

Product datasheet

Caspase 9 Assay Kit (Colorimetric) ab65608

36 References 1 Image

Overview

| | |
|---------------------------|---|
| Product name | Caspase 9 Assay Kit (Colorimetric) |
| Detection method | Colorimetric |
| Sample type | Tissue Extracts, Cell Lysate |
| Assay type | Enzyme activity |
| Assay time | 2h 00m |
| Species reactivity | Reacts with: Mammals |
| Product overview | Caspase 9 Assay Kit (Colorimetric) ab65608 provides a simple and convenient means for assaying the activity of caspases that recognize the sequence LEHD. |

The caspase 9 assay is based on spectrophotometric detection of the chromophore p-nitroanilide (p-NA) after cleavage from the labeled substrate LEHD-p-NA. The p-NA light emission can be quantified using a spectrophotometer or a microtiter plate reader at 400- or 405 nm.

Notes This product is manufactured by BioVision, an Abcam company and was previously called K119 Caspase-9 Colorimetric Assay Kit. K119-100 is the same size as the 100 test size of ab65608. Activation of ICE-family proteases/caspases initiates apoptosis in mammalian cells.

Other caspase and apoptosis assays

Review the full set of [caspase assays](#), or the [apoptosis assay and apoptosis marker guide](#).

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform Microplate reader

Properties

Storage instructions Store at -20°C. Please refer to protocols.

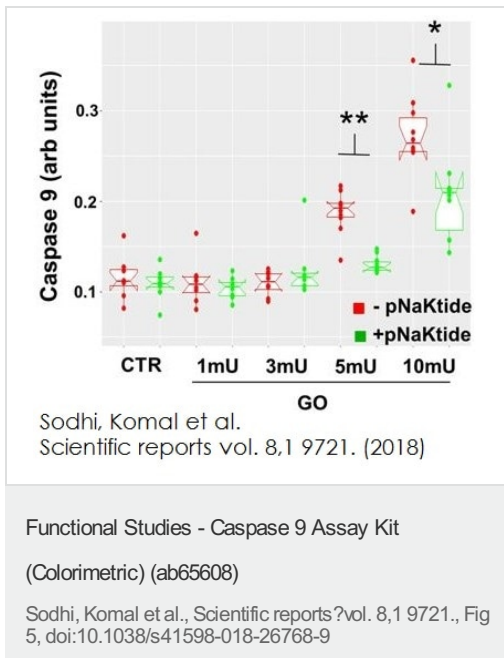
| Components | 100 tests |
|--------------------|-----------|
| 2X Reaction Buffer | 4 x 2ml |

| Components | 100 tests |
|-------------------|-----------|
| Cell Lysis Buffer | 1 x 100ml |
| Dilution Buffer | 1 x 100ml |
| DTT | 1 x 400µl |
| LEHD-pNA (4mM) | 1 x 500µl |

Relevance

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.

Images



Caspase-9 activity measured by ab65608. N = 16/group, □ p < 0.05, □□ p < 0.01 vs. corresponding GO concentration groups without pNaKtide.

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.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