

Product datasheet

Caspase-3, Caspase-8 and Caspase-9 Multiplex Activity Assay Kit (Fluorometric) ab219915

[3 References](#) [3 Images](#)

Overview

Product name	Caspase-3, Caspase-8 and Caspase-9 Multiplex Activity Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Adherent cells, Suspension cells
Assay type	Quantitative
Product overview	<p>Caspase 3, Caspase 8 and Caspase 9 Multiplex Activity Assay Kit (Fluorometric) (ab219915) provides a simple and convenient tool to quantify caspase 3, caspase 8 and caspase 9 activity in cells that are undergoing apoptosis. The assay can also be used to screen Caspase 3, caspase 8 and caspase 9 inhibitors. This product is designed to simultaneously monitor key caspases involved in apoptosis: the initiator caspases caspase 8 and caspase 9, and the executioner Caspase 3. The kit uses DEVD-ProRed, IETD-R110 and LEHD-AMC as fluorogenic indicators for Caspase 3, caspase 8 and caspase 9 activity respectively. Upon caspase cleavage, three distinct fluorophores are released: ProRed (red fluorescence), R110 (green fluorescence) and AMC (blue fluorescence), which can be readily monitored in a single assay due to their nice spectral separation.</p> <p>This product has been optimized for use in a microplate reader in 96-well plate, providing enough reagent to perform 100 tests for each caspase.</p>
Notes	<p>This product is designed to simultaneously monitor key caspases involved in apoptosis: the initiator caspases caspase 8 and caspase 9, and the executioner Caspase 3 (which share the same substrate). The kit uses DEVD-ProRed™, IETD-R110 and LEHD-AMC as fluorogenic indicators for Caspase 3, caspase 8 and caspase 9 activity respectively. Upon caspase cleavage, three distinct fluorophores are released: ProRed™ (red fluorescence), R110 (green fluorescence) and AMC (blue fluorescence), which can be readily monitored in a single assay due to their nice spectral separation.</p> <p>This product has been optimized for use in a microplate reader in 96-well plate, providing enough reagent to perform 100 tests for each caspase.</p>
Platform	Microplate reader

Properties

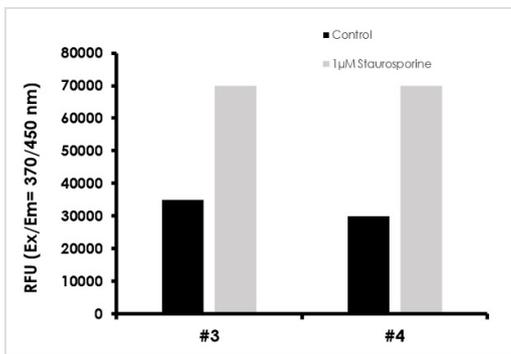
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
200X Caspase 3 Substrate	1 x 50µl
200X Caspase 8 Substrate	1 x 50µl
200X Caspase 9 Substrate	1 x 50µl
Assay Buffer	1 x 30ml

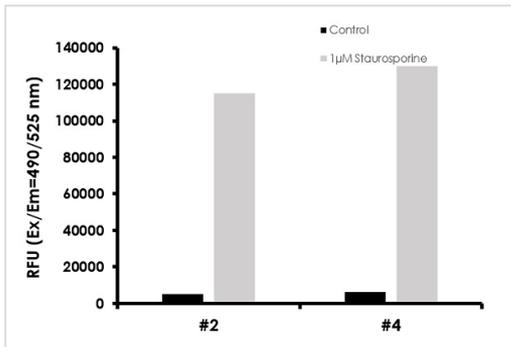
Cellular localization

Caspase 3: Cytoplasmic Caspase 8: Cytoplasmic

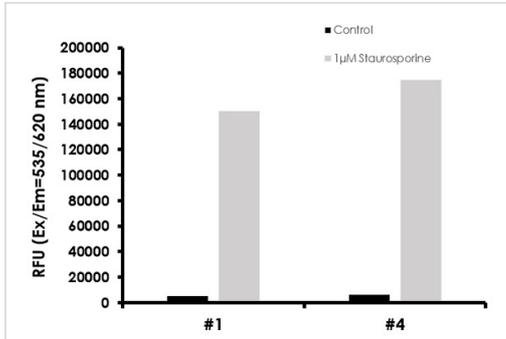
Images



Detection of Caspase activity in Jurkat cells
Caspase 9



Detection of Caspase activity in Jurkat cells
Caspase 8



Detection of Caspase activity in Jurkat cells
Caspase 3

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