

Caspase 3/7 Staining Kit (Far Red) ab270785

2 References

医薬用外劇物

製品の概要

製品名	Caspase 3/7 Staining Kit (Far Red)
製品の概要	<p>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.</p> <p>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.</p> <p>Produced using a non-baculovirus expression system.</p> <p>The target for this kit is Caspase 3/7.</p> <p>NOTE: The number of tests per kit depends on the application being tested. This kit provides 50 tests for flow and 25 for imaging.</p> <p>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.</p>

Kits are available for other caspases:

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

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

アプリケーション

適用あり: FM, Flow Cyt

製品の特性

保存方法

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

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

アプリケーション

The Abpromise guarantee

Abpromise保証は、次のテスト済みアプリケーションにおけるab270785の使用に適用されます

アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
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.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