

Caspase 2 Assay Kit (Fluorometric) ab39794

概述

产品名称	Caspase 2 Assay试剂盒(Fluorometric)
样品类型	Tissue Extracts, Cell Lysate
检测类型	Enzyme activity
检测时间	2h 00m
产品概述	<p>Abcam's Caspase 2 Assay Kit (Fluorometric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence VDVAD. The assay is based on detection of cleavage of substrate VDVAD-AFC (AFC: 7-amino-4-trifluoromethyl coumarin). VDVAD-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.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
说明	Activation of ICE-family proteases/caspases initiates apoptosis in mammalian cells.
经测试应用	适用于: Functional Studies

性能

存放说明 Store at -20°C. Please refer to protocols.

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

功能	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.
组织特异性	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.
序列相似性	Belongs to the peptidase C14A family.

Contains 1 CARD domain.

翻译后修饰

The mature protease can process its own propeptide, but not that of other caspases.

应用

Our [Abpromise guarantee](#) covers the use of **ab39794** in the following tested applications.

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

应用	Ab评论	说明
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.cn/abpromise> or contact our technical team.

Terms and conditions

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