

Generic Caspase Activity Assay Kit (Fluorometric - Green) ab112130

[7 References](#) [画像数 1](#)

製品の概要

製品名	Generic Caspase Activity Assay Kit (Fluorometric - Green)
検出方法	Fluorescent
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Direct
種交差性	交差種: Mammals, Other species
製品の概要	<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>
特記事項	<p>ab112130 should be stored dessicated.</p> <p>Other caspase and apoptosis assays</p> <p>Review the full set of caspase assays, or the apoptosis assay and apoptosis marker guide.</p>
試験プラットフォーム	Flow cytometer

製品の特性

保存方法

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

内容	100 tests
500 X TF2-VAD-FMK	1 x 100µl
500X Propidium Iodide	1 x 100µl
Assay Buffer	1 x 50ml

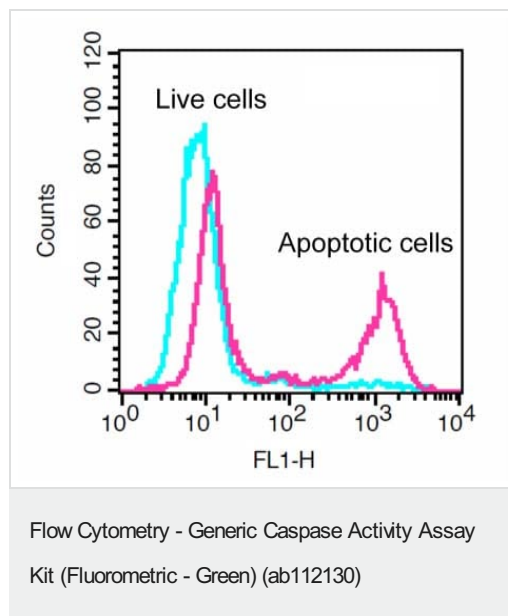
関連性

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.

細胞内局在

Cytoplasmic

画像



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. 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