

Product datasheet

pan Caspase (active) FITC Staining Kit ab65611

製品の概要

製品名	pan Caspase (active) FITC Staining Kit
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	<p>Abcam's Fluorescein Caspase Staining Kit provides a convenient and sensitive means for detecting activated caspases in living cells. The assay utilizes the caspase family inhibitor VAD-FMK conjugated to FITC (FITC-VAD-FMK) as a marker. FITC-VAD-FMK is cell permeable, nontoxic, and irreversibly binds to activated caspases in apoptotic cells.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
特記事項	Activation of caspases plays a central role in apoptosis.
アプリケーション	適用あり: FM, Functional Studies, Flow Cyt

製品の特性

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

内容	100 tests
FITC-VAD-FMK	1 x 100µl
Wash Buffer	2 x 100ml
Z-VAD-FMK	1 x 10µl

関連性 Caspases are members of the cysteine-aspartic acid protease (caspase) family. Sequential activation of caspases plays a central role in the execution-phase of cell apoptosis. Caspases exist as inactive proenzymes which undergo proteolytic processing at conserved aspartic residues to produce 2 subunits, large and small, that dimerize to form the active enzyme.

細胞内局在 Cytoplasmic

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab65611** in the following tested applications.

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

アプリケーション	Abreviews	特記事項
FM		Use at an assay dependent concentration.
Functional Studies		Use at an assay dependent dilution.
Flow Cyt		Use at an assay dependent dilution.

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.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