

Product datasheet

Generic Caspase Activity Assay Kit (Fluorometric - Green)
ab112130

2 References 1 图像

概述

产品名称	Generic Caspase Activity Assay试剂盒(Fluorometric - Green)
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Direct
种属反应性	与反应: Human
产品概述	<p>Abcam's Activity Assay kits are a set of tools for monitoring cellular functions. The activation of caspase is widely accepted as a reliable indicator for cell apoptosis. Most caspases have substrate selectivity for the peptide sequence Val-Ala-Asp (VAD). ab112130 uses TF2-VAD-FMK as a fluorescent indicator for most caspase activities. The cell permeable and nontoxic TF2-VAD-FMK irreversibly binds to activated caspase-1, -3, -4, -5, -6, -7, -8 and -9 in apoptotic cells. Once bound to caspases, the fluorescent reagent is retained inside the cell. The binding event prevents the caspases from further catalysis but will not stop apoptosis from proceeding. Within 15 minutes incubation, it starts to react with active caspase enzymes.</p> <p>ab112130 provides all the essential components with an optimized assay protocol. It is designed to detect cell apoptosis by measuring generic activation of caspases (caspase-1, -3, -4, -5, -6, -7, -8 and -9) in live cells.</p> <p>ab112130 is used for either the quantification of most activated caspase activities in apoptotic cells or screening of caspase inhibitors. TF2-VAD-FMK, the green label reagent, allows for direct detection of activated caspases in apoptotic cells by a flow cytometer at Ex/Em = 488/520 nm.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
说明	ab112130 should be stored dessicated.
经测试应用	适用于: Flow Cyt, Functional Studies
平台	Flow

性能

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

组件	100 tests
500 X TF5-VAD-FMK	1 x 100µl
500X Propidium Iodide	1 x 100µl
Assay Buffer	1 x 50ml

相关性 Caspases are members of the cysteine-aspartic acid protease (caspase) family. Sequential activation of caspases plays a central role in the execution-phase of cell apoptosis. Caspases exist as inactive proenzymes which undergo proteolytic processing at conserved aspartic residues to produce 2 subunits, large and small, that dimerize to form the active enzyme.

细胞定位 Cytoplasmic

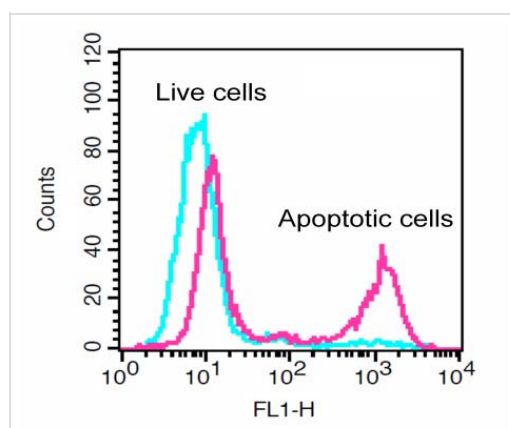
应用

Our [Abpromise guarantee](#) covers the use of **ab112130** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

应用	Ab评论	说明
Flow Cyt		Use at an assay dependent concentration.
Functional Studies		Use at an assay dependent concentration.

图片



Flow Cytometry - Generic Caspase Activity Assay Kit (Fluorometric - Green) (ab112130)

The increase in TF2-VAD-FMK (Component A) fluorescence intensity with the addition of Camptothecin in Jurkat cells.

Jurkat cells were untreated (Blue) or with 20 µM camptothecin (Pink) in a 37 °C, 5 % CO₂ incubator for 4-5 hours, and then dye loaded with TF2-VAD-FMK for 1 hour.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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