

# Mouse and Rabbit Specific HRP/AEC Detection IHC Kit ab93705

★★★★★ [1 Abreviews](#) [14 References](#) [2 图像](#)

### 概述

#### 产品名称

小鼠和兔Specific HRP/AEC Detection IHC试剂盒

#### 产品概述

Mouse and Rabbit specific HRP/AEC IHC Detection Kit ab93705 is a complete kit to detect a primary antibody raised in either mouse or rabbit, using HRP-labeled-streptavidin and a biotinylated anti-rabbit/anti-mouse secondary antibody.

It is supplied with a full protocol and with peroxidase and protein blocking reagents, biotinylated secondary antibody, streptavidin-HRP, and AEC substrate.

In the method:

- a sample is incubated with an unconjugated primary antibody specific to the target antigen
- a biotinylated secondary antibody then binds to the primary antibody
- HRP-labeled streptavidin then binds to the secondary antibody
- the HRP produces a colored substance at the site of primary antibody binding by reacting with AEC.

#### 说明

For alternative AEC streptavidin-biotin IHC detection kits see:

- anti-Rabbit HRP-AEC IHC kit [ab64260](#)
- Anti-Mouse HRP-AEC IHC Kit [ab64258](#)

For high sensitivity AEC HRP micro-polymer IHC detection kits see:

- Anti-Mouse / anti-Rabbit HRP-AEC Micro-polymer IHC Kit [ab236467](#)
- Anti-Rabbit HRP-AEC Micro-polymer IHC Kit [ab236468](#)

Alternatively, see [HRP-DAB IHC detection kits](#).

Find more kits and reagents for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the [IHC kits and reagents guide](#).

#### 经测试应用

适用于: IHC-P

### 性能

#### 存放说明

Store at +4°C. Please refer to protocols.

组件	15 ml
----	-------

组件	15 ml
AEC Single Solution	1 x 15ml
Biotinylated Goat Anti-Polyvalent	1 x 15ml
Hydrogen Peroxide Block	1 x 15ml
Protein block	1 x 15ml
Streptavidin Peroxidase	1 x 15ml

**相关性** An HRP/AEC Detection Kit is a labeled immunoenzymatic antigen detection system. Horseradish peroxidase (HRP) readily combines with hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>) and the resultant complex can oxidize a wide variety of chromogenic substrates. AEC is 3-amino-9-ethyl-carbazole, a chromogenic substrate.

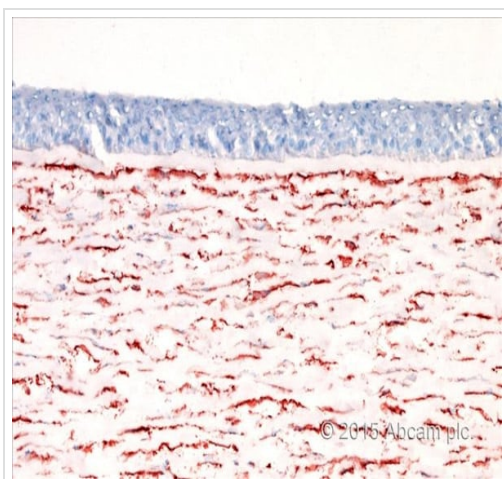
## 应用

**The Abpromise guarantee** **Abpromise™** 承诺保证使用 ab93705 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
IHC-P		Use at an assay dependent dilution.

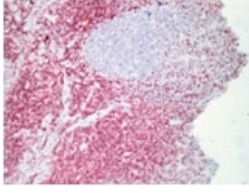
## 图片



Human cornea (cryosections fixed in cold acetone) stained with anti-CD34 antibody using ab93705 Mouse and Rabbit Specific HRP/AEC detection IHC kit.

CD34 detection using HRP/AEC Detection Kit  
(ab93705)

This image is courtesy of an Abreview submitted by  
Stanislava Reinstein Merjava



Human tonsil stained with anti-CD3 antibody using ab93705  
Polyvalent HRP/AEC Detection Kit (Ready-To-Use).

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Mouse and Rabbit Specific  
HRP/AEC detection IHC kit (ab93705)

**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