

## Product datasheet

# Caspase-3 Assay Kit (Fluorometric) ab39383

★★★★★ 1 Abreviews 27 References 2 Images

### Overview

<b>Product name</b>	Caspase-3 Assay Kit (Fluorometric)
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Tissue Extracts, Cell Lysate
<b>Assay type</b>	Enzyme activity
<b>Assay time</b>	2h 00m
<b>Product overview</b>	Caspase-3 Assay Kit (Fluorometric) (ab39383) provides a simple and convenient means for assaying DEVD-dependent caspase activity.

The Caspase-3 assay protocol is based on detection of cleavage of substrate DEVD-AFC (AFC: 7-amino-4-trifluoromethyl coumarin).

DEVD-AFC emits blue light ( $\lambda$  max = 400 nm). On cleavage of the substrate by CPP32 or related caspases, free AFC emits a yellow-green fluorescence (Ex/Em = 400/505 nm).

The signal can be quantified using a fluorometer or a fluorescence microtiter plate reader. Comparison of the fluorescence of AFC from an apoptotic sample with an uninduced control allows determination of the fold increase in Caspase-3/ CPP32 activity.

Caspase-3 assay protocol summary:

- lyse cells / homogenize and lyse tissues in lysis buffer
- incubate on ice for 10 min
- add reaction buffer and DEVD-AFC substrate and incubate for 1-2 hr at 37°C
- analyze with fluorometer or microplate reader

**Notes** Due to the nature of the substrate, this assay also detects caspase-7 activity.

#### Other caspase and apoptosis assays

Review the full set of [caspase assays](#), or the [apoptosis assay and apoptosis marker guide](#).

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

**Platform** Microplate reader

## Properties

---

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

Components	100 tests
2X Reaction Buffer	4 x 2ml
Cell Lysis Buffer	1 x 100ml
DEVD AFC	1 x 500µl
DTT	1 x 400µl

---

**Function** Involved in the activation cascade of caspases responsible for apoptosis execution. At the onset of apoptosis it proteolytically cleaves poly(ADP-ribose) polymerase (PARP) at a '216-Asp-Gly-217' bond. Cleaves and activates sterol regulatory element binding proteins (SREBPs) between the basic helix-loop-helix leucine zipper domain and the membrane attachment domain. Cleaves and activates caspase-6, -7 and -9. Involved in the cleavage of huntingtin.

**Tissue specificity** Highly expressed in lung, spleen, heart, liver and kidney. Moderate levels in brain and skeletal muscle, and low in testis. Also found in many cell lines, highest expression in cells of the immune system.

**Sequence similarities** Belongs to the peptidase C14A family.

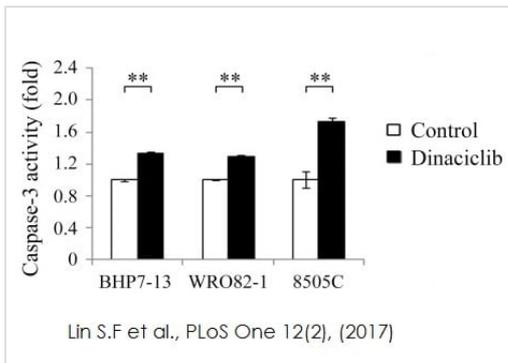
**Post-translational modifications** Cleavage by granzyme B, caspase-6, caspase-8 and caspase-10 generates the two active subunits. Additional processing of the propeptides is likely due to the autocatalytic activity of the activated protease. Active heterodimers between the small subunit of caspase-7 protease and the large subunit of caspase-3 also occur and vice versa. S-nitrosylated on its catalytic site cysteine in unstimulated human cell lines and denitrosylated upon activation of the Fas apoptotic pathway, associated with an increase in intracellular caspase activity. Fas therefore activates caspase-3 not only by inducing the cleavage of the caspase zymogen to its active subunits, but also by stimulating the denitrosylation of its active site thiol.

**Cellular localization** Cytoplasm.

---

## Images

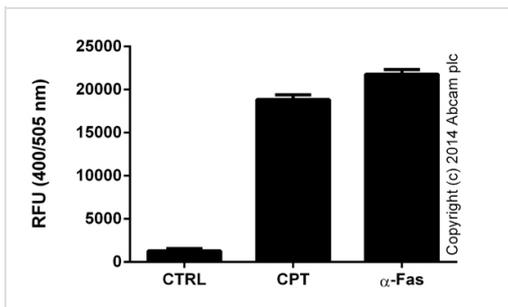
---



Functional assays: Caspase-3 assay kit (ab39383)

Image from Lin S.F et al., PLoS One 12(2), Fig4B. doi: 10.1371/journal.pone.0172315. Reproduced under the Creative Commons license <http://creativecommons.org/licenses/by/4.0/>

Caspase-3 activity was analyzed using fluorometric assay kit (ab39383). Thyroid cancer cells were plated at  $1 \times 10^6$  cells in 10 mL of media and incubated overnight. Cells were treated with Dinaciclib (25 nM) for 24 hours. Adherent cells ( $5 \times 10^5$ ) were collected, centrifuged, lysed using 50  $\mu$ L of lysis buffer on ice for 10 min, incubated with DEVD-AFC substrate and reaction buffer at 37°C for 1.5 hours. Caspase-3 activity was detected by spectrophotometry. The fluorescence intensity of the treated samples was compared with that of control samples to determine the fold-increase in caspase activity. Each condition was performed in duplicate.



Functional assays: Caspase-3 Assay Kit (Fluorometric) (ab39383)

Caspase-3 activity in Jurkat lysates ( $6.6 \times 10^5$  cells) following 20 hour exposure to 2  $\mu$ M Camptothecin (ab120115) or 10 ng/mL anti-Fas Ab (MBL). Background signal subtracted, duplicates +/- SD.

**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