

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

★★★★★ [1 Abreviews](#) [14 References](#) [画像数 2](#)

### 製品の概要

製品名	Mouse and Rabbit Specific HRP/AEC Detection IHC Kit
製品の概要	<p>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.</p> <p>It is supplied with a full protocol and with peroxidase and protein blocking reagents, biotinylated secondary antibody, streptavidin-HRP, and AEC substrate.</p> <p>In the method:</p> <ul style="list-style-type: none"><li>- a sample is incubated with an unconjugated primary antibody specific to the target antigen</li><li>- a biotinylated secondary antibody then binds to the primary antibody</li><li>- HRP-labeled streptavidin then binds to the secondary antibody</li><li>- the HRP produces a colored substance at the site of primary antibody binding by reacting with AEC.</li></ul>
特記事項	<p>For alternative AEC streptavidin-biotin IHC detection kits see:</p> <ul style="list-style-type: none"><li>- anti-Rabbit HRP-AEC IHC kit <a href="#">ab64260</a></li><li>- Anti-Mouse HRP-AEC IHC Kit <a href="#">ab64258</a></li></ul> <p>For high sensitivity AEC HRP micro-polymer IHC detection kits see:</p> <ul style="list-style-type: none"><li>- Anti-Mouse / anti-Rabbit HRP-AEC Micro-polymer IHC Kit <a href="#">ab236467</a></li><li>- Anti-Rabbit HRP-AEC Micro-polymer IHC Kit <a href="#">ab236468</a></li></ul> <p>Alternatively, see <a href="#">HRP-DAB IHC detection kits</a>.</p> <p>Find more kits and reagents for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the <a href="#">IHC kits and reagents guide</a>.</p>
アプリケーション	<b>適用あり:</b> IHC-P

### 製品の特性

**保存方法** Store at +4°C. Please refer to protocols.

内容	15 ml
----	-------

内容	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

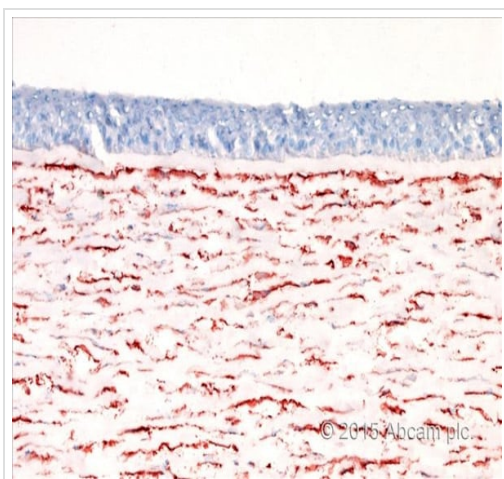
**関連性** An HRP/AEC Detection Kit is a labeled immunoenzymatic antigen detection system. Horseradish peroxidase (HRP) readily combines with hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>) and the resultant complex can oxidize a wide variety of chromogenic substrates. AEC is 3-amino-9-ethyl-carbazole, a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

### 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ab93705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検討ください。

アプリケーション	Abreviews	特記事項
IHC-P		Use at an assay dependent dilution.

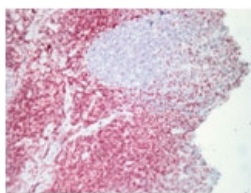
### 画像



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.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