

Product datasheet

pan Caspase (active) FITC Staining Kit ab65611

4 References

Overview

Product name	pan Caspase (active) FITC Staining Kit
Sample type	Adherent cells, Suspension cells
Assay type	Enzyme activity
Assay time	2h 00m
Product overview	<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>
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K180 CaspGLOW™ Fluorescein Active Caspase Staining Kit. K180-100 is the same size as the 100 test size of ab65611.</p> <p>Activation of caspases plays a central role in apoptosis.</p> <p>Other caspase and apoptosis assays</p> <p>Review the full set of caspase assays, or the apoptosis assay and apoptosis marker guide.</p>
Platform	Flow cytometer, Fluorescence microscope

Properties

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

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

Relevance

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.

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