

Caspase-12 Assay Kit (Fluorometric) ab65664

17 References

製品の概要

製品名	Caspase-12 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-12 Assay Kit (Fluorometric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence ATAD. The assay is based on detection of cleavage of substrate ATAD-AFC (AFC: 7-amino-4-trifluoromethyl coumarin). ATAD-AFC emits blue light ($\lambda_{max} = 400 \text{ nm}$); upon cleavage of the substrate by caspase-12 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 K139 Caspase-12 Fluorometric Assay Kit. K139-100 is the same size as the 100 test size of ab65664. 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
ATAD-AFC Substrate	1 x 500 μ l
Cell Lysis Buffer	1 x 100ml
DTT	1 x 400 μ l

関連性

Caspases are cysteine proteases that cleave C-terminal aspartic acid residues on their substrate molecules. Caspase 12 is most highly related to members of the ICE subfamily. However, human caspase 12 has no protease activity. It may reduce cytokine release in response to bacterial lipopolysaccharide during infections and reduces activation of NF-kappa-B in response to TNF.

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