

# Cleaved Caspase-3 Staining Kit (Red) ab65617

5 References 画像数 1

### 製品の概要

製品名	Cleaved Caspase-3 Staining Kit (Red)
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Enzyme activity
全工程の試験時間	2h 00m
製品の概要	Cleaved Caspase-3 Staining Kit (Red) ab65617 provides a convenient means for detecting activated Caspase-3 in living cells. The assay utilizes the Caspase-3 inhibitor DEVD-FMK conjugated to sulfo-rhodamine (Red-DEVD-FMK) as the fluorescent in situ marker. Red-DEVD-FMK is cell permeable, nontoxic, and irreversibly binds to activated caspase-3 in apoptotic cells.
特記事項	This product is manufactured by BioVision, an Abcam company and was previously called K193 CaspGLOW™ Red Active Caspase-3 Staining Kit. K193-100 is the same size as the 100 test size of ab65617.
試験プラットフォーム	Microplate reader, Fluor. microscope, Flow cyt.

### 製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
Red-DEVD-FMK	1 x 100µl
Wash Buffer IV	2 x 100ml
Z-VAD-FMK	1 x 10µl

**関連性**

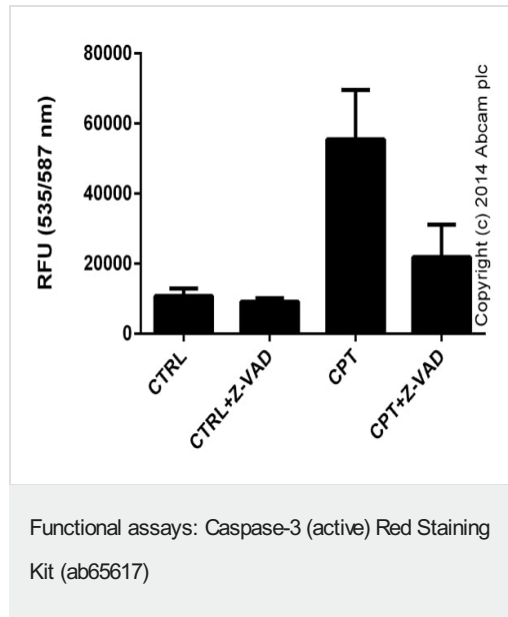
The caspase family of cysteine proteases play a key role in apoptosis. Caspase 3 (also known as CPP32, YAMA and apopain) is the most extensively studied apoptotic protein among caspase family members. Caspase 3 is synthesized as an inactive pro enzyme that is processed in cells undergoing apoptosis by self proteolysis and/or cleavage by other upstream proteases (e.g. Caspases 8, 9 and 10). The processed form of Caspase 3 consists of large (17kD) and small (12kD) subunits which associate to form an active enzyme. Caspase 3 is cleaved at Asp28 - Ser29 and Asp175 - Ser176. The active Caspase 3 proteolytically cleaves and activates other caspases (e.g. Caspases 6, 7 and 9), as well as relevant targets in the cells (e.g. PARP and DFF). Alternative splicing of this gene results in two transcript variants which encode the same

protein. In immunohistochemical studies Caspase 3 expression has been shown to be widespread but not present in all cell types (e.g. commonly reported in epithelial cells of skin, renal proximal tubules and collecting ducts). Differences in the level of Caspase 3 have been reported in cells of short lived nature (eg germinal centre B cells) and those that are long lived (e.g. mantle zone B cells). Caspase 3 is the predominant caspase involved in the cleavage of amyloid beta 4A precursor protein, which is associated with neuronal death in Alzheimer's disease.

細胞内局在

Cytoplasmic

## 画像



Caspase-3 activity in Jurkat cells ( $3 \times 10^5$  cells) following 24 hour exposure to 2  $\mu$ M Camptothecin ([ab120115](#)) with or without 50  $\mu$ M caspase inhibitor Z-VAD(OMe)-FMK ([ab120487](#)). Background signal subtracted, duplicates;  $\pm$  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.co.jp/abpromise> or contact our technical team.

## Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors