

**ab102498**

# **Caspase 10 Inhibitor Drug Detection Kit**

## **Instructions for Use**

For the rapid, sensitive and accurate detection of Caspase 10 inhibition by various compounds

This product is for research use only and is not intended for diagnostic use.

# Table of Contents

---

Table of Contents	1
1. Overview	2
2. Protocol Summary	3
3. Components and Storage	4
4. Assay Protocol	6
5. Factors to consider for caspase activity assays	8
6. Troubleshooting	10

# 1. Overview

---

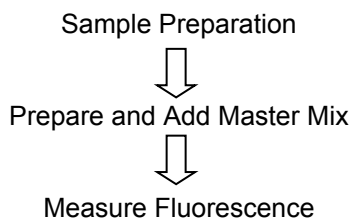
Caspases have been shown to play a crucial role in apoptosis induced by various deleterious and physiologic stimuli. Inhibition of caspases can delay apoptosis, implicating a potential role in drug screening efforts.

Abcam's Caspase 10 Inhibitor Drug Detection Kit provides an effective means for screening caspase inhibitors using fluorometric methods. The assay utilizes synthetic peptide substrate AEVD-AFC (AFC, 7-amino-4-trifluoromethyl coumarin). Active caspase 10 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 10 activity can be determined by comparison of the fluorescence intensity in samples with and without the testing inhibitors.

## 2. Protocol Summary

---



### 3. Components and Storage

---

#### A. Kit Components

Item	Quantity
2X Reaction Buffer	10 mL
Caspase Substrate AEVD-AFC (1 mM)	0.5 mL
DTT (1 M)	100 $\mu$ L
Active Caspase 10 (Lyophilized)	100 units
Caspase Inhibitor, Z-VAD-FMK (2 mM)	10 $\mu$ L

\* Store kit at -20°C. All components are stable for 6 months under proper storage conditions.

Protect AEVD-AFC from light.

REACTION BUFFER: After thawing, store the 2X Reaction Buffer at +4°C. Aliquot enough 2X Reaction Buffer for the number of assays to be performed. Add DTT to the 2X Reaction Buffer immediately before use (10 mM final concentration: add 10  $\mu$ l of 1.0 M DTT stock per 1 ml of 2X Reaction Buffer).

ACTIVE CASPASE 10: Reconstitute in 550  $\mu$ l 2X Reaction Buffer. Aliquot and immediately store at -80°C.

## **B. Additional Materials Required**

- Microcentrifuge
- Pipettes and pipette tips
- Fluorescent microplate reader
- 96-well plate
- Orbital shaker

## 4. Assay Protocol

---

1. Prepare test samples in dH<sub>2</sub>O up to a final volume of 50 µl/well. Add 5 µl of Active Caspase 10. Mix well.

Prepare a **background control** by omitting the Active Caspase 10 from the reaction mixture.

Prepare a **positive inhibition control** by adding 1 µl of the Caspase 10 Inhibitor (provided with the kit) instead of your testing inhibitor.

**Note:**

This product detects proteolytic activity. Do not use protease inhibitors in the sample preparation step as it might interfere with the assay.

2. Prepare a Master Mix for each assay containing the following:

2X Reaction Buffer	45 µL
(containing 10 mM DTT)	
1 mM AEVD-AFC substrate	5 µL
(50 µM final concentration)	
3. Mix well and add 50 µl of the Master Mix to each well to start the reaction. Incubate at 37°C for 0.5-1 hour.
4. Read samples in a fluorescence plate reader equipped with a 400-nm excitation filter and 505-nm emission filter.

Compare the fluorescence intensity of the test samples with samples containing no inhibitors to determine the inhibition efficiency of the testing inhibitors.

## 5. Factors to consider for caspase activity assays

Three major factors need to be taken into account when using caspase activity assays:

1. The substrate in a particular assay is not necessarily specific to a particular caspase.

Cleavage specificities overlap so reliance on a single substrate/assay is not recommended. Other assays, such as Western blot, use of fluorescent substrates e.g. FRET assays should be used in combination with caspase activity assays.

