

Caspase-2 Inhibitor Assay Kit ab102490

[1 Image](#)

Overview

| | |
|-------------------------|--|
| Product name | Caspase-2 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 2 Inhibitor Drug Detection Kit provides an effective means for screening caspase inhibitors using fluorometric methods. The assay utilizes synthetic peptide substrate VDVAD-AFC (AFC, 7-amino-4-trifluoromethyl coumarin). Active caspase 2 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 2 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>This product is manufactured by BioVision, an Abcam company and was previously called K152 Caspase-2 Inhibitor Drug Screening Kit (Fluorometric). K152-100 is the same size as the 100 test size of ab102490.</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> |
| Platform | Microplate reader |

Properties

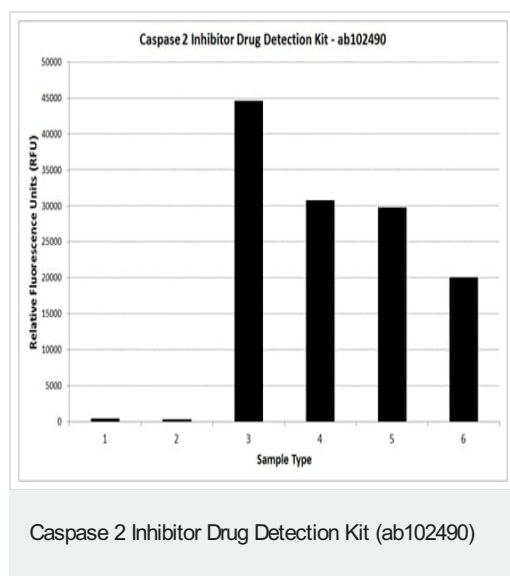
Storage instructions Store at -20°C. Please refer to protocols.

| Components | 100 tests |
|------------------------------|---------------|
| 2X Reaction Buffer | 1 x 10ml |
| Active Caspase-2 | 1 x 100 units |
| Caspase Inhibitor, Z-VAD-FMK | 1 x 10µl |

| Components | 100 tests |
|-----------------------------|-----------|
| Caspase Substrate VDVAD-AFC | 1 x 0.5ml |
| DTT | 1 x 100µl |

| | |
|---|---|
| Function | Involved in the activation cascade of caspases responsible for apoptosis execution. Might function by either activating some proteins required for cell death or inactivating proteins necessary for cell survival. |
| Tissue specificity | Expressed at higher levels in the embryonic lung, liver and kidney than in the heart and brain. In adults, higher level expression is seen in the placenta, lung, kidney, and pancreas than in the heart, brain, liver and skeletal muscle. |
| Sequence similarities | Belongs to the peptidase C14A family. Contains 1 CARD domain. |
| Post-translational modifications | The mature protease can process its own propeptide, but not that of other caspases. |

Images



Caspase activity (RFU) in presence of 0µM - 40µM of z-VAD-FMK (generic caspase inhibitor), assessed using VDVAD-AFC as caspase 2 substrate and following Caspase 2 Inhibitor Drug Detection Kit (ab102490) protocol. Lanes:

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