

# Cytochrome c Apoptosis ICC Antibody Kit ab110417

[9 References](#) [画像数 1](#)

### 製品の概要

---

製品名	Cytochrome c Apoptosis ICC Antibody Kit
検出方法	Fluorescent
サンプルの種類	Adherent cells
アッセイタイプ	Direct
種交差性	交差種: Mouse, Rat, Human
製品の概要	<p>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.</p> <p>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.</p> <p>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.</p>
特記事項	<p>This product was previously called ApoTrack™ Cytochrome c Apoptosis ICC Antibody Kit.</p> <p><b>Other apoptosis assays</b></p> <p>For more apoptosis assays, review the <a href="#">apoptosis assay and apoptosis marker guide</a>.</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. 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>
アプリケーション	適用あり: ICC/IF

### 製品の特性

---

保存方法 Please refer to protocols.

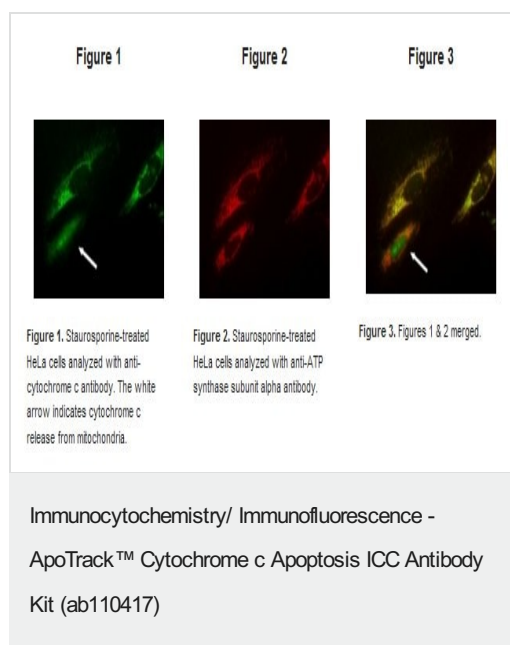
内容	ラベル	200 µg
100% normal goat serum		1 x 20ml
10X Buffer (10X PBS, pH 7.4)		1 x 100ml
Anti-Complex Va monoclonal antibody	clone 15H4C4	1 x 50µg
Anti-Cytochrome c monoclonal antibody	clone 37BA11	1 x 50µg
Detergent (0.1% (v/v) Triton® X-100 in PBS)		1 x 30ml
Goat anti-mouse IgG2a - FITC secondary antibody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

## アプリケーション

**The Abpromise guarantee** **Abpromise保証は、** 次のテスト済みアプリケーションにおけるab110417の使用に適用されます  
 アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご確認ください。

アプリケーション	Abreviews	特記事項
ICC/IF		Use at an assay dependent concentration.

## 画像



**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