

Caspase-3/7, Caspase-8 and Caspase-9 Multiplex Activity Assay Kit (Fluorometric) ab219915

[26 References](#) [3 图像](#)

概述

产品名称	Caspase-3/7, Caspase-8 and Caspase-9 Multiplex Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Semi-quantitative
产品概述	Caspase 3/7, Caspase 8 and Caspase 9 Multiplex Activity Assay Kit (Fluorometric) (ab219915) provides a simple and convenient tool to monitor caspase 3/7, caspase 8 and caspase 9 activity in cells that are undergoing apoptosis.

This product is designed to simultaneously monitor key caspases involved in apoptosis: the initiator caspases caspase 8 and caspase 9, and the executioner Caspase 3/7. The kit uses DEVD-ProRed™, IETD-R110 and LEHD-AMC as fluorogenic indicators for Caspase 3/7, 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.

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.

平台	Microplate reader
----	-------------------

性能

存放说明	Store at -20°C. Please refer to protocols.
------	--

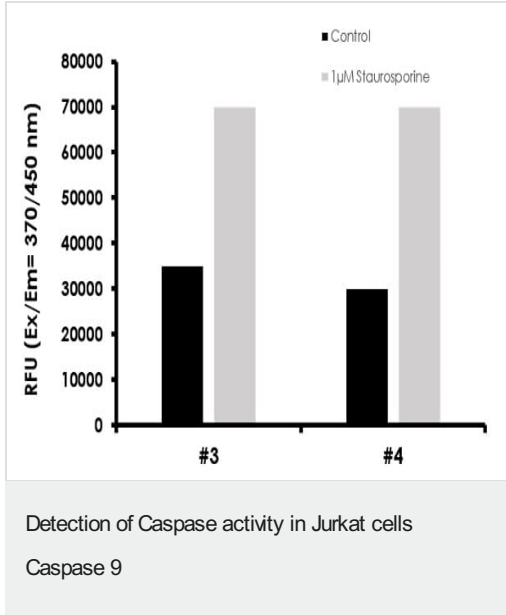
组件	100 tests
200X Caspase 3/7 Substrate	1 x 50µl
200X Caspase 8 Substrate	1 x 50µl
200X Caspase 9 Substrate	1 x 50µl

组件	100 tests
Assay Buffer	1 x 30ml

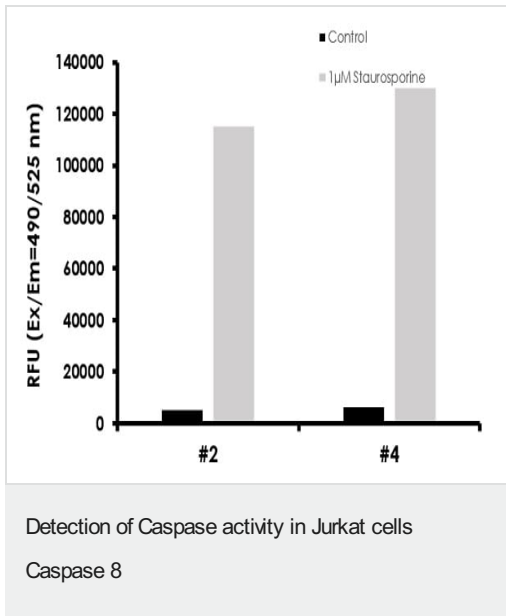
细胞定位

Caspase-7: Cytoplasm. Caspase-3: Cytoplasmic Caspase-8: Cytoplasmic

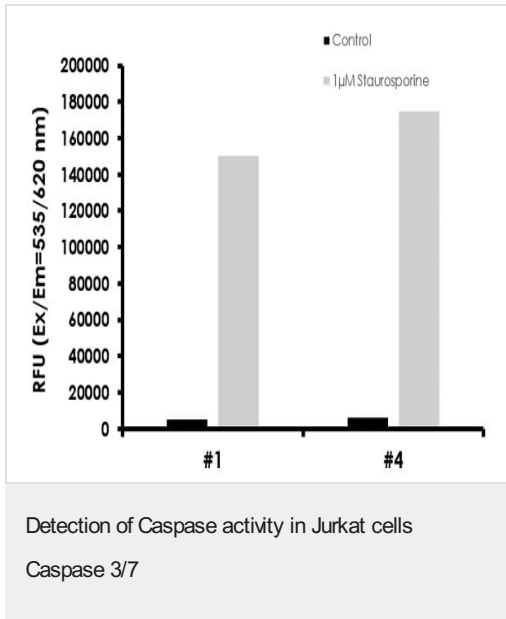
图片



Caspase 3/7, Caspase 8 and Caspase 9 Multiplex Activity Assay Kit (Fluorometric) (ab219915). Detection of Caspase activity in Jurkat cells graphs shown Caspase 9 activity. Jurkat cells were seeded on the same day at 2×10^5 cells/well in a Costar black wall/clear bottom 96-well plate. Cells were either left untreated (black bar) or treated with 1 μM staurosporine for 4 hours (gray bar). Single-caspase assay loading solution (100 μL/well; #1 for Caspase 3/7, #2 for caspase 8 or #3 for caspase 9) or Triple-caspase assay loading solution (100 μL/well; #4 for Caspase 3/7, 8 and 9 together) was added to cells, followed by an incubation at RT for 1 hour. The fluorescence intensity was measured with FlexStation fluorescence microplate reader at the indicated wavelength. Caspase 3/7, 8 and 9 activities can be detected in a single assay without interferences from other caspases.



Caspase 3/7, Caspase 8 and Caspase 9 Multiplex Activity Assay Kit (Fluorometric) (ab219915). Detection of Caspase activity in Jurkat cells graphs shown Caspase 8 activity. Jurkat cells were seeded on the same day at 2×10^5 cells/well in a Costar black wall/clear bottom 96-well plate. Cells were either left untreated (black bar) or treated with 1 μM staurosporine for 4 hours (gray bar). Single-caspase assay loading solution (100 μL/well; #1 for Caspase 3, #2 for caspase 8 or #3 for caspase 9) or Triple-caspase assay loading solution (100 μL/well; #4 for Caspase 3/7, 8 and 9 together) was added to cells, followed by an incubation at RT for 1 hour. The fluorescence intensity was measured with FlexStation fluorescence microplate reader at the indicated wavelength. Caspase 3/7, 8 and 9 activities can be detected in a single assay without interferences from other caspases.



Caspase 3/7, Caspase 8 and Caspase 9 Multiplex Activity Assay Kit (Fluorometric) (ab219915). Detection of Caspase activity in Jurkat cells graphs shown Caspase 3/7 activity. Jurkat cells were seeded on the same day at 2×10^5 cells/well in a Costar black wall/clear bottom 96-well plate. Cells were either left untreated (black bar) or treated with 1 µM staurosporine for 4 hours (gray bar). Single-caspase assay loading solution (100 µL/well; #1 for Caspase 3/7, #2 for caspase 8 or #3 for caspase 9) or Triple-caspase assay loading solution (100 µL/well; #4 for Caspase 3/7, 8 and 9 together) was added to cells, followed by an incubation at RT for 1 hour. The fluorescence intensity was measured with FlexStation fluorescence microplate reader at the indicated wavelength. Caspase 3/7, 8 and 9 activities can be detected in a single assay without interferences from other caspases.

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors