

## Product datasheet

# Caspase 9 Assay Kit (Colorimetric) ab65608

### 8 References

#### 製品の概要

製品名	Caspase 9 Assay Kit (Colorimetric)
サンプルの種類	Tissue Extracts, Cell Lysate
アッセイタイプ	Enzyme activity
全工程の試験時間	2h 00m
製品の概要	<p>Abcam's Caspase 9 Assay Kit (Colorimetric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence LEHD. The 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.</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p>
特記事項	Activation of ICE-family proteases/caspases initiates apoptosis in mammalian cells.

#### 製品の特性

**保存方法** Store at -20°C. Please refer to protocols.

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

**関連性** 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 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 <https://www.abcam.co.jp/abpromise> or contact our technical team.

### **Terms and conditions**

---

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