

## Product datasheet

# Caspase 6 Assay Kit (Fluorometric) ab39707

### Overview

<b>Product name</b>	Caspase 6 Assay Kit (Fluorometric)
<b>Sample type</b>	Tissue Extracts, Cell Lysate
<b>Assay type</b>	Enzyme activity
<b>Assay time</b>	2h 00m
<b>Product overview</b>	<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 (<math>\lambda_{max} = 400 \text{ nm}</math>); upon cleavage of the substrate by caspase 6 or related caspases, free AFC emits a yellow-green fluorescence (<math>\lambda_{max} = 505 \text{ nm}</math>), which can be quantified using a fluorometer or a fluorescence microtiter plate reader.</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p>

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

### Properties

**Storage instructions** Store at  $-20^{\circ}\text{C}$ . Please refer to protocols.

Components	100 tests
2X Reaction Buffer	4 x 2ml
Cell Lysis Buffer	1 x 100ml
DTT	1 x 400 $\mu\text{l}$
VEID-AFC	1 x 500 $\mu\text{l}$

**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.

**Cellular localization** Cytoplasmic

**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 <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