

Cytochrome c Apoptosis ICC Antibody Kit ab110417

[9 References](#) [1 Image](#)

Overview

Product name	Cytochrome c Apoptosis ICC Antibody Kit
Detection method	Fluorescent
Sample type	Adherent cells
Assay type	Direct
Species reactivity	Reacts with: Mouse, Rat, Human
Product overview	Apoptotic cells which have released mitochondrial cytochrome c into the cytosol can be differentiated from non apoptotic cells which still retain cytochrome c in their mitochondria by fluorescence microscopy.

ab110417 (MSA07) contains a cytochrome c monoclonal antibody (50 µg) and an isotype-specific secondary antibody (50 µg) which has a fluorescein fluorophore conjugated to it. The kit also contains an ATP synthase V subunit alpha monoclonal antibody (50 µg) and the appropriate isotype-specific secondary antibody (50 µg) which has a Texas Red fluorophore conjugated to it.

ATP synthase is an inner mitochondrial membrane protein and serves as a mitochondrial marker since it is not released into the cytoplasm during apoptosis, unlike cytochrome c. Thus one is able to observe cytochrome c and a mitochondrial marker (ATP synthase subunit alpha) simultaneously by fluorescence microscopy using ab110417. Buffers and blocking solution for immunocytochemical analysis are also included in the kit.

Notes This product was previously called ApoTrack™ Cytochrome c Apoptosis ICC Antibody Kit.

Other apoptosis assays

For more apoptosis assays, review the [apoptosis assay and apoptosis marker guide](#).

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.

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.

Tested applications **Suitable for:** ICC/IF

Properties

Storage instructions Please refer to protocols.

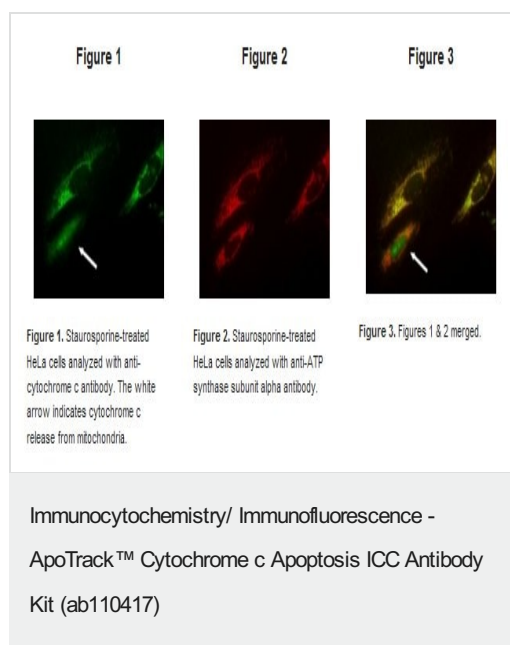
Components	Identifier	200 µg
100% normal goat serum		1 x 20ml
10X Buffer (10X PBS, pH 7.4)		1 x 100ml
Anti-Complex Va monoclonal antibody	clone 15H4C4	1 x 50µg
Anti-Cytochrome c monoclonal antibody	clone 37BA11	1 x 50µg
Detergent (0.1% (v/v) Triton® X-100 in PBS)		1 x 30ml
Goat anti-mouse IgG2a - FITC secondary antibody		1 x 50µg
Goat anti-mouse IgG2b - TXRD secondary antibody		1 x 50µg

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab110417 in the following tested applications. The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
ICC/IF		Use at an assay dependent concentration.

Images



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