

Caspase-2 Assay Kit (Colorimetric) ab39830

Overview

Product name	Caspase-2 Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Tissue Extracts, Cell Lysate
Assay type	Enzyme activity
Assay time	2h 00m
Product overview	Abcam's Caspase-2 Assay Kit (Colorimetric) provides a simple and convenient means for assaying the activity of caspases that recognize the sequence VDVAD. The assay is based on spectrophotometric detection of the chromophore p-nitroanilide (p-NA) after cleavage from the labeled substrate VDVAD-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.

Notes This product is manufactured by BioVision, an Abcam company and was previously called K117 Caspase-2 Colorimetric Assay Kit. K117-100 is the same size as the 100 test size of ab39830. 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](#).

Platform 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
VDVAD-pNA	1 x 500µl

Function	Involved in the activation cascade of caspases responsible for apoptosis execution. Might function by either activating some proteins required for cell death or inactivating proteins necessary for cell survival.
Tissue specificity	Expressed at higher levels in the embryonic lung, liver and kidney than in the heart and brain. In adults, higher level expression is seen in the placenta, lung, kidney, and pancreas than in the heart, brain, liver and skeletal muscle.
Sequence similarities	Belongs to the peptidase C14A family. Contains 1 CARD domain.
Post-translational modifications	The mature protease can process its own propeptide, but not that of other caspases.

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