

Product datasheet

Caspase Assay Kit (orange, VAD-FMK) ab270770

1 References

Overview

Product name Caspase Assay Kit (orange, VAD-FMK)

Product overview Caspases play important roles in apoptosis and inflammation. Caspase Assay Kit (orange, VAD-FMK) (ab270770) is used by researchers seeking to quantitate apoptosis via caspase activity in cultured cells and tissues.

The FLICA reagent SR-VAD-FMK enters each cell and irreversibly binds to activated caspases. Because the SR-VAD-FMK FLICA reagent becomes covalently coupled to the active enzymes, it is retained within the cell, while any unbound SR-VAD-FMK FLICA reagent diffuses out of the cell and is washed away. The remaining red fluorescent signal is a direct measure of the active caspase enzyme activity present in the cell at the time the reagent was added. Cells that contain the bound FLICA can be analyzed by a fluorescence plate reader, fluorescence microscopy or flow cytometry. Cells labeled with the FLICA reagent may be read immediately or preserved for 16 hours using the fixative included in the kit. Unfixed samples may also be analyzed with Hoechst 33342 to detect changes in nuclear morphology.

Produced using a non-baculovirus expression system.

Tested applications **Suitable for:** Flow Cyt, FM

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	25 tests	100 tests
10X Apoptosis Wash Buffer	1 x 15ml	1 x 60ml
Caspase inhibitor reagent SR-VAD-FMK	1 vial	4 vials
Fixative	1 x 6ml	1 x 6ml
Hoechst 33342, 200 µg/mL	1 x 1ml	1 x 1ml

Applications

The Abpromise guarantee Our [Abpromise guarantee](#) covers the use of ab270770 in the following tested applications. The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Flow Cyt		Use at an assay dependent concentration.
FM		Use at an assay dependent concentration.

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