

# Rabbit specific HRP/AEC IHC Detection Kit - Micro-polymer ab236468

4 References 1 图像

概述

产品名称	兔specific HRP/AEC IHC Detection试剂盒- Micro-polymer
检测方法	Colorimetric
检测类型	Micro-Polymer
产品概述	<p>This IHC detection kit is used to detect rabbit primary antibodies in immunohistochemistry (IHC). It uses an HRP micro-polymer secondary antibody for highly sensitive detection.</p> <p>HRP micro-polymer methods are more sensitive than the traditional streptavidin-biotin method, and they do not require a blocking step for endogenous biotin.</p> <p>The kit includes ready-to-use hydrogen peroxide blocking reagent, protein blocking reagent, HRP micro-polymer conjugated goat anti-rabbit secondary antibody, and AEC chromogen.</p>
说明	<p>For alternative kits, see Anti-Mouse / anti-Rabbit HRP-AEC Micro-polymer IHC Kit <a href="#">ab236467</a>, and <a href="#">HRP-DAB micro-polymer IHC 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>
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
AEC Single Solution	1 x 15ml
Goat anti-rabbit HRP Conjugate	1 x 15ml
Hydrogen Peroxide Block	1 x 15ml
Protein block	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.

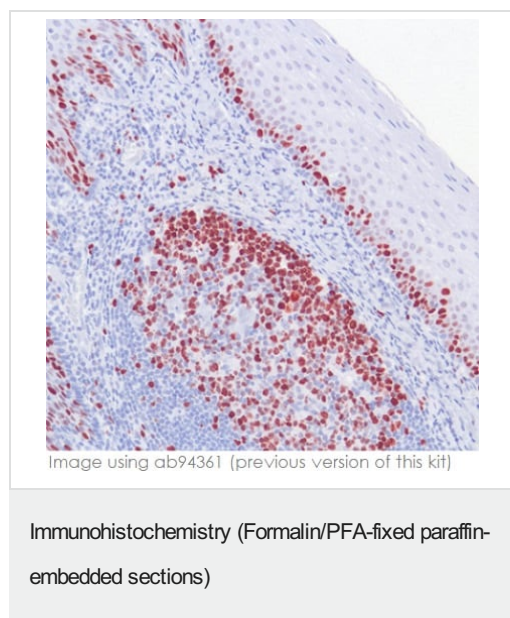
## 应用

**The Abpromise guarantee** Abpromise™ 承诺保证使用 ab236468 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 concentration.

## 图片



Immunohistochemical analysis of Human Tonsil tissue labeling Ki-67 using the [ab94361](#) anti-EXPOSE Rabbit specific HRP-AEC detection IHC kit (this kit is identical to [ab94361](#), except that it uses the HRP micro-polymer secondary antibody used by Abcam's R&D team for antibody validation in IHC).

**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