

Product datasheet

Caspase 3/7 Staining Kit (Far Red) ab270785

Overview

Product name Caspase 3/7 Staining Kit (Far Red)

Product overview Caspases play important roles in apoptosis and inflammation. Caspase 3/7 Staining Kit (Far Red) are used by researchers seeking to quantitate apoptosis via caspase activity in cultured cells. The fluorescent 660-DEVD-FMK caspase-3/7 inhibitor reagent allows researchers to assess caspase activation.

The fluorescent reagent 660-DEVD-FMK enters each cell and irreversibly binds to activated Caspase 3/7. Because the 660-DEVD-FMK reagent becomes covalently coupled to the active enzymes, it is retained within the cell, while any unbound 660-DEVD-FMK reagent diffuses out of the cell and is washed away. The remaining far 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 660-DEVD-FMK can be analyzed by fluorescence microscopy or flow cytometry. Cells labeled with the 660-DEVD-FMK reagent may be read immediately or preserved for 16 hours using the fixative. Unfixed samples may also be analyzed with Hoechst stain to detect changes in nuclear morphology.

Produced using a non-baculovirus expression system.

The target for this kit is Caspase 3/7.

NOTE: The number of tests per kit depends on the application being tested. This kit provides 50 tests for flow and 25 for imaging.

FLICA 660-DEVD-FMK caspase-3/7 inhibitor reagent contains the preferred binding sequence, Asp-Glu-Val-Asp (DEVD) for executioner caspases -3 and -7. This preferred caspase-3 and -7 binding sequence, DEVD, is labeled at the amino terminus end with a far-red fluorescent 660 dye and linked at the carboxyl end to a fluoromethyl ketone (FMK) reactive entity. The resulting cell permeant, fluorescent molecule, 660-DEVD-FMK, optimally excites at 660 nm and emits between 685-690 nm. A conventional red HeNe laser with a 633 nm excitation provides excellent excitation efficiency, enabling cells labeled with 660-DEVD-FMK to be analyzed with most flow cytometers, as well as fluorescence microscopes equipped with electronic grey scale image capabilities.

Kits are available for other caspases:

Poly Caspase Staining Kit (Far Red) ([ab270783](#))

Caspase 1 Staining Kit (Far Red) ([ab27084](#))

Tested applications

Suitable for: FM, Flow Cyt

Properties

Storage instructions

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

Components	50 tests
10X Apoptosis Wash Buffer	1 x 15ml
Fixative	1 x 6ml
660-DEVD-FMK caspase-3/7 inhibitor reagent	1 vial

Applications

The Abpromise guarantee

Our [Abpromise guarantee](#) covers the use of ab270785 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
FM		Use at an assay dependent concentration.
Flow Cyt		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