

Product datasheet

Caspase 6 Assay Kit (Fluorometric) ab39707

Overview

Product name	Caspase 6 Assay Kit (Fluorometric)
Sample type	Tissue Extracts, Cell Lysate
Assay type	Enzyme activity
Assay time	2h 00m
Product overview	<p>Abcam's Caspase 6 Assay Kit (Fluorometric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence VEID. The assay is based on detection of cleavage of substrate VEID-AFC (AFC: 7-amino-4-trifluoromethyl coumarin). VEID-AFC emits blue light ($\lambda_{max} = 400 \text{ nm}$); upon cleavage of the substrate by caspase 6 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.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
Notes	Activation of ICE-family proteases/caspases initiates apoptosis in mammalian cells.
Tested applications	Suitable for: Functional Studies

Properties

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

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

Relevance Caspase 6 is involved in the activation cascade of caspases responsible for apoptosis execution. Cleaves poly(ADP-ribose) polymerase in vitro, as well as lamins. Overexpression promotes programmed cell death. Caspase 3 and Caspase 6 are the major active caspases in apoptotic cells, and are activated in response to distinct apoptosis inducing stimuli and in all cell lines analyzed. Both Caspase 3 and 6 are present in apoptotic cells as multiple active species. Caspase 6 cleaves nuclear mitotic apparatus protein (NuMA) and mediates the shrinkage and fragmentation of nuclei.

Applications

Our [Abpromise guarantee](#) covers the use of **ab39707** in the following tested applications.

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

Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration.

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.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