

# Caspase 3 (active) FITC Staining Kit ab65613

6 References 1 图像

## 概述

|       |   |
|-------|---|
| 产品名称  | Caspase 3 (active) FITC Staining试剂盒   |
| 样品类型  | Adherent cells, Suspension cells  |
| 检测类型  | Enzyme activity   |
| 检测时间  | 2h 00m  |
| 产品概述  | <p>Abcam's Caspase 3 (active) FITC Staining Kit provides a convenient means for sensitive detection of activated caspase 3 in living cells. The assay utilizes the caspase 3 inhibitor, DEVD-FMK, conjugated to FITC (FITC-DEVD-FMK) as a marker. FITC-DEVD-FMK is cell permeable, non-toxic, and irreversibly binds to activated caspase 3 in apoptotic cells. The FITC label allows detection of activated caspase-3 in apoptotic cells directly by fluorescence microscopy, flow cytometry, or fluorescence plate reader.</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p> |
| 说明    | Activation of caspases plays a central role in apoptosis.   |
| 经测试应用 | 适用于: FM, Functional Studies, Flow Cyt   |

## 性能

存放说明 Store at -20°C. Please refer to protocols.

| 组件            | 100 tests |
|---------------|-----------|
| FITC-DEVD-FMK | 1 x 100µl |
| Wash Buffer   | 2 x 100ml |
| Z-VAD-FMK     | 1 x 10µl  |

**相关性**

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.

**细胞定位**

Cytoplasmic

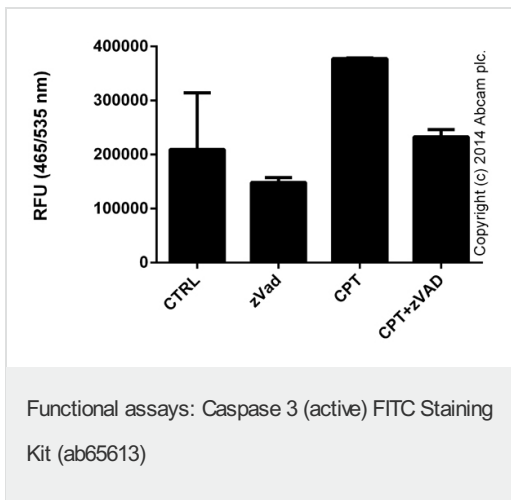
**应用**

Our [Abpromise guarantee](#) covers the use of **ab65613** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

| 应用                 | Ab评论 | 说明                                       |
|--------------------|------|--|
| FM                 |      | Use at an assay dependent concentration. |
| Functional Studies |      | Use at an assay dependent dilution.      |
| Flow Cyt           |      | Use at an assay dependent dilution.      |

**Caspase 3 (active) FITC Staining Kit 图像**



Caspase 3 in Jurkat cells following 24 hour exposure to 2 uM Camptothecin (CPT; [ab120115](#)) with or without 20 nM caspase inhibitor Z-VAD(OMe)-FMK (zVAD; [ab120487](#)). Background signal subtracted, duplicates; +/- SD.

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

**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 <http://www.abcam.cn/abpromise> or contact our technical team.

## Terms and conditions

---

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