

# Cytochrome c Apoptosis ICC Antibody Kit ab110417

**9 References**   [1 图像](#)

### 概述

产品名称	Cytochrome c Apoptosis ICC抗体试剂盒
检测方法	Fluorescent
样品类型	Adherent cells
检测类型	Direct
种属反应性	与反应: Mouse, Rat, Human
产品概述	<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>
说明	<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.</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>
经测试应用	适用于: ICC/IF

### 性能

**存放说明** Please refer to protocols.

组件	标识符	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

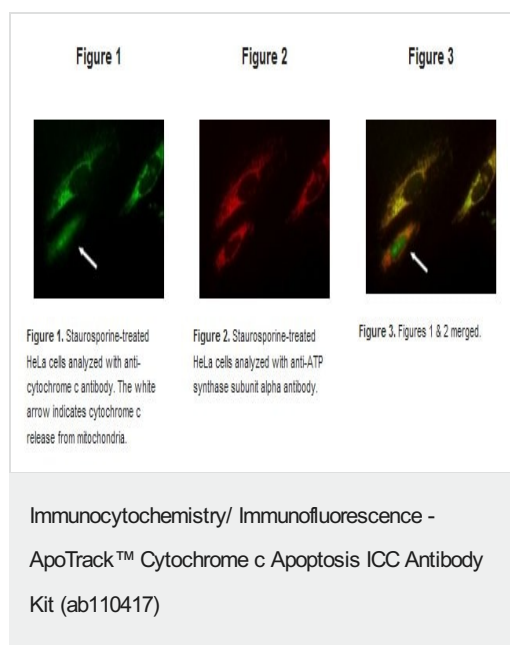
## 应用

**The Abpromise guarantee** **Abpromise™**承诺保证使用ab110417于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
ICC/IF		Use at an assay dependent concentration.

## 图片



**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.cn/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors