

## Product datasheet

# pan Caspase (active) FITC Staining Kit ab65611

### 概述

<b>产品名称</b>	pan Caspase (active) FITC 荧光 Staining 试剂盒
<b>样品类型</b>	Adherent cells, Suspension cells
<b>检测类型</b>	Enzyme activity
<b>检测时间</b>	2h 00m
<b>产品概述</b>	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. Visit our <a href="#">FAQs page</a> for tips and troubleshooting.
<b>说明</b>	Activation of caspases plays a central role in apoptosis.
<b>经测试应用</b>	<b>适用于:</b> FM, Functional Studies, Flow Cyt

### 性能

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

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

**相关性** 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.

**细胞定位** Cytoplasmic

### 应用

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.

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.

**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