

Caspase 5 Assay Kit (Colorimetric) ab65661

製品の概要

製品名 Caspase 5 Assay Kit (Colorimetric)

検出方法 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 Abcam's Caspase 5 Assay Kit (Colorimetric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence WEHD. The assay is based on spectrophotometric detection of the chromophore p-nitroanilide (p-NA) after cleavage from the labeled substrate WEHD-p-NA. The p-NA light emission can be quantified using a spectrophotometer or a microtiter plate reader at 400- or 405 nm. Visit our [FAQs page](#) for tips and troubleshooting.

特記事項 This product is manufactured by BioVision, an Abcam company and was previously called K123 Caspase-5 Colorimetric Assay Kit. K123-100 is the same size as the 100 test size of ab65661. 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](#).

試験プラットフォーム Microplate reader

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 II	1 x 100ml
DTT	1 x 400µl
WEHD-pNA (4mM)	1 x 500µl

関連性

Caspase 5 belongs to the Ced3/ICE family of cysteine proteases. Caspase 5 exists as an inactive proenzyme which undergoes proteolytic processing at conserved aspartic residues, leading to p20 and p10 subunits that dimerize to form the active enzyme. Sequential activation of caspases plays a central role in the execution-phase of cell apoptosis.

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.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