

### Caspase 4 Assay Kit (Colorimetric) ab65659

#### 2 References

#### Overview

<b>Product name</b>	Caspase 4 Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Tissue Extracts, Cell Lysate
<b>Assay type</b>	Enzyme activity
<b>Assay time</b>	2h 00m
<b>Product overview</b>	<p>Abcam's Caspase 4 Assay Kit (Colorimetric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence LEVD. The assay is based on spectrophotometric detection of the chromophore p-nitroanilide (p-NA) after cleavage from the labeled substrate LEVD-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>
<b>Notes</b>	<p>This product is manufactured by BioVision, an Abcam company and was previously called K127 Caspase-4 Colorimetric Assay Kit. K127-100 is the same size as the 100 test size of ab65659.</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>
<b>Platform</b>	Microplate reader

#### Properties

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

Components	100 tests
2X Reaction Buffer	4 x 2ml
Cell Lysis Buffer	1 x 100ml
Dilution Buffer II	1 x 100ml
DTT	1 x 400µl

Components	100 tests
LEVD-pNA (4mM)	1 x 500µl

#### Relevance

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.

#### Cellular localization

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.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