

Caspase 9 (active) Red Staining Kit ab65619

1 References

概述

| | |
|------|--|
| 产品名称 | Caspase 9 (active) Red Staining试剂盒 |
| 检测方法 | Fluorescent |
| 样品类型 | Adherent cells, Suspension cells |
| 检测类型 | Enzyme activity |
| 检测时间 | 2h 00m |
| 产品概述 | Abcam's Caspase 9 (active) Red Staining Kit provides a convenient means for detecting activated caspase 9 in living cells. The assay utilizes the caspase 9 inhibitor LEHD-FMK conjugated to sulfo-rhodamine (Red-LEHD-FMK) as the fluorescent in situ marker. Red-LEHD-FMK is cell permeable, nontoxic, and irreversibly binds to activated caspase 9 in apoptotic cells. Visit our FAQs page for tips and troubleshooting. |
| 说明 | <p>This product is manufactured by BioVision, an Abcam company and was previously called K199 CaspGLOW™ Red Active Caspase-9 Staining Kit. K199-100 is the same size as the 100 test size of ab65619.</p> <p>Activation of caspases plays a central role in apoptosis.</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, Fluor. microscope, Flow cyt. |

性能

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

| 组件 | 100 tests |
|----------------|-----------|
| Red-LEHD-FMK | 1 x 100µl |
| Wash Buffer IV | 2 x 100ml |
| Z-VAD-FMK | 1 x 10µl |

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.

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