

Product datasheet

Caspase 8 Assay Kit (Red) ab270786

Overview

Product name Caspase 8 Assay Kit (Red)

Product overview Caspases play important roles in apoptosis and inflammation. Caspase 8 Assay Kit (Red) (ab270786) is used by researchers seeking to quantitate apoptosis via caspase activity in cultured cells and tissues.

The reagent SR-LETD-FMK enters each cell and irreversibly binds to activated caspase-8. Because the SR-LETD-FMK reagent becomes covalently coupled to the active enzymes, it is retained within the cell, while any unbound SR-LETD-FMK reagent diffuses out of the cell and is washed away. The remaining red fluorescent signal is a direct measure of the active caspase-8 enzyme activity present in the cell at the time the reagent was added. Cells that contain the bound SR-LETD-FMK can be analyzed by a fluorescence plate reader, fluorescence microscopy, or flow cytometry. Cells labeled with the SR-LETD-FMK 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.

There is no interference from pro-caspases or inactive forms of the enzymes. If the treatment is causing cell death via apoptosis, apoptotic cells will have an elevated level of caspase activity relative to nonapoptotic or negative control cells and fluoresce with SR-LETD-FMK.

Produced using a non-baculovirus expression system.

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 SR-LETD-FMK	1 vial	4 vials
Fixative	1 x 6ml	1 x 6ml
Hoechst 33342, 200 µg/mL	1 vial	1 vial

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