

Product datasheet

Caspase-8 (active) FITC Staining Kit ab65614

1 References 3 Images

Overview

Product name	Caspase-8 (active) FITC Staining Kit
Sample type	Adherent cells, Suspension cells
Assay type	Enzyme activity
Assay time	2h 00m
Product overview	<p>Caspase 8 (active) FITC Staining Kit (ab65614) provides a convenient means for sensitive detection of activated caspase 8 in living cells. The assay utilizes the caspase 8 inhibitor, IETD-FMK, conjugated to FITC (FITC-IETD-FMK) as a marker. FITC-IETD-FMK is cell permeable, non-toxic, and irreversibly binds to activated caspase 8 in apoptotic cells. The FITC label allows detection of activated caspase-8 in apoptotic cells directly by fluorescence microscopy, flow cytometry, or fluorescence plate reader.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
Notes	<p>Activation of caspases plays a central role in apoptosis.</p> <p>Other caspase and apoptosis assays</p> <p>Review the full set of caspase assays, or the apoptosis assay and apoptosis marker guide.</p>
Platform	Microplate reader, Fluor. microscope, Flow cyt.

Properties

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

Components	100 tests
FITC-IETD-FMK	1 x 100µl
Wash Buffer	2 x 100ml
Z-VAD-FMK	1 x 10µl

Function Most upstream protease of the activation cascade of caspases responsible for the TNFRSF6/FAS mediated and TNFRSF1A induced cell death. Binding to the adapter molecule FADD recruits it to either receptor. The resulting aggregate called death-inducing signaling complex (DISC) performs CASP8 proteolytic activation. The active dimeric enzyme is then liberated from the DISC and free to activate downstream apoptotic proteases. Proteolytic fragments of the N-terminal propeptide (termed CAP3, CAP5 and CAP6) are likely retained in the

DISC. Cleaves and activates CASP3, CASP4, CASP6, CASP7, CASP9 and CASP10. May participate in the GZMB apoptotic pathways. Cleaves ADPRT. Hydrolyzes the small-molecule substrate, Ac-Asp-Glu-Val-Asp-

-AMC. Likely target for the cowpox virus CRMA death inhibitory protein. Isoform 5, isoform 6, isoform 7 and isoform 8 lack the catalytic site and may interfere with the pro-apoptotic activity of the complex.

Tissue specificity

Isoform 1, isoform 5 and isoform 7 are expressed in a wide variety of tissues. Highest expression in peripheral blood leukocytes, spleen, thymus and liver. Barely detectable in brain, testis and skeletal muscle.

Involvement in disease

Defects in CASP8 are the cause of caspase-8 deficiency (CASP8D) [MIM:607271]. CASP8D is a disorder resembling autoimmune lymphoproliferative syndrome (ALPS). It is characterized by lymphadenopathy, splenomegaly, and defective CD95-induced apoptosis of peripheral blood lymphocytes (PBLs). It leads to defects in activation of T-lymphocytes, B-lymphocytes, and natural killer cells leading to immunodeficiency characterized by recurrent sinopulmonary and herpes simplex virus infections and poor responses to immunization.

Sequence similarities

Belongs to the peptidase C14A family.

Contains 2 DED (death effector) domains.

Domain

Isoform 9 contains a N-terminal extension that is required for interaction with the BCAP31 complex.

Post-translational modifications

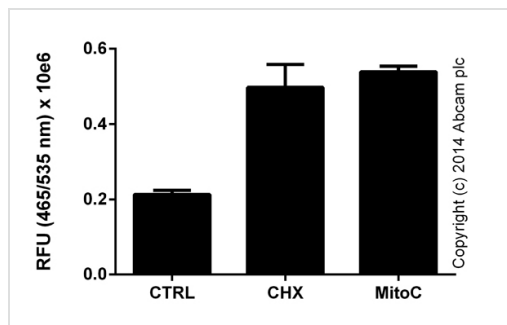
Generation of the subunits requires association with the death-inducing signaling complex (DISC), whereas additional processing is likely due to the autocatalytic activity of the activated protease. GZMB and CASP10 can be involved in these processing events.

Phosphorylated upon DNA damage, probably by ATM or ATR.

Cellular localization

Cytoplasm.

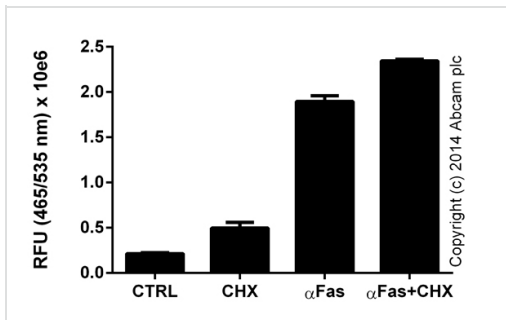
Images



Active caspase 8 in control Jurkat cells (10e6/mL) or cells treated for five hours with 10 ug/mL Cyclohexamide (CHX) ([ab120093](#)) or four hours with 25 ug/mL Mitomycin C (MitoC) ([ab120797](#)).

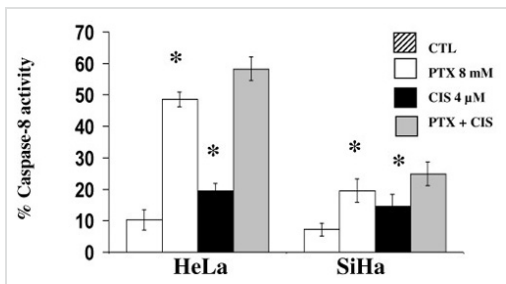
Background signal subtracted, duplicates; +/- SD.

Functional assays: Caspase 8 (active) FITC Staining Kit (ab65614)



Active caspase 8 in Jurkat cells following four hours exposure to 50 ng/mL anti-Fas Ab (α Fas) (MBL), five hours with 10 μ g/mL cyclohexamide (CHX) ([ab120093](#)), or one hour pretreatment with CHX followed by four hours with α Fas. Background signal subtracted, duplicates; +/- SD.

Functional assays: Caspase 8 (active) FITC Staining Kit (ab65614)



Caspase 8 activation of HeLa (left) and SiHa (right) cells after *in vitro* treatment with pentoxyllyne (PTX) or cisplatin (CIS) either alone or in combination. Results represent the mean \pm SD of three independent experiments carried out in triplicate. (*) $p < 0.001$ vs CTL.

Image obtained from Hernandez-Flores G et al; BMC Cancer, 2011 Nov 11; 11:483

Flow Cytometry - Caspase 8 (active) FITC Staining Kit (ab65614)

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.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