

Caspase 9 Assay Kit (Fluorometric) ab65607

20 References 1 图像

概述

产品名称	Caspase 9 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 9 Assay Kit (Fluorometric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence LEHD. The assay is based on detection of cleavage of substrate LEHD-AFC (AFC: 7-amino-4-trifluoromethyl coumarin). LEHD-AFC emits blue light ($\lambda_{\text{max}} = 400 \text{ nm}$); upon cleavage of the substrate by caspase-9 or related caspases, free AFC emits a yellow-green fluorescence ($\lambda_{\text{max}} = 505 \text{ nm}$), which can be quantified using a fluorometer or fluorescence microtiter plate reader.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K118 Caspase-9 Fluorometric Assay Kit. K118-100 is the same size as the 100 test size of ab65607.</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>
平台	Microplate reader

性能

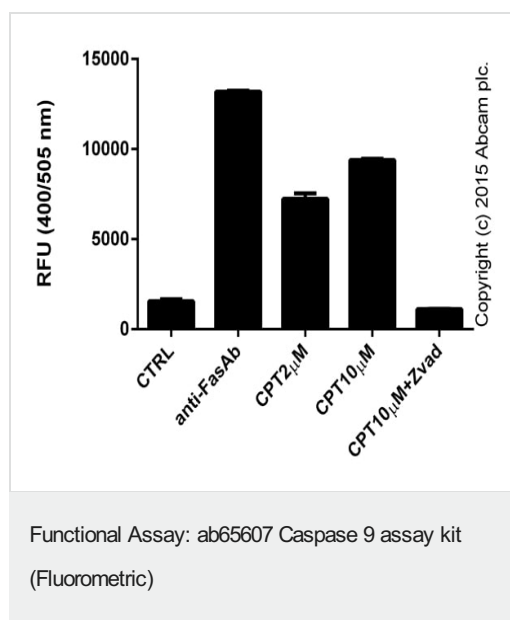
存放说明 Store at -20°C. Please refer to protocols.

组件	100 tests
2X Reaction Buffer	4 x 2ml
Cell Lysis Buffer	1 x 100ml
DTT	1 x 400μl
LEHD-AFC (1mM)	1 x 500μl

相关性

Caspases are cysteine proteases, expressed as inactive precursors, that mediate apoptosis by proteolysis of specific substrates. Caspases have the ability to cleave after aspartic acid residues. There are two classes of caspases involved in apoptosis; initiators (activation by receptor cluster) and effectors (activation by mitochondrial permeability transition). Proapoptotic signals autocatalytically activate initiator caspases, such as Caspase 8 and Caspase 9. Activated initiator caspases then process effector caspases, such as Caspase 3 and Caspase 7, which in turn cause cell collapse. Caspase 9 (also known as ICE like apoptotic protease 6 (ICE LAP6), apoptotic protease Mch6, and apoptotic protease activating factor 3 (Apaf3)) is a member of the peptidase family C14 that contains a CARD domain. It is active as a heterotetramer and has been reported to have two isoforms. ProCaspase 9 is approximately 47 kD. It is present in the cytosol and, upon activation, translocates to the mitochondria. Caspase 9 is involved in the caspase activation cascade responsible for apoptosis execution and cleaves/activates Caspase 3 and Caspase 6. It becomes activated when recruited to the Apaf1/cytochrome c complex.

图片



Caspase 9 activity in Jurkat cells following 24 hour exposure to 50 ng x mL⁻¹ anti-Fas Ab (MBL), 2 uM or 10 uM Camptothecin (CPT; [ab120115](#)) with or without 20 nM caspase inhibitor Z-VAD(OMe)-FMK (Zvad; [ab120487](#)). Sample input 5-25 uL, background signal subtracted, , duplicates; +/- SD.

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