	1 x 0.5ml
Z-VAD-FMK	1 x 10µl

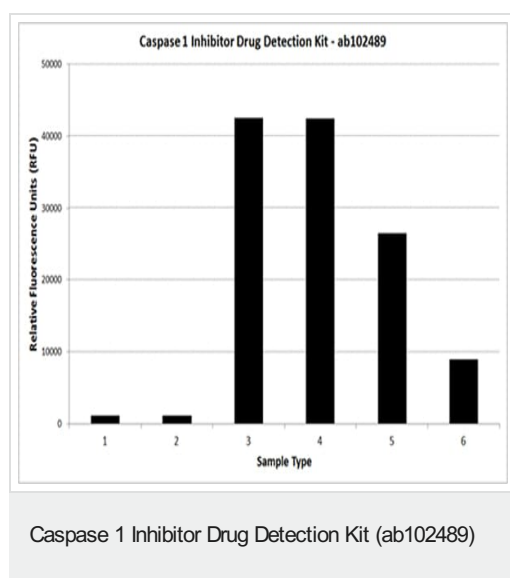
## 相关性

Caspases are a family of cysteine proteases that are key mediators of programmed cell death or apoptosis. The precursor form of all caspases is composed of a prodomain, and large and small catalytic subunits. The active forms of caspases are generated by several stimuli including ligand-receptor interactions, growth factor deprivation and inhibitors of cellular functions. All known caspases require cleavage adjacent to aspartates to liberate one large and one small subunit, which associate into a2b2 tetramer to form the active enzyme. Caspase 1 is similar to the cell death gene CED3 of *C. elegans* and regulates multiple proinflammatory cytokines, including Interleukin 1b and interferon-gamma-inducing factor. Caspase 1 plays a role in down stream of Caspase 8 which is involved in Fas-mediated apoptosis.

## 细胞定位

Cytoplasmic

图片



Caspase activity (RFU) in presence of 0 $\mu$ M - 40 $\mu$ M of z-VAD-FMK (generic caspase inhibitor), assessed using YVAD-AFC as caspase 1 substrate and following Caspase 1 Inhibitor Drug Detection Kit (ab102489) protocol. Lanes:

- 1.- Background Control: no Caspase 1, no z-VMAD-FMK
- 2.- Positive Inhibition Control: no Caspase 1, +z-VAD-FMK
- 3.- Positive Control: + Caspase 1, no z-VAD-FMK
- 4.- Caspase 1 + 2.5µM z-VAD-FMK
- 5.- Caspase 1 + 10µM z-VAD-FMK
- 6.- Caspase 1 + 40µM z-VAD-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