

Caspase-2 Assay Kit (Fluorometric) ab39794

製品の概要

製品名 Caspase-2 Assay Kit (Fluorometric)

検出方法 Fluorescent

サンプルの種類 Tissue Extracts, Cell Lysate

アッセイタイプ Enzyme activity

全工程の試験時間 2h 00m

製品の概要

Abcam's Caspase-2 Assay Kit (Fluorometric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence VDVAAD. The assay is based on detection of cleavage of substrate VDVAAD-AFC (AFC: 7-amino-4-trifluoromethyl coumarin). VDVAAD-AFC emits blue light ($\lambda_{max} = 400 \text{ nm}$); upon cleavage of the substrate by caspase-2 or related caspases, free AFC emits a yellow-green fluorescence ($\lambda_{max} = 505 \text{ nm}$), which can be quantified using a fluorometer or a fluorescence microtiter plate reader. Visit our [FAQs page](#) for tips and troubleshooting.

特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K116 Caspase-2 Fluorometric Assay Kit. K116-100 is the same size as the 100 test size of ab39794.

Activation of ICE-family proteases/caspases initiates apoptosis in mammalian cells.

Other caspase and apoptosis assays

Review the full set of [caspase assays](#), or the [apoptosis assay and apoptosis marker guide](#).

試験プラットフォーム

Microplate reader

製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
2X Reaction Buffer	4 x 2ml
Cell Lysis Buffer	1 x 100ml
DTT	1 x 400 μ l
VDVAAD-AFC	1 x 500 μ l

機能	Involved in the activation cascade of caspases responsible for apoptosis execution. Might function by either activating some proteins required for cell death or inactivating proteins necessary for cell survival.
組織特異性	Expressed at higher levels in the embryonic lung, liver and kidney than in the heart and brain. In adults, higher level expression is seen in the placenta, lung, kidney, and pancreas than in the heart, brain, liver and skeletal muscle.
配列類似性	Belongs to the peptidase C14A family. Contains 1 CARD domain.
翻訳後修飾	The mature protease can process its own propeptide, but not that of 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.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