

# Caspase 4 Assay Kit (Fluorometric) ab65658

## 1 References

### 概述

产品名称	Caspase 4 Assay试剂盒(Fluorometric)
检测方法	Fluorescent
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 4 Assay Kit (Fluorometric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence LEVD. The assay is based on detection of cleavage of substrate LEVD-AFC (AFC: 7-amino-4-trifluoromethyl coumarin). LEVD-AFC emits blue light (<math>\lambda_{\text{max}} = 400 \text{ nm}</math>); upon cleavage of the substrate by caspase 4 or related caspases, free AFC emits a yellow-green fluorescence (<math>\lambda_{\text{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>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K126 Caspase-4 Fluorometric Assay Kit. K126-100 is the same size as the 100 test size of ab65658.</p> <p>Activation of ICE-family proteases/caspases initiates apoptosis in mammalian cells.</p> <p><b>Other caspase and apoptosis assays</b></p> <p>Review the full set of <a href="#">caspase assays</a>, or the <a href="#">apoptosis assay and apoptosis marker guide</a>.</p>
平台	Microplate reader

### 性能

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

组件	100 tests
2X Reaction Buffer I	4 x 2ml
DTT I	1 x 400 $\mu$ l
LEVD-AFC	1 x 500 $\mu$ l
Lysis Buffer IV	1 x 100ml

<b>相关性</b>	Caspase 4 is a member of the cysteine-aspartic acid protease (caspase) family. It is involved in the activation cascade of caspases responsible for apoptosis execution. Caspases exist as inactive proenzymes. Activation of caspases requires proteolytic processing at conserved internal aspartic residues to generate a heterodimeric enzyme consisting of large and small subunits. Caspase 4 is able to cleave and activate its own precursor protein, as well as the caspase 1 precursor. Caspase 4 mRNA is found in most tissues examined with the exception of the brain. Highest expression is found in spleen and lung, with moderate expression in heart and liver. Low expression is observed with skeletal muscle, kidney, and testis.
<b>细胞定位</b>	Cytoplasmic and Mitochondrial

**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.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