

Product datasheet

Caspase-8 Inhibitor Assay Kit ab102496

1 Image

Overview

| | |
|-------------------------|---|
| Product name | Caspase-8 Inhibitor Assay Kit |
| Detection method | Fluorescent |
| Sample type | Adherent cells, Suspension cells |
| Assay type | Quantitative |
| Assay time | 1h 30m |
| Product overview | <p>Abcam's Caspase 8 Inhibitor Drug Detection Kit provides an effective means for screening caspase inhibitors using fluorometric methods. The assay utilizes synthetic peptide substrate IETD-AFC (AFC, 7-amino-4-trifluoromethyl coumarin). Active caspase 8 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 8 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> |
| Notes | <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> |
| Platform | Microplate reader |

Properties

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

| Components | 100 tests |
|------------------------------|---------------|
| Active Caspase 8 | 1 x 100 units |
| Caspase Inhibitor, Z-VAD-FMK | 1 x 10µl |
| Caspase Substrate IETD-AFC | 1 x 0.5ml |
| DTT | 1 x 100µl |
| Reaction buffer | 1 x 10ml |

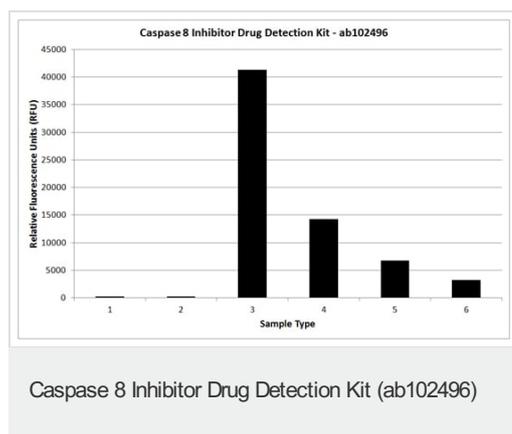
Relevance

Caspases are cysteine proteases, expressed as inactive precursors, that mediate apoptosis by proteolysis of specific substrates. Caspases have the ability to cleave after aspartic acid residues. There are two classes of caspases involved in apoptosis; initiators (activation by receptor cluster) and effectors (activation by mitochondrial permeability transition). Proapoptotic signals autocatalytically activate initiator caspases, such as Caspase 8 and Caspase 9. Activated initiator caspases then process effector caspases, such as Caspase 3 and Caspase 7, which in turn cause cell collapse.

Cellular localization

Cytoplasmic

Images



Caspase activity (RFU) in presence of 0µM - 40µM of z-VAD-FMK (generic caspase inhibitor), assessed using IETD-AFC as caspase 8 substrate and following Caspase 8 Inhibitor Drug Detection Kit (ab102496) protocol. Lanes:

- 1.- Background Control: no Caspase 8, no z-VAD-FMK
- 2.- Positive Inhibition Control: no Caspase 8, +z-VAD-FMK
- 3.- Positive Control: + Caspase 8, no z-VAD-FMK
- 4.- Caspase 8 + 2.5µM z-VAD-FMK
- 5.- Caspase 8 + 10µM z-VAD-FMK
- 6.- Caspase 8 + 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.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