

Product datasheet

Caspase-3 Immunoassay Kit (Fluorometric) ab234045

Overview

Product name	Caspase-3 Immunoassay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Cell Lysate
Assay type	Quantitative
Product overview	The Caspase 3 Immunoassay Kit (Fluorometric) (ab234045) provides an effective immunosorbent enzyme assay for specific, quantitative detection of caspase 3 activity in microtiter plates. The assay utilizes caspase 3 polyclonal antibody to capture activated caspase 3 from cell lysates. Caspase substrate DEVD-AFC is then added and is cleaved proportionally to the amount of activated caspase 3 in the cell lysate. The cleavage generates free AFC which can be analyzed fluorometrically (Ex/Em = 400 nm/505 nm) using a fluorescence plate reader. The assay ensures absolute specific detection of caspase 3. Other known caspases and non-specific proteases are not detected.

Notes	<p>Activation of caspase 3 plays a key role in initiation of cellular events during apoptosis.</p> <p>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.</p> <p>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.</p>
--------------	---

Platform	Microplate reader
-----------------	-------------------

Properties

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

Components	100 tests
Cell Lysis Buffer	1 x 25ml
Coating Buffer	1 x 10ml
Anti-Caspase 3 Antibody (20X)	1 x 0.5ml
Blocking Buffer	1 x 15ml

Components	100 tests
Incubation Buffer	1 x 100ml
DTT (1M)	1 x 400µl
DEVD-AFC Substrate (1 mM)	1 x 500µl
Positive Control (rh-Caspase 3)	1 x 10 units
Microtite Plate	1 unit
Adhesive Plate Cover	1 x 2 units

Function	Involved in the activation cascade of caspases responsible for apoptosis execution. At the onset of apoptosis it proteolytically cleaves poly(ADP-ribose) polymerase (PARP) at a '216-Asp-Gly-217' bond. Cleaves and activates sterol regulatory element binding proteins (SREBPs) between the basic helix-loop-helix leucine zipper domain and the membrane attachment domain. Cleaves and activates caspase-6, -7 and -9. Involved in the cleavage of huntingtin. Triggers cell adhesion in sympathetic neurons through RET cleavage.
Tissue specificity	Highly expressed in lung, spleen, heart, liver and kidney. Moderate levels in brain and skeletal muscle, and low in testis. Also found in many cell lines, highest expression in cells of the immune system.
Sequence similarities	Belongs to the peptidase C14A family.
Post-translational modifications	Cleavage by granzyme B, caspase-6, caspase-8 and caspase-10 generates the two active subunits. Additional processing of the propeptides is likely due to the autocatalytic activity of the activated protease. Active heterodimers between the small subunit of caspase-7 protease and the large subunit of caspase-3 also occur and vice versa. S-nitrosylated on its catalytic site cysteine in unstimulated human cell lines and denitrosylated upon activation of the Fas apoptotic pathway, associated with an increase in intracellular caspase activity. Fas therefore activates caspase-3 not only by inducing the cleavage of the caspase zymogen to its active subunits, but also by stimulating the denitrosylation of its active site thiol.
Cellular localization	Cytoplasm.

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