

Product datasheet

Caspase 3 Inhibitor Drug Detection Kit ab102491

3 图像

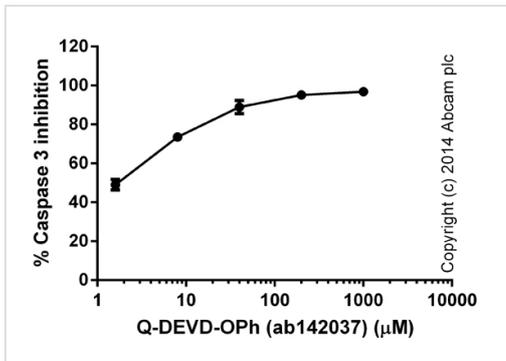
概述

产品名称	Caspase 3抑制剂Drug Detection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 3 Inhibitor Drug Detection Kit provides an effective means for screening caspase inhibitors using fluorometric methods. The assay utilizes synthetic peptide substrate DEVD-AFC (AFC, 7-amino-4-trifluoromethyl coumarin). Active caspase 3 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 3 activity can be determined by comparison of the fluorescence intensity in samples with and without the testing inhibitors.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
说明	<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>
经测试应用	适用于: Functional Studies
平台	Microplate

性能

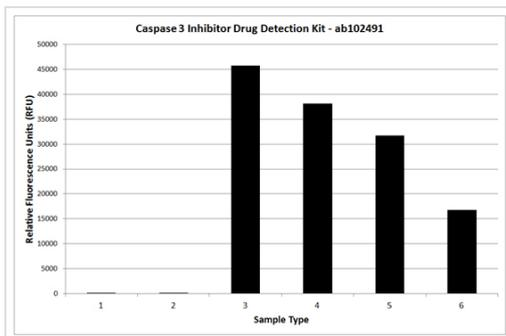
存放说明 Store at -20°C. Please refer to protocols.

组件	100 tests
2X Reaction Buffer	1 x 10ml
Active Caspase-3	1 x 100 units
Caspase Substrate DEVD-AFC	1 x 0.5ml
Caspase-3 Inhibitor	1 x 10µl



Titration of the Caspase 3 specific inhibitor Q-DEVD-Oph (ab142037) (duplicates; +/- SD).

Functional assays: Caspase 3 Inhibitor Drug Detection Kit (ab102491)



Caspase 3 Inhibitor Drug Detection Kit (ab102491)

Caspase activity (RFU) in presence of 0μM - 40μM of z-VAD-FMK (generic caspase inhibitor), assessed using DEVD-AFC as caspase 3 substrate and following Caspase 3 Inhibitor Drug Detection Kit (ab102491) protocol. Lanes:

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

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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