2. The expression and abundance of each caspase in a particular cell type and cell line will vary.
3. As the activation and cleavage of caspases in the cascade will change over time, you should consider when particular caspase will be at its peak concentration e.g. after 3 hours, after 20 hours etc.

The table below shows the known cross-reactivities with other caspases.

Classification of caspases based on synthetic substrate preference, does not reflect the real caspase substrate preference *in vivo* and may provide inaccurate information for discriminating amongst caspase activities. Thus, caution is advised in applying the intrinsic

tetrapeptide preferences to predict the targets of individual caspases.

### Apoptotic Initiator Caspases

Caspase	Cleavage motif	Inhibitor motif	Cross-reactivity with other caspase:											
			1	2	3	4	5	6	7	8	9	10		
<b>Caspase 2</b>	VDVAD				Y				Y					
<b>Caspase 8</b>	IETD	IETD, LETD			Y			Y						Y
<b>Caspase 9</b>	LEHD				Y			Y		Y				Y
<b>Caspase 10</b>	AEVD				Y				Y	Y	?			

## 6. Troubleshooting

<b>Problem</b>	<b>Reason</b>	<b>Solution</b>
Assay not working	Reaction buffer at wrong temperature	Assay buffer must be at 4°C
	Protocol step missed	Re-read and follow the protocol exactly
	Plate read at incorrect wavelength	Ensure you are using appropriate reader and filter settings (refer to datasheet)
	Unsuitable microtiter plate for assay	Fluorescence: Black plates (clear bottoms); Luminescence: White plates; Colorimetry: Clear plates. If critical, datasheet will indicate whether to use flat- or U-shaped wells
Unexpected results	Measured at wrong wavelength	Use appropriate reader and filter settings described in datasheet
	Samples contain impeding substances	Troubleshoot and also consider deproteinizing samples
	Unsuitable sample type	Use recommended samples types as listed on the datasheet

General Issues	Improperly thawed components	Thaw all components completely and mix gently before use
	Incorrect incubation times or temperatures	Refer to datasheet & verify the correct incubation times and temperatures
	Incorrect volumes used	Use calibrated pipettes and aliquot correctly
	Air bubbles formed in the well/tube	Pipette gently against the wall of the well/tubes
	Substituting reagents from older kits/ lots	Use fresh components from the same kit

**For further technical questions please do not hesitate to contact us by email ([technical@abcam.com](mailto:technical@abcam.com)) or phone (select “*contact us*” on [www.abcam.com](http://www.abcam.com) for the phone number for your region).**







**UK, EU and ROW**

Email: [technical@abcam.com](mailto:technical@abcam.com) | Tel: +44-(0)1223-696000

**Austria**

Email: [wissenschaftlicherdienst@abcam.com](mailto:wissenschaftlicherdienst@abcam.com) |  
Tel: 019-288-259

**France**

Email: [supportscientifique@abcam.com](mailto:supportscientifique@abcam.com) | Tel: 01-46-94-62-96

**Germany**

Email: [wissenschaftlicherdienst@abcam.com](mailto:wissenschaftlicherdienst@abcam.com) | Tel: 030-896-779-154

**Spain**

Email: [soportecientifico@abcam.com](mailto:soportecientifico@abcam.com) | Tel: 911-146-554

**Switzerland**

Email: [technical@abcam.com](mailto:technical@abcam.com)  
Tel (Deutsch): 0435-016-424 | Tel (Français): 0615-000-530

**US and Latin America**

Email: [us.technical@abcam.com](mailto:us.technical@abcam.com) | Tel: 888-77-ABCAM (22226)

**Canada**

Email: [ca.technical@abcam.com](mailto:ca.technical@abcam.com) | Tel: 877-749-8807

**China and Asia Pacific**

Email: [hk.technical@abcam.com](mailto:hk.technical@abcam.com) | Tel: 400 921 0189 / +86 21 2070 0500

**Japan**

Email: [technical@abcam.co.jp](mailto:technical@abcam.co.jp) | Tel: +81-(0)3-6231-0940

[www.abcam.com](http://www.abcam.com) | [www.abcam.cn](http://www.abcam.cn) | [www.abcam.co.jp](http://www.abcam.co.jp)