

Product datasheet

Caspase-12 Assay Kit (Fluorometric) ab65664

8 References

Overview

Product name	Caspase-12 Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Tissue Extracts, Cell Lysate
Assay type	Enzyme activity
Assay time	2h 00m
Product overview	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.

Notes	<p>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.</p> <p>Activation of ICE-family proteases/caspases initiates apoptosis in mammalian cells.</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> <p>Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.</p>
--------------	--

Platform	Microplate reader
-----------------	-------------------

Properties

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

Components	100 tests
2X Reaction Buffer	4 x 2ml

Components	100 tests
ATAD-AFC Substrate	1 x 500µl
Cell Lysis Buffer	1 x 100ml
DTT	1 x 400µl

Relevance

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.com/abpromise> or contact our technical team.

Terms and conditions

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